

EpiCypher® CUTANA™ CUT&RUN Protocol

For histone PTMs, transcription factors (TFs), and chromatin regulators

This protocol has been validated for genomic profiling of:

- Histone PTMs (e.g. lysine methylation, acetylation and ubiquitylation)
- Transcription factors (e.g. CTCF, FOXA1)
- Chromatin remodelers (e.g. ATPase subunits of SWI/SNF, ISWI, INO80, CHD)
- Chromatin writers & readers (e.g. MLL1, BRD4)
- Nuclear hormone receptor signaling factors (e.g. Estrogen Receptor)
- Epitope-tagged proteins (e.g. HA, FLAG tags)

Table of Contents

1. Overview	2
2. CUTANA™ Products & Services: Advantages.....	3
3. Outline of CUT&RUN Workflow	4
4. Experimental Design & Key Protocol Notes.....	6
5. Buffers, Reagents & Materials Needed.....	9
6. EpiCypher CUTANA™ CUT&RUN Protocol	13
Section I: ConA Bead Activation (~30 min)	13
Section II: Binding Cells to Activated Beads (~30 min)	13
Section III: Binding of Antibodies (~30 min + overnight).....	14
Section IV: Binding of pAG-MNase (~30 min)	16
Section V: Targeted Chromatin Digestion and Release (~3 hrs)	16
Section VI: Library Preparation (~4 hrs)	17
Section VII: Agilent 2100 Bioanalyzer System (~1 hr).....	18
Section VIII: Illumina Sequencing	18
7. Frequently Asked Questions (FAQs)	20
8. References	27
Appendix I: Nuclei Isolation & Cryopreservation Protocol for CUT&RUN.....	28
Appendix II: Sample Integrity & Bead Conjugation Quality Control Checks	30
Appendix III: SNAP-CUTANA™ Spike-in Controls.....	32
Appendix IV: Use of <i>E. coli</i> Spike-in DNA for Experimental Normalization	47

*EpiCypher® and SNAP-ChIP® are registered trademarks of EpiCypher, Inc.
All other trademarks and trade names in this document are property of their respective corporations.*

1. Overview

Cleavage Under Targets & Release Using Nuclease (CUT&RUN) is a revolutionary genomic mapping strategy developed by the group of Dr. Steven Henikoff¹. It builds on Chromatin ImmunoCleavage (ChIC) from Dr. Ulrich Laemmli², whereby a fusion of Proteins A and/or G to Micrococcal Nuclease (pAG-MNase) is recruited to selectively cleave antibody-bound chromatin *in situ*³. In CUT&RUN, cells or nuclei are immobilized to a solid support, with pAG-MNase cleaved target DNA fragments isolated from solution. The workflow is compatible with next-generation sequencing (NGS) to provide high resolution genome-wide profiles of histone post-translational modifications (PTMs) and chromatin-associated proteins (e.g. TFs and chromatin remodelers; **Figure 1**).

The following protocol describes detailed recommendations for performing CUT&RUN, based on extensively optimized workflows developed by EpiCypher scientists. Check back at epicypher.com/protocols for regular protocol updates.

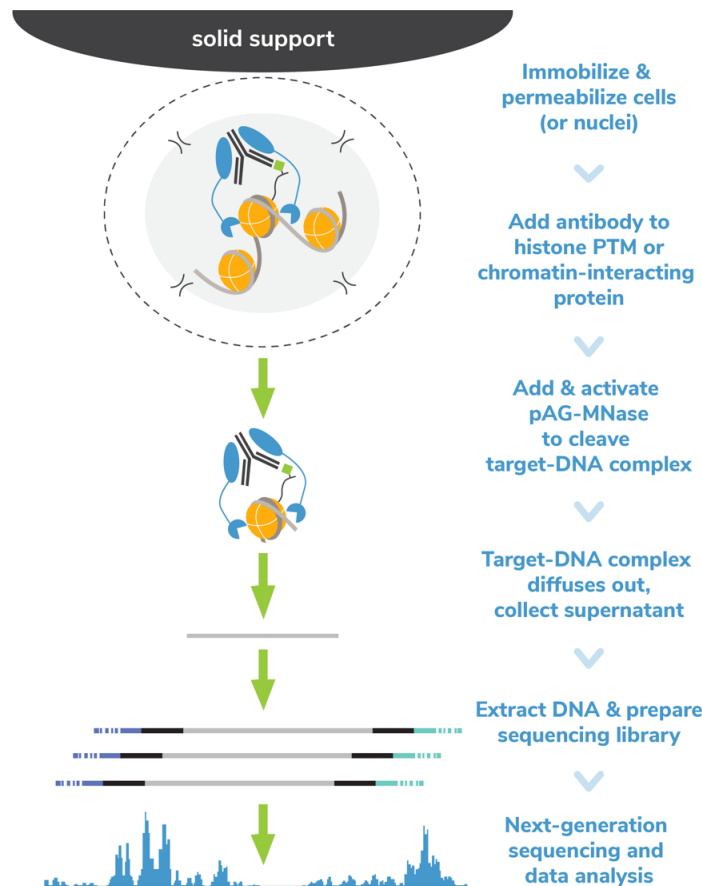


Figure 1. Overview of the CUTANA™ CUT&RUN protocol.

2. CUTANA™ Products & Services: Advantages

ChIC and CUT&RUN have revolutionized the study of chromatin regulation. Compared to ChIP-seq (the current leading approach for genome-wide mapping of histone PTMs and chromatin-associated proteins), CUT&RUN offers the following advantages:

Platform Comparison	ChIP-seq	CUTANA™ CUT&RUN
Required cells	> 1 million	5,000 – 500,000
Cell Input	Fragmented chromatin	Intact cells or nuclei
Compatible Targets	Histone PTMs & chromatin-associated proteins	Histone PTMs & chromatin-associated proteins, including difficult ChIP targets
Sequencing Depth (Reads)	> 30 million	3-8 million
Experimental throughput	Low	High
Signal : Noise	Low	High
Assay Automation	Difficult	Yes

EpiCypher now offers a suite of products to support CUT&RUN workflows under our CUTANA™ assay platform, including:

- **pAG-MNase** ([EpiCypher 15-1016](#)), the essential reagent and first-in-class commercial product for ChIC/CUT&RUN assays. The EpiCypher enzyme contains an optimized fusion of Proteins A and G with Micrococcal Nuclease (pAG-MNase) to enable compatibility with a broad range of antibody isotypes.
- **ChIC/CUT&RUN Kit** ([EpiCypher 14-1018](#)) with all reagents included to go from cells to purified CUT&RUN DNA.
- **CUT&RUN Compatible Antibodies**, rigorously lot-to-lot validated directly in CUT&RUN. See epicypher.com/cut-and-run-compatible-antibodies for more information.
- **CUT&RUN Spike-in Controls**
 - *E. coli* Spike-in DNA ([EpiCypher 18-1401](#)) for data normalization.
 - **SNAP-CUTANA™ Spike-in Controls** are DNA-barcoded semi-synthetic/recombinant nucleosome spike-in panels that control for all aspects of CUT&RUN workflows, including antibody specificity and assay success, guide troubleshooting experiments, and enable quantitative normalization (**Appendix III**). Now available for histone lysine methylation PTMs (SNAP-CUTANA K-MetStat Panel, [EpiCypher 19-1002](#)).
- **CUT&RUN supporting reagents** selected and validated for optimal performance in the EpiCypher CUT&RUN protocol. See epicypher.com/cut-and-run-assays for more info.
- **CUTANA CUT&Tag** reagents for ultra-low input applications. Go to epicypher.com/CUT&Tag for more information.
- Inquire for more information or to connect with EpiCypher scientists: info@epicypher.com

3. Outline of CUT&RUN Workflow

Description: EpiCypher’s in-house optimized protocol for CUTANA™ CUT&RUN assays (**Figure 2**). Before starting, we strongly recommend reading this section, the **Experimental Design & Key Protocol Notes** section and the **Protocol** to carefully plan your experiment and familiarize yourself with the assay.

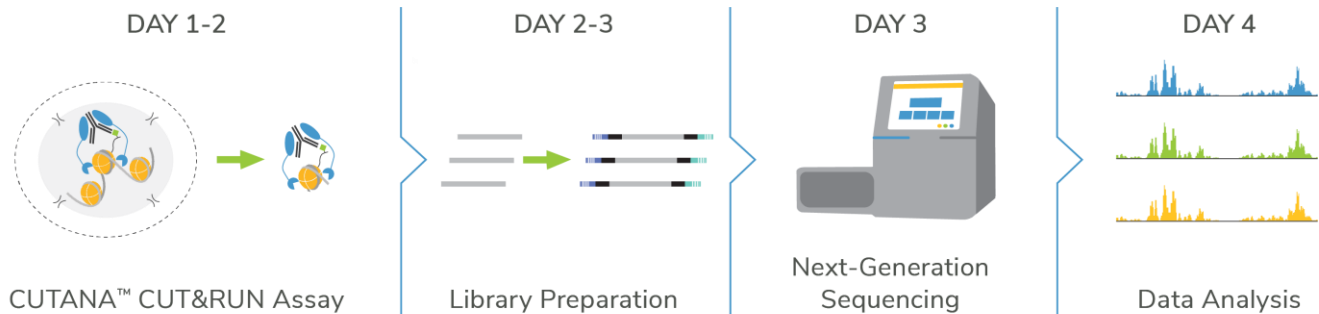


Figure 2: Timeline of CUT&RUN assay using EpiCypher’s CUTANA™ CUT&RUN protocol.

Section I: ConA Bead Activation (~30 min)

Description: This section “activates” Concanavalin A coated beads (**Con A beads**) to bind and immobilize cell/nuclei samples. If preparing for multiple CUT&RUN reactions, it is recommended to batch process the full volume of beads needed for all reactions in a single 1.5 mL tube. This helps ensure homogeneity across reactions. For these steps, beads can be washed using a 1.5 mL magnetic rack (e.g. [EpiCypher 10-0012](#)).

Section II: Binding Cells to Activated Beads (~30 min)

Description: In this section, cell/nuclei samples are prepared and immobilized to ConA beads in bulk, and then separated into 8-strip PCR tubes for individual CUT&RUN reactions. This protocol section was developed using 500,000 unfixed (*i.e.* native) K562 cells per CUT&RUN reaction and is specifically designed for batch processing of cell samples for multiple CUT&RUN reactions (see **Protocol Notes**). Once the cell-coupled beads are transferred to PCR tubes, it is recommended to use multi-channel pipettes and a compatible magnetic rack ([EpiCypher 10-0008](#); **Figure 3**), which helps increase experimental throughput and reproducibility.

Importantly, we have protocol adaptations for many types of inputs (adherent vs. suspension cells, nuclei, cryopreservation, and cross-linking) noted within **Section II** of the protocol and our **FAQs**. To ensure that cells are immobilized to ConA beads, we have also developed simple quality control checks (**Appendix II**) that we strongly recommend using in every experiment.

Section III: Binding of Antibodies (~30 min + overnight)

Description: After coupling to activated ConA beads, the cell – bead mixture is resuspended in cold **Antibody Buffer**, and a target-specific antibody is added for an overnight incubation. Note that antibody selection in CUT&RUN is critical to success; see **FAQs** for more information.

Importantly, **Antibody Buffer** is the first to contain digitonin (see **Buffer Recipes**), which permeabilizes cells and allows the antibody to bind its target *in situ* (histone PTM or chromatin-associated protein). The digitonin concentration required for CUT&RUN assays varies by sample (e.g. cell type, fixation) and must be optimized for every unique input, as described in the **Protocol Notes**. It is crucial to use the minimum amount of digitonin needed for efficient permeabilization to avoid cell lysis and digitonin precipitation during overnight incubations.

A second feature of this section is the addition of the **SNAP-CUTANA™ K-MetStat Panel** ([EpiCypher 19-1002](#)) to CUT&RUN reactions using a methyl-lysine antibody (e.g. H3K4me3) or IgG control. This spike-in panel comprises highly pure, DNA-barcoded semi-synthetic/recombinant nucleosomes carrying defined lysine methylation PTMs, and is useful for in-assay antibody validation, quantitative normalization, and experimental troubleshooting. We outline its appropriate use in the **Section III** of the protocol (*i.e.* added immediately before antibody addition; also see **Table 5**). **Appendix III** includes detailed information about how the spike-ins work and how they can be incorporated as critical controls to master CUT&RUN.

Section IV: Binding of pAG-MNase (~30 min)

Description: At this stage of the protocol, **pAG-MNase** ([EpiCypher 15-1016](#)) is added to each reaction and incubated briefly at room temperature to allow binding to antibody-labelled chromatin. This incubation is performed in the **Digitonin Buffer**, which is free of CaCl₂ to avoid premature activation of MNase. ConA bead – cell suspensions often become “clumpy” at this point and are difficult to pipette. Taking care to minimize this clumping is key for quality results. In this case, we recommend using a P200 pipette to gently disperse beads in buffer. Cut-off pipette tips can be used if beads are clogging pipette tips or if cells/nuclei are easily damaged.

Section V: Targeted Chromatin Digestion and Release (~3 hrs)

Description: During this part of the experiment, MNase is activated by addition of CaCl₂ to cleave and release antibody-bound chromatin. The fragmented chromatin containing the histone PTM or protein of interest diffuses into the supernatant, where it can be easily separated from bead-coupled cells. The fragmented chromatin is purified, quantified using a Qubit™ fluorometer, and used for next-generation sequencing (NGS) library preparation. Of note, the **CUTANA™ DNA Purification Kit** ([EpiCypher 14-0050](#)) is designed to capture CUT&RUN DNA fragments as small as 50 bp, which is adequate for most chromatin targets, including TFs.

We have also included instructions for the addition of exogenous ***E. coli* Spike-in DNA** ([EpiCypher 18-1401](#)), which can be added to CUT&RUN reactions following MNase activation as a component of the **Stop Buffer** and used to normalize downstream NGS data. For further instructions on optimizing *E. coli* DNA for CUT&RUN experiments, see **Appendix IV**.

IMPORTANT: At this step, many researchers are tempted to assess fragment size distribution of CUT&RUN purified DNA and/or analyze enrichment of known targets by qPCR. We do not recommend either of these analyses, which are derived from ChIP-seq and are not robust

indicators of CUT&RUN success (see **FAQs** for a detailed explanation). CUT&RUN is distinct from CHIP-seq, and the quality controls applied in one assay cannot be transferred to the other. In fact, raw CUT&RUN DNA yields are often below the sensitivity of Bioanalyzer/TapeStation approaches used to examine fragment distribution. Furthermore, because CUT&RUN is performed *in situ*, there is no chromatin Input, complicating enrichment analysis by qPCR.

The best indicator of CUT&RUN success at this step is that raw DNA yields from positive control reaction(s) are greater than yields from the IgG negative control. For low abundance targets, such as H3K4me3, this difference may be slight, while high abundance targets (e.g. H3K27me3) will display a more pronounced difference. Proceed directly to library preparation.

Section VI: Library Preparation (~4 hrs)

Description: Here, we provide guidance on CUT&RUN sequencing library preparation using the NEBNext® Ultra™ II Library Prep Kit for Illumina®. This includes PCR steps specifically optimized to amplify short CUT&RUN library fragments compatible with Illumina paired-end sequencing (100 bp – 700 bp including universal adapters; see **FAQs** for further guidance).

Section VII: Agilent 2100 Bioanalyzer System (~1 hr)

Description: Prior to sequencing, CUT&RUN libraries should be examined using the Bioanalyzer, TapeStation, or equivalent approach to assess fragment size, library concentration, and DNA quality. Predominant enrichment of mononucleosome-sized fragments (~300 bp including universal adapters) is the best indicator of CUT&RUN experimental success prior to NGS (e.g. see **Figure 5**; also see **FAQs** for expanded discussion on quality control checks for CUT&RUN). Here we describe the use of Bioanalyzer and the Agilent High Sensitivity DNA Chip for CUT&RUN library analysis along with expected results.

Section VIII: Illumina Sequencing

Description: The final step of the protocol is loading and sequencing your CUT&RUN libraries on an Illumina sequencing system. In this section, we provide guidance on pooling indexed CUT&RUN libraries for multiplexed sequencing. We also offer recommendations for loading pooled libraries onto MiniSeq and NextSeq systems. These suggestions are based on the fact that CUT&RUN only requires 3-8 million reads per sample to generate high signal-to-noise data (vs. 30 million or more with CHIP-seq). This allows researchers to pool >48 samples per sequencing run, if using the Illumina NextSeq, or use a benchtop sequencer (e.g. Illumina MiniSeq) for smaller projects.

For guidelines on NGS analysis see the **FAQ** section. For guidance on SNAP-CUTANA K-MetStat spike-in analysis, see **Appendix III**; for *E. coli* spike-in DNA, see **Appendix IV**.

4. Experimental Design & Key Protocol Notes

Description: This section is considered essential reading for CUTANA™ CUT&RUN assays. For CUT&RUN workflows to be successful you must include proper controls and optimize key steps

for your unique cell input (e.g. digitonin permeabilization, number of cells) as detailed in this section. We also offer tips on common problems with the protocol and explain our rationale for using 8-strip PCR tubes in the CUTANA CUT&RUN protocol.

1. Include controls in every experiment. We suggest the following controls (at minimum):
 - Negative control antibody (e.g. IgG negative control antibody: [EpiCypher 13-0042](#))
 - Positive control antibody (e.g. for PTMs, H3K4me3: [EpiCypher 13-0041](#); for chromatin-associated protein targets, BRD4: [EpiCypher 13-2003](#) or CTCF [EpiCypher 13-2014](#))
 - SNAP-CUTANA™ Spike-in Controls ([EpiCypher 19-1002](#); see **Appendix III**). They should be included with H3K4me3 and IgG control antibodies in every experiment, as well as any assays targeting histone methyl-lysine PTMs.

These controls are especially critical when optimizing CUT&RUN for new experimental conditions (e.g. new cell types, reduced inputs, drug treatments, fixation method), but should also be included as standard controls for continuous monitoring of assay success.

2. Optimize conditions and become familiar with the CUT&RUN workflow using a control cell line (e.g. K562 cells) before attempting different sample types (see **FAQs**).
 - This protocol was optimized using 500,000 human K562 cells per reaction. However, without any further modifications, this protocol has been validated on as few as 5,000 cells with antibodies against H3K4me3 ([EpiCypher 13-0041](#)), H3K27me3 ([Thermo Fisher MA5-11198](#)), and BRD4 ([EpiCypher 13-2003](#)).
3. We strongly recommend performing the **Quality Control Checks** outlined in **Appendix II**, which include evaluating the integrity of starting cells/nuclei and binding to Concanavalin A (**ConA**) beads.
4. While the standard protocol is written for native (unfixed) suspension cells, we have included specific sample preparation instructions for working with nuclei (**Appendix I**), cryopreserved cells/nuclei (**Appendix I**), adherent cells (**CUT&RUN Protocol, Section II**), and cross-linked material (see **Cross-linking Protocol** at epicypher.com/protocols).
5. We recommend a 5% digitonin stock solution be prepared in DMSO (as opposed to heated H₂O), as this improves detergent solubility and protocol reproducibility.
6. To avoid digitonin precipitation and cell lysis, use the minimal concentration of digitonin required for efficient permeabilization in the **Digitonin Buffer** and **Antibody Buffer**. This step should be optimized for every type of cell/nuclei input used in CUT&RUN workflows. A detailed protocol for digitonin optimization is described in the CUTANA™ ChIC/CUT&RUN Kit manual ([EpiCypher 14-1048](#)). In general, perform a digitonin titration (e.g. 3-fold dilutions from 0.1% down to 0.001%) and monitor cell lysis using Trypan blue staining (see **Appendix II**). Find the minimum digitonin concentration needed to achieve >95% permeabilized cells.

7. This protocol has been adapted to 8-strip PCR tubes (vs. 1.5 mL tubes) for rapid “batch processing” of multiple CUT&RUN reactions from bulk cell samples. The first few steps, *i.e.* cell harvest and conjugation to ConA beads, are performed in 1.5 mL tubes, and then the reactions are split into 8-strip PCR tubes for the remainder of the assay. These steps:
 - Minimize beads sticking to tubes
 - Enable more rapid workflows with multichannel pipettes
 - Provide more consistent sample handling
 - Enable high-throughput sample preparation
8. ConA beads dry out easily, which can result in sample loss. To avoid this problem in the CUT&RUN assay, take caution to prevent ConA beads sticking to the sides/caps of tubes.
 - To avoid ConA beads sticking to tube sides/caps and drying out, it is essential to use a **nutator** rather than a **rotator** (see **Table 3**), since nutators gently agitate by shaking or rocking tubes rather than rotating end-over-end.
 - Take note of steps that indicate when to pipette or vortex to disperse clumps and keep ConA beads in an even suspension.
9. Although protocols with shortened antibody and/or CaCl₂ incubation times have been published³, in our hands such changes adversely impact yields and reproducibility.
10. **IMPORTANT:** Since CUT&RUN has lower background and is compatible with fewer cells compared to ChIP-seq, **it is not recommended to assess fragment size distribution using agarose gel or capillary electrophoresis prior to library preparation.** Such analysis is not indicative of the success of a CUT&RUN experiment, and further the amount of DNA recovered is often below the sensitivity of detection for these approaches (see **FAQs**). Instead, assess DNA yield compared to positive (*e.g.* H3K4me3, BRD4, CTCF) and negative (IgG) controls, determine fragment size distribution of sequencing-ready libraries (**Figure 5**), and evaluate peak structure and expected genome-wide distribution in sequencing data.

5. Buffers, Reagents & Materials Needed

Table 1: Buffer components

Components	Source	Cat #
HEPES	Sigma-Aldrich	H3375
KCl	Sigma-Aldrich	P3911
CaCl ₂	Sigma-Aldrich	C1016
MnCl ₂	Sigma-Aldrich	203734
Molecular biology grade H ₂ O (RNase, DNase free)	VWR	VWRV02-0201-0500
NaCl	Sigma-Aldrich	S5150-1L
EDTA (prepare 0.5 M stock at pH 8.0)	Sigma-Aldrich	E5134
EGTA (prepare 0.5 M stock at pH 8.0)	Sigma-Aldrich	E3889
RNase A	Thermo Fisher Scientific	EN0531
Glycogen	Sigma-Aldrich (Roche)	10930193001
Spermidine trihydrochloride*	Sigma-Aldrich	S2501
Digitonin (store aliquots of 5% stock in DMSO at -20°C)	Millipore Sigma	300410
DMSO	Sigma-Aldrich	D8418-100ml
cOmplete™, Mini, EDTA-free Protease Inhibitor Cocktail	Roche	11836170001
Trypan blue	Thermo Fisher Scientific	T10282

*1M spermidine preparation: Dissolve 1 gram spermidine (MW = 254.63) in 3.93 mL molecular grade H₂O. Store in single-use aliquots at -20°C for 6 months.

Buffer recipes

Bead Activation Buffer

20 mM HEPES, pH 7.9
10 mM KCl
1 mM CaCl₂
1 mM MnCl₂
Filter sterilize. Store at 4°C for up to 6 months.

Pre-Wash Buffer

20 mM HEPES, pH 7.5
150 mM NaCl
Filter sterilize. Store at 4°C for up to 6 months.

Wash Buffer

Pre-Wash Buffer (recipe above)
0.5 mM Spermidine*
1x Roche cOmplete™, Mini, EDTA-free Protease Inhibitor (CPI-mini, 1 tab/10mL)
Filter sterilize. Store at 4°C for up to 1 week.

Digitonin Buffer

Wash Buffer + 0.01% digitonin**
Prepare fresh each day and store at 4°C.

Antibody Buffer

Digitonin Buffer*** + 2 mM EDTA
Prepare fresh each day and store at 4°C.

Stop Buffer

340 mM NaCl
20 mM EDTA
4 mM EGTA
50 µg/mL RNase A
50 µg/mL Glycogen
Filter sterilize. Store at 4°C for up to 6 months.

Buffer Preparation Notes

* Spermidine is added to compensate for the removal of Mg²⁺ from the buffer. Mg²⁺ can cause DNA degradation and is typically omitted from CUT&RUN buffers.

** Optimal [digitonin] for each cell type should be determined empirically, as described in **Protocol Notes**. Starting concentration validated for K562, MCF7, and A549 cells is 0.01% digitonin.

*** See **Protocol Notes** for guidance on optimizing [digitonin] for a given cell type.

Table 2: Reagents

Item	Vendor	Catalog No.	Notes
Concanavalin A (ConA) Conjugated Paramagnetic Beads	EpiCypher	21-1401	ConA is a lectin, which can cause immune cell activation. For technical support re. immune cell studies, see the FAQs section or contact info@epicypher.com .
CUTANA™ pAG-MNase	EpiCypher	15-1016 or 15-1116	50 & 250 reaction pack sizes available. Supplied as 20X stock.
SNAP-CUTANA™ K-MetStat Panel	EpiCypher	19-1002	<p>Spike-in controls for Sample Normalization & Antibody Profiling (SNAP Spike-in Controls): The K-MetStat panel comprises fifteen semi-synthetic/recombinant nucleosomes carrying unique methyl-lysine modifications (me1/2/3 at H3K4, H3K9, H3K27, H3K36, and H4K20), plus an unmodified control, immobilized to magnetic beads. Each histone octamer is wrapped with two different barcoded DNA templates, providing an internal technical replicate for each histone PTM.</p> <p>Add spike-ins to CUT&RUN reactions targeting one of the PTMs in the panel as well as in CUT&RUN samples designated for H3K4me3 positive and IgG negative control antibodies. For more information about using SNAP-CUTANA Spike-ins, see Appendix III.</p>
Rabbit IgG Negative Control Antibody	EpiCypher	13-0042	See datasheet for application-specific dilutions.
SNAP-ChIP® Certified, CUTANA Compatible H3K4me3 Positive Control Antibody	EpiCypher	13-0041	See datasheet for application-specific dilutions.
SNAP-ChIP® Certified H3K27me3 Positive Control Antibody	Thermo Fisher	MA5-11198	Use 0.5 µg per reaction.
CUTANA Compatible CTCF Positive Control Antibody	EpiCypher	13-2014	See datasheet for application-specific dilutions.
Antibody to target of interest (histone PTM, TF, or chromatin regulator)	User-dependent		EpiCypher continues to conduct <u>extensive</u> antibody characterization studies ⁴ . For antibodies directly validated in CUT&RUN, visit: epicypher.com/cut-and-run-compatible-antibodies . For other targets not found on our webpage, contact us for recommendations: info@epicypher.com .
<i>E. coli</i> Spike-in DNA	EpiCypher	18-1401	Use as an exogenous spike-in control for experimental normalization. See Appendix IV for detailed instructions.
CUTANA™ DNA Purification Kit	EpiCypher	14-0050	Designed to maximize yield from low [DNA] samples/low cell inputs; is optimized for low elution volume and retention of >50bp fragments. Suitable for both nucleosomal (PTMs) and subnucleosomal (TFs) fragments.
Agencourt AMPure XP Magnetic Beads	Beckman Coulter	A63880	For PCR cleanup and size selection (<i>i.e.</i> adapter dimer and primer removal).
Qubit™ 1x dsDNA HS Assay Kit	Thermo Fisher Scientific	Q33230	For DNA quantification.
NEBNext® Ultra™ II DNA Library Prep Kit for Illumina®	New England Biolabs	E7645S	For DNA sequencing applications.

Table 3: Equipment

Item	Vendor	Catalog No.	Notes
1.5 mL Magnetic Separation Rack	EpiCypher	10-0012	For bulk or “batch” processing of ConA beads and ConA bead-conjugated cells in Section I and II of CUT&RUN protocol; see Figure 3A .
8-strip PCR tube Magnetic Separation Rack	EpiCypher	10-0008	For processing of individual CUT&RUN reactions in Section III onward; see Figure 3B . Enables streamlined sample handling for higher experimental throughput and improved reproducibility.
8-strip 0.2 mL PCR tubes	EpiCypher	10-0009	Compatible with the magnetic stand.
Qubit™ 4 Fluorometer	Thermo Fisher Scientific	Q33226	For DNA quantification.
Agilent 2100 Bioanalyzer	Agilent	G2939A	For analysis of purified CUT&RUN sequencing libraries. May substitute comparable capillary electrophoresis instrument (e.g. Agilent TapeStation).
High Performance Multi-Channel Pipettors, 8-Channel	VWR	76169-252	For performing CUT&RUN in 8-strip PCR tubes e.g. for aspiration and wash steps. May substitute comparable multi-channel pipettor.
Tube Nutator	VWR	82007-202	For bead incubation steps (overnight antibody incubation, pAG-MNase digest reaction). It is critical to use a tube nutator rather than a rotator for these steps.
Vortex-Genie	Scientific Industries	SI-0236	For bead mixing steps.

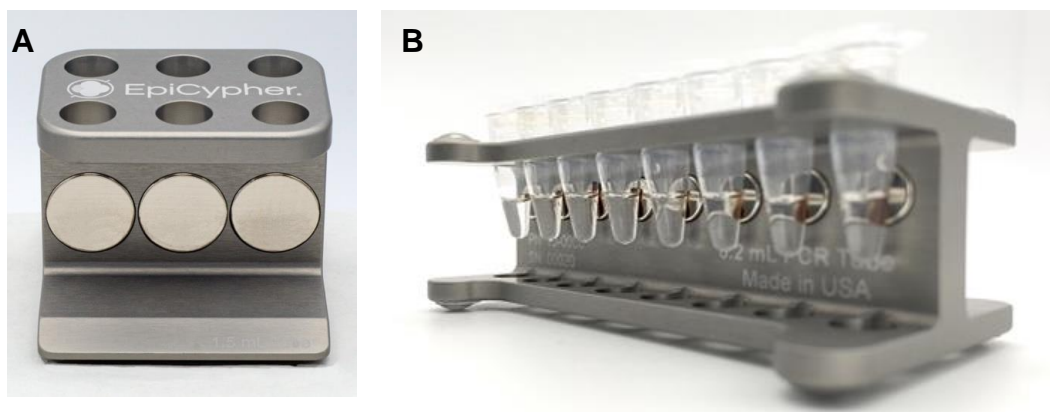


Figure 3: Magnetic racks for CUT&RUN assays. **(A)** For batch processing of ConA beads, use a 1.5 mL magnetic rack (e.g. EpiCypher 10-0012, pictured). **(B)** For processing samples in 8-strip PCR tubes we recommend using a multi-channel pipette and compatible magnetic rack (e.g. EpiCypher 10-0008, pictured).

6. EpiCypher CUTANA™ CUT&RUN Protocol

Essential Reading: Before starting, we strongly recommend reading the **Outline of CUT&RUN Workflow** and the **Experimental Design & Key Protocol Notes** for important information about assay controls and optimization. The **FAQs section** also provides guidance on adapting the protocol for unique inputs, targets, and other cell preparation methods.

6.1. CUT&RUN Protocol (~5hrs)

---Day 1---

Section I: ConA Bead Activation (~30 min)

1. Gently resuspend the **ConA beads** and transfer 11 μL per planned CUT&RUN reaction to a 1.5 mL tube for batch processing.
2. Place the tube on a 1.5 mL magnetic separation rack until slurry clears (30 s – 2 min) and pipette to remove supernatant.
3. Immediately add 100 μL /reaction cold **Bead Activation Buffer**, remove from magnet, and pipette to mix. Place back on magnet until slurry clears and pipette to remove supernatant.
4. Repeat the previous step for total of two washes.
5. Resuspend beads in 11 μL /reaction cold **Bead Activation Buffer**.
6. Keep activated **ConA beads** on ice until needed.

Section II: Binding Cells to Activated Beads (~30 min)

7. Harvest 500,000 cells per planned CUT&RUN reaction. If more than one CUT&RUN reaction is planned for the cell sample, cells can be processed together at this step (see **Table 4** for guidance). For suspension cell cultures, spin for 3 min at 600 x g at room temperature (**RT**) in a 1.5 mL tube. Remove supernatant by decanting or pipetting.

Notes and alternative cell preparation protocols:

- It is recommended to prepare excess cells (~10% excess if working in batch [preferred] or one extra sample if preparing individual reactions) for Quality Control (QC) Checks described in **Appendix II**. See **Table 4** for guidance.
- To harvest adherent cells, we recommend a very mild trypsin treatment: 0.05% trypsin, incubated at 37°C, for the minimal time necessary to detach cells (see **FAQs**).
- Freshly isolated nuclei, frozen nuclei, and frozen cells are compatible with CUT&RUN; for instructions, see **Appendix I**.
- For samples that require cross-linking (e.g. formaldehyde) to preserve labile marks, such as histone lysine acetylation, see our cross-linking protocol at epicypher.com/protocols.
- See **FAQs “Sample Input Compatibility”** section for special considerations when using adherent, cryopreserved and cross-linked cells, immune cells, and tissue.

Cell Type	# Reactions	Cells/Reaction	Total Cells*	Vol. Wash Buffer*	Vol. ConA Beads*
K562 cells	8	500,000	4.4 million	880 µL	88 µL
K562 cells	1	500,000	1 million	200 µL	20 µL

Table 4: Preparation of cells for CUT&RUN reactions with extra volume allotted for key Quality Control Checks (**Appendix II**). **Total cell number and volumes include a 10% excess for batch processing cells, or one extra sample volume for preparing individual CUT&RUN reactions. Extra rows are provided to customize suggestions for user workflows.*

8. Resuspend cells in 100 µL/reaction RT **Wash Buffer**, spin for 3 min at 600 x g at RT, and remove supernatant by decanting or pipetting.
9. Repeat the previous step for total of two washes.
10. Resuspend cells in 100 µL/reaction in RT **Wash Buffer**. Pipette to mix.
11. Proceed with **ConA bead** binding.
 - **Note 1:** If batch processing cell samples for multiple CUT&RUN reactions (as in **Table 4, row 1**), process bulk cells + ConA beads together in 1.5 mL tubes to ensure homogeneity across reactions. Add 10 µL of activated **ConA beads** per 100 µL washed cells. Gently vortex (setting #7) to mix.
 - **Note 2:** If working with individual cell samples at this step (e.g. each CUT&RUN reaction uses a distinct cell type or sample input; see **Table 4, row 2**), add 10 µL activated **ConA beads** per 100 µL cells, and transfer individual samples (110 µL) to 8-strip PCR tubes for subsequent steps.
12. **Incubate** cell – bead slurry for **10 min at RT**. Cells will adsorb to the activated ConA beads.

Section III: Binding of Antibodies (~30 min + overnight)

13. Place the tube(s) on a magnet until slurry clears (30 s – 2 min). Pipette to remove and discard supernatant.
14. Add 50 µL cold **Antibody Buffer** per reaction quickly, to avoid bead drying. Gently vortex immediately and thoroughly to an even resuspension.
 If still working in 1.5 mL tubes for batch processing:
 - Take into account 10% excess volume of Antibody Buffer as in **Table 4**. Can save an aliquot of cell – bead slurry for QC checks described in **Appendix II**.
 - Transfer individual reactions (50 µL) into 8-strip PCR tubes.
 - Continue the remaining steps using an 8-strip PCR tube magnetic rack.
15. For reactions designated for positive control histone PTM (e.g. H3K4me3) and negative control (IgG) antibodies, as well as samples assigned a target in the K-MetStat Panel (me1, me2, and me3 at H3K4, H3K9, H3K27, H3K36 and H4K20): Add 2 µL **SNAP-CUTANA™ K-MetStat Panel** per 500,000 cells. If using less than 500,000 cells, decrease the amount of

SNAP-CUTANA Spike-ins linearly by preparing a “working stock” dilution of the panel in **Antibody Buffer**. General starting recommendations are provided in **Table 5**.

- **Note 1:** SNAP-CUTANA Spike-ins must be added before addition of Primary Antibody.
- **Note 2:** Aim for spike-ins to comprise ~1% of total sequencing reads; adjust amount added as needed. It is acceptable for this bandwidth to be higher for low abundance targets and negative controls (e.g. IgG = 10-20%) and lower for high abundance targets (e.g. H3K27me3 = 0.1-1%).
- See **Appendix III** for more information about SNAP-CUTANA Spike-ins.

Starting # Cells in CUT&RUN	Working Stock Dilution in Antibody Buffer [use <i>FRESH</i> the day of preparation]	Volume added to reaction	Final dilution in reaction
500,000	Stock	2 µL	1:25
250,000	1:2	2 µL	1:50
100,000	1:5	2 µL	1:125
50,000 or less*	1:10	2 µL	1:250

Table 5: Recommended SNAP-CUTANA Spike-in amounts for varying numbers of starting cells in CUT&RUN. ***NOTE:** *additional dilutions of the SNAP-CUTANA Panels may be added for lower cell inputs. However, dilution of spike-ins beyond 1:250 increases risk of experimental variation and may require end user optimization.*

16. Add manufacturer’s recommended amount (or 0.5 µg if application is untested) of **Primary Antibody** to each reaction and gently vortex immediately and thoroughly.

- **Note:** Antibodies stored in glycerol solution may be viscous. Take care to ensure accurate pipetting by aspirating slowly, check tip for accuracy, and pipette up and down ~3x times into CUT&RUN samples to clear remaining glycerol from tip.

17. **Incubate** 8-strip PCR tubes on nutator **overnight at 4°C**.

- **Critical step:** To keep beads in solution, slightly elevate cap side of 8-strip PCR tubes on nutator to ensure bead solution remains in bottom of conical tube (**Figure 4**). **DO NOT USE** a rotator or turn tubes over end-to-end for this step.



Bead slurry before O/N incubation at 4°C | Nutator O/N | Bead slurry after O/N incubation at 4°C

Figure 4. ConA beads should appear homogenous and not clumpy throughout the procedure. Vortex or pipette as needed to disperse clumps. Ensure tube caps are elevated during incubation steps on tube nutator to keep beads in solution.

---Day 2---

Section III (continued)

18. Place the 8-strip PCR tubes on a magnet until slurry clears (30 s – 2 min) and pipette to remove and discard supernatant.
19. Keeping tubes on the magnet, add 250 µL cold **Digitonin Buffer** directly onto beads of each reaction, and then pipette to remove supernatant.
20. Repeat the previous step for total of two washes, keeping tubes + beads on magnet the entire time.
21. After second wash, remove supernatant and discard. Add 50 µL cold **Digitonin Buffer** per reaction, and gently vortex the 8-strip PCR tubes to resuspend bead – cell slurry in buffer. Continue to the addition of pAG-MNase.
 - Note: Beads are often clumpy at this point but can easily be dispersed by gentle pipetting with a P200 pipette. A slightly cut-off pipette tip may be used to aid in resuspension and/or preserve delicate cells/nuclei.

Section IV: Binding of pAG-MNase (~30 min)

22. Add 2.5 µL **CUTANA pAG-MNase** (20x stock) to each CUT&RUN reaction, and gently vortex to mix.
 - Critical step: To evenly distribute pAG-MNase across cells/nuclei, ensure beads are thoroughly resuspended by gentle vortexing and/or pipetting using a P200 pipette (cut-off pipette tip optional).
23. **Incubate** reactions for **10 min at RT**.
24. Return 8-strip PCR tubes to magnet, allow slurry to clear (30 s – 2 min). Pipette to remove and discard supernatant.
25. Keeping tubes on the magnet, add 250 µL cold **Digitonin Buffer** directly onto beads of each reaction, and then pipette to remove supernatant.
26. Repeat the previous step for total of two washes, keeping tubes + beads on magnet the entire time.
27. After second wash, remove supernatant and discard. Remove 8-strip PCR tubes from the magnet and add 50 µL cold **Digitonin Buffer** to each reaction. Gently vortex to mix.
 - Note: If beads are clumpy, gently pipette to mix using a P200 pipette (cut-off pipette tip optional).

Section V: Targeted Chromatin Digestion and Release (~3 hrs)

28. Place 8-strip PCR tubes on ice and add 1 µL **100 mM CaCl₂** to each reaction. Gently vortex to mix.
 - Critical step: For efficient pAG-MNase digestion, ensure the beads are thoroughly resuspended by vortexing and/or pipetting using a P200 pipette (cut-off pipette tip optional).
29. **Incubate** 8-strip PCR tubes on nutator for **2 hours at 4°C**, with tubes slightly elevated as in

Figure 4. Note that this is the key step wherein MNase tethered to antibody-bound chromatin is activated by calcium to cleave target chromatin.

30. Add 33 μ L **Stop Buffer** to each reaction, and gently vortex to mix. This buffer stops MNase enzymatic activity by chelating Ca^{2+} ions.
 - Note: If using *E. coli* spike-in DNA for NGS normalization, prepare a **Stop Buffer Master Mix** by adding the optimized amount of ***E. coli* DNA (Appendix IV)** to the Stop Buffer immediately before adding to reactions. Make enough Master Mix for all reactions plus 10% excess volume. Mix thoroughly and add 33 μ L **Stop Buffer Master Mix** per reaction.
31. **Incubate** 8-strip PCR tubes for **10 min at 37°C** in a thermocycler. This step releases cleaved chromatin to supernatant and degrades RNA.
32. Perform a quick spin of 8-strip PCR tubes in benchtop microfuge to collect beads/buffer. Place 8-strip PCR tubes on a magnet stand until slurry clears (30 s – 2 min).
33. Transfer supernatant containing CUT&RUN-enriched DNA to new 1.5 mL tubes. The ConA beads can now be discarded.
34. Purify DNA from supernatant using the **CUTANA™ DNA Purification Kit** and provided instructions. This kit captures DNA fragments >50 bp, which is adequate for most targets.
35. Elute DNA in 12 μ L **Elution Buffer** and use 1 μ L to quantify the CUT&RUN-enriched DNA using the **Qubit™ fluorometer** per manufacturer's instructions.
 - Note 1: Yields are influenced by a variety of factors, including cell type, antibody, and target abundance. The best indicator of experimental success at this stage is that the CUT&RUN DNA yield of the target(s) of interest are greater than IgG negative control (even if slightly so, *i.e.* for low abundance targets such as H3K4me3).
 - Note 2: EpiCypher's typical CUT&RUN DNA yields from 500,000 native K562 cells (pre-PCR) are noted below:
 - IgG \rightarrow ~ 2-5 ng
 - H3K4me3 (low abundance target) \rightarrow ~5-10 ng
 - H3K27me3 (high abundance target) \rightarrow ~20-50 ng
36. **PROCEED DIRECTLY TO LIBRARY PREPARATION.** DO NOT assess fragment size distribution prior to library preparation, as the yields may be below the limit of detection for this analysis and is not indicative of experimental success. See **FAQs** for more information.

6.2. Library prep (4 hrs), Bioanalyzer (1 hr), & Illumina Sequencing

Section VI: Library Preparation (~4 hrs)

37. Use 5-10 ng purified CUT&RUN-enriched DNA to prepare Illumina NGS libraries using the **NEBNext® Ultra™ II Library Prep Kit for Illumina®** per manufacturer's instructions.
 - Note: The NEB Ultra™ II Kit contains hot-start Q5 DNA polymerase.
38. Overview of library prep workflow, according to manufacturer's instructions:
 - a. End repair
 - b. Universal adapter ligation

- c. DNA cleanup using 1.1x **AMPure beads** : reaction volume (retains fragments >150 bp)
 - d. PCR and primer indexing according to cycling parameters below
39. CUT&RUN-specific PCR cycling parameters. See **FAQs** for additional information and suggestions.
- a. 45 sec at 98°C → activation of hot-start Q5 polymerase
 - b. 15 sec at 98°C → DNA melting
 - c. 10 sec at 60°C → hybrid primer annealing & short extension (<700 bp)
 - d. Repeat Steps b-c for a total of 14 cycles
 - e. 1 min at 72°C → final extension
40. DNA cleanup using 1.1x **AMPure beads** : reaction volume (e.g. 55 µL beads : 50 µL PCR).
41. Elute DNA in 15 µL **0.1x TE buffer**. Use 1 µL to quantify the purified PCR product using the **Qubit fluorometer** per manufacturer's instructions.
- Note: Typical yield of purified PCR DNA library: ~500-750 ng (30-50 ng/µL in 15 µL)

Section VII: Agilent 2100 Bioanalyzer System (~1 hr)

42. For each purified CUT&RUN library, including your IgG negative control, prepare 5 µL at 10 ng/µL for loading on the Agilent Bioanalyzer. 1 µL of each diluted library will be loaded onto the **Agilent High Sensitivity DNA Chip** (Cat# 5067-4626).
- Note the dilution factor, which will be needed to calculate the library molarity for desired DNA size range (100-1,000 bp) using Bioanalyzer results.
43. Prepare and run the **Agilent High Sensitivity DNA Chip** per the manufacturer's instructions.
44. Typical Bioanalyzer results for CUT&RUN sequencing libraries are shown in **Figure 5**.
- Note 1: Confirm that positive control histone PTM antibodies (e.g. H3K4me3) enriched for predominantly mononucleosome fragments (~275 bp peak with nucleosomes + sequencing adapters).
 - Note 2: The typical molarity for a 15 µL purified CUT&RUN sequencing library (100-1,000 bp region) is 100-200 nM.

Section VIII: Illumina Sequencing

45. Based on Bioanalyzer molarity calculations, pool libraries at the desired ratios.
- Note: Only 3-8 million paired-end (PE) reads are needed for good CUT&RUN coverage. For lower abundance targets (e.g. H3K4me3), 3-5 million reads are adequate. For higher abundance targets (e.g. H3K27me3), aim for 5-8 million reads.
46. To obtain 3-8M PE reads for 8 samples, we typically load 0.8 pM (500 µL) into a cartridge for a MiniSeq High Output Reagent Kit, 150-cycles (catalog # FC-420-1002).
- Note: Flow cells using the MiniSeq High Output Kit typically yield 25-40 million pass-filter clusters (i.e. 25-40 million PE reads).
47. To obtain 3-8M PE reads for >48 samples, we typically load 0.8 pM (1,500 µL) into a cartridge for a NextSeq 500/550 High Output Kit v2.5, 150-cycles (catalog # 20024907).
- Note: Flow cells using the NextSeq High Output Kit v 2.5 typically yield 400 million pass-

filter clusters (*i.e.* 400 million PE reads).

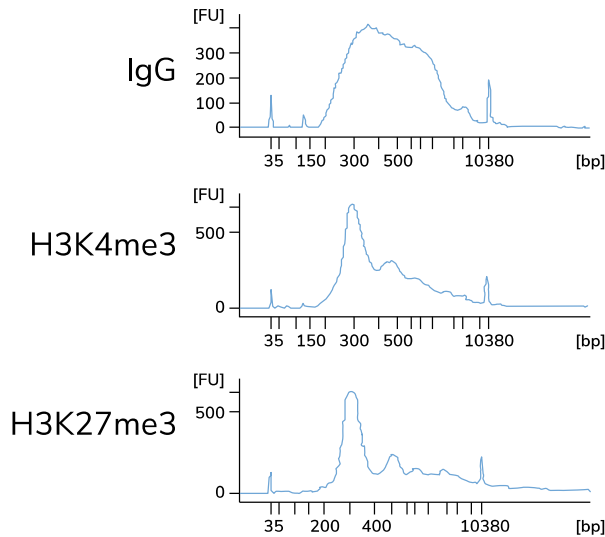


Figure 5: Typical Bioanalyzer traces for IgG negative control and H3K4me3 and H3K27me3 positive control CUT&RUN sequencing libraries (5 ng PCR amplified libraries loaded on Bioanalyzer). H3K4me3 and H3K27me3 libraries are predominantly enriched for mononucleosomes as indicated by the peak at 275 bp (~150 bp mononucleosomes + 125 bp sequence adapters). Quality Bioanalyzer (or TapeStation or other equivalent approach) traces are the best indicator of success prior to sequencing. However, this analysis should not be performed prior to NGS library preparation, since the amount of CUT&RUN DNA is likely to be below the limit of detection for these platforms.

7. Frequently Asked Questions (FAQs)

7.1 General

1. What is the best way to know if a CUT&RUN experiment worked prior to sequencing?

Results from challenging cell inputs/targets may be ambiguous, so EpiCypher recommends including positive and negative controls in every experiment. We have included a series of quality control (QC) checks to analyze sample quality, confirm ConA bead binding, optimize cell permeabilization, assess DNA yields at various stages, and confirm fragment size distribution (see **Appendix II** and **the CUTANA™ CUT&RUN Protocol, Sections V - VII**). If the QC checks and positive and negative controls perform as expected, then proceeding to sequencing with all samples is recommended. If sequencing results for challenging cell inputs/targets are not satisfactory, further optimization may be necessary (e.g. cell type and/or number, digitonin permeabilization, antibody concentration/alternate vendors, etc.).

Recommended experimental design:

- a. Always include control antibodies and control cells
 - i. Start with 500,000 K562 cells
 - ii. Positive control antibody (e.g. H3K4me3 for PTMs; CTCF for TFs)
 - iii. Negative control antibody (IgG)
- b. Experimental sample antibodies and cell types
- c. Use spike-in controls whenever possible (e.g. the SNAP-CUTANA™ K-MetStat Panel can be added into the positive and negative control reactions as well as any sample with an antibody to a methyl-lysine histone PTM target; see **Appendix III**).

Quality control checks before decision to sequence:

- a. Confirm integrity of cells/nuclei before and after permeabilization (**Appendix II**). Samples should not contain clumps and starting counts should be accurate.
- b. Optimize digitonin permeabilization conditions (>95% cells permeabilized).
- c. Confirm sample binding to ConA beads (**Appendix II**).
- d. Ideally, the yield of CUT&RUN enriched DNA for the positive control antibody (e.g. H3K4me3) should be greater than negative control antibody (e.g. IgG). See **Section V**. Note that at low cell inputs, this difference may not be observed, but good quality sequencing data can still be obtained.
- e. Yields from PCR amplified sequencing libraries (~750 ng from 5-10 ng library input). See **Section VI**.
- f. Bioanalyzer or TapeStation traces of sequence-ready libraries for positive control histone PTMs (e.g. H3K4me3) and other PTMs should be enriched for mononucleosome size fragments (~275 bp = ~150 bp nucleosome + 125 bp adapters, see **Section VII**).

- **IMPORTANT:** It is not recommended to assess fragment size distribution of CUT&RUN purified DNA (from **Section V**) prior to library preparation (see further explanation below). This analysis is not indicative of the success of a CUT&RUN experiment, and the amount of purified CUT&RUN DNA is often below the sensitivity of detection for Bioanalyzer/TapeStation approaches. For these reasons, we recommend assessment of CUT&RUN fragment size distribution after library preparation.
- Examples of library distribution in our protocol (**Figure 5**) are based on CUT&RUN analysis of histone PTMs. However, EpiCypher routinely maps TFs using CUT&RUN, and generally recovers 100-120 bp DNA fragments, similar to PTMs. These data provide reliable motif analysis and expected peak structures.
- Nevertheless, some DNA-bound TFs may have smaller size distributions due to processive MNase cleavage, which can generate sub-nucleosomal (<120 bp) DNA fragments. Our DNA purification kit is designed to recover fragments as small as 50 bp, and PCR conditions can also be adjusted to enrich for smaller library fragments if desired (see **FAQs** below).

2. Can I use Bioanalyzer or TapeStation traces to evaluate the success of CUT&RUN prior to library preparation?

We do not advise assessing the fragment size distribution of raw CUT&RUN DNA (from **Section V**) before library preparation. This habit is derived from ChIP-seq methods, which requires careful optimization of chromatin shearing to achieve mononucleosome-sized fragments in Input chromatin. Fragment size distribution in Input chromatin and immunoprecipitated (IP'd) DNA is monitored by agarose gel or capillary electrophoresis (e.g. Bioanalyzer) to confirm assay success at multiple steps.

However, CUT&RUN and ChIP-seq are not the same method, and thus do not entail the same quality control steps. ChIP-DNA yields are typically high due to high cell inputs and elevated background, which are inherent to the IP process. CUT&RUN, on the other hand, is a ChIP-less *in situ* technique that bypasses bulk chromatin fragmentation steps. CUT&RUN also has low background and requires ~10-fold fewer cells vs. ChIP-seq. As a result, raw CUT&RUN DNA yields are below the limit of sensitivity for fragment size distribution using the Bioanalyzer or TapeStation.

The best indicator of CUT&RUN success at this step is that raw DNA yields (e.g. from Qubit) from positive control reaction(s) are greater than yields from the IgG negative control. For low abundance targets, such as H3K4me3, this difference may be slight, while high abundance targets (e.g. H3K27me3) will display a more pronounced difference.

3. Can I use qPCR to evaluate the success of a CUT&RUN experiment?

We do not recommend qPCR for validation of CUT&RUN enrichment. To understand why, it is important to consider why qPCR is a go-to method for ChIP-seq, and how ChIP-seq approaches differ from CUT&RUN.

ChIP involves antibody-mediated enrichment of targets (histone PTMs or chromatin-associated proteins) from bulk fragmented chromatin, or “Input.” qPCR is traditionally used to verify the enrichment of a known on-target region following a ChIP experiment as a readout for experimental success. To determine enrichment, qPCR for the same region must also be performed on the bulk chromatin Input. Regions not enriched by the ChIP reaction are also included as negative controls for comparison.

Not only are these steps unnecessary for CUT&RUN assays, they are also incredibly challenging. CUT&RUN is an *in situ* assay, meaning it is performed in intact cells. Antibody-bound chromatin is cleaved by pAG-MNase, diffuses outside the cell, and is collected for sequencing. There is no immunoprecipitation step and no bulk chromatin Input for enrichment comparisons. In fact, due to the processive nature of MNase, the CUT&RUN fragments may be 50-100 bp (for some TFs), challenging primer design and qPCR analysis.

Instead, we recommend analysis of post-library prepped, purified CUT&RUN sequencing libraries (using the Agilent Bioanalyzer or TapeStation) to confirm DNA size distribution and molarity (see **Section VII**) and proceed to sequencing without qPCR. As you will need only 3-8 million reads per sample, you can multiplex samples, save on sequencing costs, and obtain the genomic data regarding your target. Due to this cost savings, it is much more effective to obtain full sequencing data rather than spending time and money on attempted qPCR optimization.

7.2 Spike-in Controls

4. Can residual *E. coli* in the pAG-MNase prep be used for sample input normalization? What spike-in DNA control does EpiCypher recommend?

Carry-over *E. coli* DNA is present at very low levels in EpiCypher’s pAG-MNase preps. However, at a typical sequencing depth of 3-8 million reads, too few *E. coli* DNA fragments (~hundreds) are recovered for reliably computing sample normalization. Thus, EpiCypher now offers *E. coli* Spike-in DNA to provide sufficient read depth for experimental normalization (**Appendix IV**).

5. Does EpiCypher offer spike-in nucleosome controls for CUT&RUN?

EpiCypher recently launched SNAP Spike-in Controls for CUTANA™ CUT&RUN assays, and now offers the SNAP-CUTANA™ K-MetStat Panel for CUT&RUN reactions against histone lysine methylation targets (see **Appendix III**). The panel comprises highly pure, modified semi-synthetic/recombinant nucleosomes wrapped with DNA containing a PTM-

specific barcode, allowing detection in next-generation sequencing (NGS). SNAP-CUTANA Spike-ins are the ideal physiological control because they replicate the natural substrate of histone PTM antibodies in CUT&RUN, and thus can provide a direct readout on assay success. Our SNAP-CUTANA Spike-ins offer multiple advantages, including sample normalization and in-assay antibody validation. SNAP-CUTANA Spike-ins also provide essential information on the quality of sample inputs, pAG-MNase activity, DNA purification, and library preparation, making them an essential tool for developing, optimizing, and troubleshooting CUT&RUN.

7.3 Sample Input Compatibility

6. What types of cell inputs are compatible with CUT&RUN?

EpiCypher has confirmed that CUT&RUN is compatible with whole cells and nuclei derived from mammalian suspension and adherent cell lines. Protocols to isolate nuclei for CUT&RUN are provided in **Appendix I**. A number of groups have successfully performed CUT&RUN on human and mouse primary tissue⁵⁻⁸, FACS isolated cells⁹, and immune cells^{10,11}. Below are general recommendations for specific sample types:

Adherent cells: Adherent cells present a special challenge for CUT&RUN, as the process must be strong enough to detach cells from culture plates and disaggregate cell clumps, yet gentle enough to preserve cell and nuclear membranes for binding to ConA beads. A major concern is that strong or prolonged enzymatic treatment (e.g. trypsin) can degrade glycoproteins, thereby impairing cell adsorption to the ConA beads.

EpiCypher has compared different methods of collecting adherent cells for CUT&RUN, including cell scraping, trypsin, and accutase. We analyzed the impact of each method on the cell adsorption rate onto ConA beads using MCF-7 cells. Based on these optimization experiments, we recommend a very mild trypsin treatment (0.05% trypsin at 37°C), for the minimal incubation time needed to detach cells as optimized for cell type. Collect cells by centrifugation for ~3 min at 600 x g at RT. Discard supernatant, and then proceed directly to cell counting and CUT&RUN wash steps as outlined in **Section II**. Trypsin is washed away by subsequent washes that are a standard part of the CUT&RUN protocol. This method detaches and monodisperses cells, resulting in >95% adsorption onto ConA beads.

Tissue samples: While EpiCypher does not have tissue-specific protocols available at this time, the primary requirement is that tissue is processed to a monodispersed cell suspension (typically by mechanical maceration or douncing). Enzymatic digestion (e.g. collagenase/dispase) can be used for connective tissue and trypsin (as described above; monitoring dissolution to single cells) may be used for macro-dissected tissue.

Working with immune cells: Lectins (e.g. ConA) play a known role in the innate immune system, and immune cell types may be inadvertently stimulated via binding to ConA beads.

To circumvent this potential problem in CUT&RUN, EpiCypher recommends using nuclei⁹ (see **Appendix I**) or a cross-linking strategy¹² (see detailed cross-linking protocol at epicypher.com/protocols).

7. Is CUT&RUN compatible with frozen or cross-linked cell preparations?

Yes. General guidelines are noted below; our detailed **CUTANA CUT&RUN Cross-linking Protocol** is provided at epicypher.com/protocols, and additional protocols are available upon request (email info@epicypher.com):

Cryopreservation: EpiCypher has confirmed that freeze/thawed cells (e.g. slow freeze in 10% DMSO/media, and then quick thaw at 37°C) generate data of indistinguishable quality to fresh material. This cryopreservation method is optimized to preserve native physiological interactions, minimize cell lysis, and reduce background signal in CUT&RUN. We use similar methods for freeze/thawing isolated nuclei for CUT&RUN; see **Appendix I**.

Cross-linking: It is recommended to first try native samples in CUT&RUN, since this works well for most targets. Of note, EpiCypher has tested previously reported CUT&RUN cross-linking conditions and recommended wash buffers¹². Although yields are significantly lower than from native cells, the resulting data tracks display similar quality (*i.e.* signal-to-noise). Furthermore, for labile targets or highly transient chromatin binding proteins, CUT&RUN signal may be improved by light cross-linking.

When using native CUT&RUN, histone deacetylase activity may contribute to incomplete or low-resolution genomic profiles for certain targets. For example, EpiCypher has observed that certain acetyl PTMs such as H3K27ac and H3K18ac may show enhanced signal after light (e.g. 0.1% formaldehyde, 1 min) to moderate (e.g. 1% formaldehyde, 1 min) cross-linking even though total yields are reduced. However, heavy cross-linking such as that typically used for ChIP-seq (e.g. 1% formaldehyde, 10 min) has been shown to damage histone acetylation signal from K562 cells in CUT&RUN. Therefore, optimal cross-linking conditions for profiling histone acetylation and histone acetylation binding proteins (e.g. some bromodomain-containing proteins) should be empirically determined in the model system of interest.

Importantly, not all acetyl-PTM and readers require cross-linking for high quality data (e.g. H3K9ac, BRD4). Therefore, cross-linking should only be used as a last resort to improve signal-to-noise, since (1) cross-linking reduces yield; (2) cross-linking can decrease specificity and increase artifacts; and (3) key protocol adaptations are required when applying CUT&RUN to cross-linked samples (supplementing Wash, Digitonin, and Antibody Buffers with Triton X-100 and SDS detergents). EpiCypher continues to optimize cross-linking protocols and determine cases where this approach may improve signal. Our current **CUTANA CUT&RUN Cross-linking Protocol** is available at epicypher.com/protocols; similar protocols can be found in the literature¹².

7.4 Antibodies and Targets

8. Does EpiCypher's CUT&RUN protocol work on non-PTM targets?

Yes. The current protocol has been used at EpiCypher to generate CUT&RUN data for numerous non-PTM targets, including CTCF, FOXA1/HNF3A, BRD4, and SMARCA4 (BRG1). Antibodies to chromatin-associated protein targets validated for use in CUT&RUN can be found at epicypher.com/cut-and-run-compatible-antibodies; for targets not yet on this list, contact info@epicypher.com for recommendations. No protocol modifications were necessary to generate these data since the columns used to purify the CUT&RUN-enriched chromatin fragments retain >50 bp DNA, allowing for NGS library distributions down to ~175 bp (~50bp DNA + 125 bp sequencing adapters).

For TFs, which may generate sub-nucleosomal size (<120 bp) DNA fragments, modifications to the library protocol have been reported^{7,13} to improve the representation of these smaller fragments during library preparation. Briefly, to enrich for CUT&RUN library fragments <120 bp, the authors altered the MNase inactivation step (after end repair) from 65°C for 30 min to 50°C for 1 hr. Also, after the ligation reaction, AMPure bead size-selection ratio was increased from 1.1x volume to 1.75x volume (retains library fragments >100 bp). For TF CUT&RUN fragment sizes, both nucleosomal (~150 bp) and sub-nucleosomal (<120 bp) reflect TF chromatin occupancy locations. However, the sub-nucleosomal fragments provide higher resolution mapping.

9. What antibodies does EpiCypher recommend for CUT&RUN? Will reliable ChIP antibodies work for CUT&RUN?

Through our extensive development of CUT&RUN assays to various targets, EpiCypher has found that antibodies that work well in ChIP may not always work in CUT&RUN. To address this unmet need for researchers, EpiCypher has begun screening antibodies for high quality performance in CUT&RUN and its sister technology CUT&Tag. Below we outline our criteria for antibodies to chromatin-associated proteins and histone PTMs.

Chromatin-associated protein targets: EpiCypher has recently launched **CUTANA™ Compatible Antibodies** for chromatin associated protein targets (see the growing list of antibodies: epicypher.com/cut-and-run-compatible-antibodies). Every lot of a CUTANA Compatible Antibody is validated directly in CUT&RUN and/or CUT&Tag mapping assays. To pass testing, the antibody must generate CUT&RUN (and/or CUT&Tag) genomic distribution profiles consistent with the reported function of the target protein (for example, DNA binding motif analysis for TFs). Our scientists are continuously testing antibodies to novel targets. For targets of interest not on this list, contact us for antibody recommendations: techsupport@epicypher.com.

Histone PTM targets: Antibodies to histone PTMs are particularly susceptible to off-target cross-reactivity which can compromise biological interpretations⁴. EpiCypher has conducted extensive studies of histone PTM antibodies in ChIP-seq by using our exclusive SNAP-ChIP® nucleosome spike-in control technology. Although the majority of antibodies exhibit major problems such as poor specificity and/or low efficiency, best-in-class reagents for ChIP-seq do exist, and have been identified using SNAP-ChIP Spike-ins (see ref ⁴ and chromatinantibodies.com *Maryanski et al., In preparation*). As noted above, EpiCypher has recently launched the first SNAP Spike-in Controls for CUTANA CUT&RUN assays (*i.e.* the SNAP-CUTANA K-MetStat Panel). We are currently using these novel spike-ins for histone PTM antibody validation in CUT&RUN and CUT&Tag, and plan to launch the first set of validated histone lysine methylation antibodies in the coming months.

For more information or for antibody recommendations, please contact techsupport@epicypher.com.

8. References

- 1 Skene, P. J. & Henikoff, S. An efficient targeted nuclease strategy for high-resolution mapping of DNA binding sites. *eLife* **6**, doi:10.7554/eLife.21856 (2017).
- 2 Schmid, M., Durussel, T. & Laemmli, U. K. ChIC and ChEC; genomic mapping of chromatin proteins. *Molecular cell* **16**, 147-157, doi:10.1016/j.molcel.2004.09.007 (2004).
- 3 Skene, P. J., Henikoff, J. G. & Henikoff, S. Targeted in situ genome-wide profiling with high efficiency for low cell numbers. *Nature protocols* **13**, 1006-1019, doi:10.1038/nprot.2018.015 (2018).
- 4 Shah, R. N. *et al.* Examining the Roles of H3K4 Methylation States with Systematically Characterized Antibodies. *Molecular cell* **72**, 162-177.e167, doi:10.1016/j.molcel.2018.08.015 (2018).
- 5 de Bock, C. E. *et al.* HOXA9 Cooperates with Activated JAK/STAT Signaling to Drive Leukemia Development. *Cancer Discov* **8**, 616-631, doi:10.1158/2159-8290.CD-17-0583 (2018).
- 6 Janssens, D. H. *et al.* Automated in situ chromatin profiling efficiently resolves cell types and gene regulatory programs. *Epigenetics & chromatin* **11**, 74, doi:10.1186/s13072-018-0243-8 (2018).
- 7 Liu, N. *et al.* Direct Promoter Repression by BCL11A Controls the Fetal to Adult Hemoglobin Switch. *Cell* **173**, 430-442.e417, doi:10.1016/j.cell.2018.03.016 (2018).
- 8 Uyehara, C. M. & McKay, D. J. Direct and widespread role for the nuclear receptor EcR in mediating the response to ecdysone in *Drosophila*. *Proceedings of the National Academy of Sciences of the United States of America* **116**, 9893-9902, doi:10.1073/pnas.1900343116 (2019).
- 9 Hainer, S. J., Boskovic, A., McCannell, K. N., Rando, O. J. & Fazio, T. G. Profiling of Pluripotency Factors in Single Cells and Early Embryos. *Cell* **177**, 1319-1329.e1311, doi:10.1016/j.cell.2019.03.014 (2019).
- 10 Mathsyaraja, H. *et al.* Max deletion destabilizes MYC protein and abrogates Emicro-Myc lymphomagenesis. *Genes & development*, doi:10.1101/gad.325878.119 (2019).
- 11 Roth, T. L. *et al.* Reprogramming human T cell function and specificity with non-viral genome targeting. *Nature* **559**, 405-409, doi:10.1038/s41586-018-0326-5 (2018).
- 12 Zheng, X. Y. & Gehring, M. Low-input chromatin profiling in *Arabidopsis* endosperm using CUT&RUN. *Plant reproduction* **32**, 63-75, doi:10.1007/s00497-018-00358-1 (2019).
- 13 Zhu, Q., Liu, N., Orkin, S. H. & Yuan, G. C. CUT&RUNTools: a flexible pipeline for CUT&RUN processing and footprint analysis. *Genome Biol* **20**, 192, doi:10.1186/s13059-019-1802-4 (2019).

Appendix I: Nuclei Isolation & Cryopreservation Protocol for CUT&RUN

Description: This protocol gives instructions on fresh nuclei isolation, cryopreservation of nuclei, and thawing nuclei – all of which are compatible with CUT&RUN assays. Note that this protocol is designed similarly to the CUTANA CUT&RUN protocol, *i.e.* to batch process cells for multiple CUT&RUN reactions, and was developed using 500,000 K562 cells per CUT&RUN reaction. Adjust volumes and cell numbers for your experiments as needed.

Reagents & Buffers Needed

Item	Vendor	Catalog No.
cOmplete™, Mini, EDTA-free Protease Inhibitor Cocktail Tablets	Roche	11836170001
Spermidine*	Sigma Millipore	S2501
0.4% Trypan blue	Any	
1X Phosphate Buffered Saline (PBS)	Any	

*1M spermidine preparation: Dissolve 1 gram spermidine (MW = 254.63) in 3.93 mL molecular grade H₂O. Store in single-use aliquots at -20°C for 6 months.

Nuclear Extraction (NE) Buffer

NE Buffer base:

20 mM HEPES, pH 7.9

10 mM KCl

0.1% Triton X-100

20% Glycerol

Filter sterilize. Store at 4°C for up to 6 months.

Add to NE Buffer on day of experiment:

1X cOmplete Mini-Tablet (1 tablet)

0.5 mM Spermidine

Keep at 4°C throughout experiment. Discard after 1 day.

Protocol: Nuclei Harvest, from Beginning of Section II of CUT&RUN Protocol

- Harvest 500,000 cells per planned CUT&RUN reaction by spinning for 3 min at 600 x g at RT in 1.5 mL tube. Pipette or aspirate to discard supernatant.
 - Note 1:** For all steps, scale buffer volumes with number of cells, *e.g.* use 1 mL buffer for 5 million cells. Adjust volumes and cell numbers for your experiment as needed.
 - Note 2:** Prepare ~10% excess cells to confirm cell integrity prior to nuclear isolation and nuclei integrity at the end of the protocol (see **Protocol Notes**, below).
- Resuspend cells in 100 µL/reaction RT **1X PBS**.
 - Note:** Set 10 µL aside for confirmation of cell integrity.
- Spin for 3 min at 600 x g at RT. Decant or pipette to remove and discard supernatant.
- Resuspend cell pellet in 100 µL/reaction cold **NE Buffer**.

5. **Incubate** reactions for **10 min on ice**.
6. Spin for 3 min at 600 x g at **4°C**. Pipette or aspirate to discard supernatant.
 - Note: The pellet should change in appearance from a sticky, pale yellow pellet (cells) to a white, fluffy pellet (nuclei).
7. Resuspend nuclei in 100 µL/reaction cold **NE Buffer**.
8. Freeze (see below) or proceed to ConA bead conjugation step (**Section II Step 11**). Nuclei in NE Buffer can be directly added to activated ConA beads.
 - Note: Set 10 µL aside for confirmation of isolated nuclei integrity

Protocol Notes (also see Appendix II)

To examine the efficiency of nuclear isolation and ensure intact nuclei:

1. Remove 10 µL aliquots from washed cells/final nuclei.
2. Combine 10 µL aliquots with 10 µL 0.4% Trypan blue dye, mix, and load onto cell counter or hemacytometer slide.
3. Examine under brightfield or phase microscope (**Figure 6, Appendix II**). Cells should not take up Trypan blue and appear clear/white, whereas nuclei should take up Trypan and appear blue.

Protocol: Cryopreservation and thawing of nuclei

1. To cryopreserve nuclei, slowly freeze aliquots in an isopropanol-filled chiller in a -80°C freezer.
 - Note: If necessary, nuclei can be shipped on dry ice in this state.
2. To avoid nuclear lysis and chromatin fragmentation, thaw nuclei quickly by placing tubes on 37°C block until thawed.
3. Proceed to ConA bead conjugation step, **Section II Step 11**. Nuclei in NE Buffer can be directly added to activated ConA beads.

Protocol: Cryopreservation and thawing of cells

1. To cryopreserve cells, supplement cell culture media with a cryoprotective agent (e.g. 10% DMSO in media) and slowly freeze aliquots in an isopropanol-filled chiller in a -80°C freezer.
2. When ready to perform CUT&RUN, quickly and completely thaw samples at 37°C and then prepare cells as described in **Section II Step 7**.

Appendix II: Sample Integrity & Bead Conjugation Quality Control Checks

Description: This Appendix provides detailed instructions on checking the quality of cells/nuclei before and after permeabilization + conjugation to ConA beads. Of note, these Quality Control (QC) checks should also be used to check the integrity of cells before nuclei isolation, as described in **Appendix I**. All QC checks were developed using batch processed K562 cells and nuclei (*i.e.* for multiple CUT&RUN reactions).

We recommend checking the quality of starting materials (cells or nuclei) prior to every CUT&RUN experiment. For new users or for working with new cell types/sample inputs, we strongly recommend performing the ConA bead conjugation checks. This is a key step in the CUT&RUN protocol, and if cells are not successfully bound and permeabilized, CUT&RUN yields will be dramatically reduced.

Reagents, Materials & Equipment Needed

Item	Vendor
0.4% Trypan blue	Any
Hemocytometer	Any
Cell counting slides	Any
Brightfield/phase contrast microscope or automated cell counter	Any

Protocol: Sample Integrity and ConA Bead Conjugation QC Checks, Starting from Beginning of Section II of CUT&RUN Protocol

- In **Section II Step 7**, process excess cells (or nuclei) to have leftover enough leftover material (*e.g.* prepare ~10% excess volume for batch processing or one extra sample).
- Prior to ConA bead conjugation, take 10 μ L from washed cells (or prepared nuclei) and evaluate sample integrity as follows:
 - Add 10 μ L of 0.4% **Trypan blue**.
 - Mix 10 times by pipetting.
 - Transfer 10 μ L to a counting slide.
 - View under brightfield/phase microscope or cell counter.
 - Confirm sample integrity or troubleshoot as needed (see **Figure 6** and **Table 6**):
 - Cells/nuclei should not be clumped and should show the expected morphology
 - Intact cells will not absorb trypan blue and appear white (**Figure 6A**)
 - Nuclei will take up Trypan and appear blue (**Figure 6B**)
 - Troubleshoot as necessary (**Table 6**)
- Proceed with bead binding (working in batch if possible) by adding 10 μ L activated **ConA Beads** per 100 μ L washed cells/nuclei in a 1.5 mL tube. Again, make sure to prepare ~10% extra volume.
- Gently vortex and/or pipette to mix ConA beads with cells/nuclei.
- Incubate** bead slurry for **10 min at RT**. Cells/nuclei will adsorb to the activated ConA beads.

6. Place 1.5 mL tube(s) on magnet until slurry clears. Transfer 10 μ L supernatant into a fresh 1.5 mL tube (**Unbound fraction**) and set aside for bead binding integrity check. Pipette to remove remaining supernatant and discard.
7. Add 50 μ L/reaction cold **Antibody Buffer** quickly, to avoid bead drying. If working in bulk, scale volume for total number of planned reactions. Gently vortex immediately and thoroughly to an even resuspension.
8. Remove 10 μ L and transfer to a fresh tube (**Bead-Bound Fraction**). Place remaining sample on ice.
9. Perform Trypan blue staining and bead binding integrity check as described in **Steps 2a-e** (above), comparing **Unbound Fraction** and **Bead-Bound Fraction**.
 - Successful binding will show Trypan positive cells/nuclei surrounded by ConA beads with little leftover in the Unbound Fraction; compare **Figures 6C and 6D**. For troubleshooting approaches, see **Table 6**.
9. Continue with the **CUT&RUN Protocol, Section III Steps 15-16** (Antibody Binding).

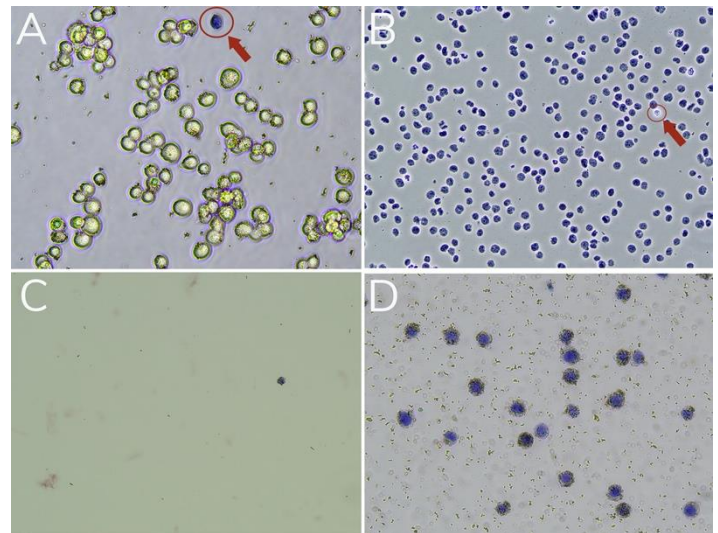


Figure 6: Representative images of cell and nuclei samples for ConA Bead binding. Samples were stained with Trypan blue and visualized under brightfield microscope. **(A) Cells** before bead binding. A dead cell (blue; Trypan positive) is circled in red, whereas the remaining live cells are bright white and round. **(B) Nuclei** before bead binding. An intact cell (Trypan negative) is circled in red, whereas isolated nuclei are Trypan stained (see **Appendix I** for preparation of nuclei). **(C) Unbound Fraction** shows little to no material leftover after ConA Bead binding. **(D) Representative Bead-Bound Fraction** showing nuclei (blue) successfully bound to activated ConA Beads (brown specks).

Samples	Figure	Success Metric (Trypan Blue staining)	Troubleshooting Tips
Cells	Fig. 6A	Cells should be bright (Trypan blue excluded), round, unclumped, and ideally show >90% viability. Over 80% minimum viability is recommended, as excess dead cells increase background in CUT&RUN.	Optimize cell culture conditions; use fresh media, evaluate potential contamination issues.
Nuclei	Fig. 6B	Nuclei should be >95% Trypan blue positive and unclumped.	See Appendix I for a detailed nuclei preparation protocol.
Unbound Fraction	Fig. 6C	Little to no material should be present if binding to beads occurred.	Ensure that ConA Beads were never frozen, cells/nuclei were not clumped, beads did not become clumped or dried out, and all buffers were correctly prepared.
Bead-Bound Fraction	Fig. 6D	Successful ConA bead binding will show Trypan blue positive permeabilized cells/nuclei surrounded by beads.	

Table 6: Quality Control testing of cell/nuclei integrity and ConA bead binding steps.

Appendix III: SNAP-CUTANA™ Spike-in Controls

Description: This Appendix describes EpiCypher’s quantitative nucleosome spike-in technology, or SNAP (Sample Normalization and Antibody Profilng) Spike-in Controls for CUTANA™ CUT&RUN/CUT&Tag assays. In **Section I**, we provide an overview of SNAP-CUTANA™ Spike-ins and how they should be incorporated into CUT&RUN/Tag workflows. **Section II** contains more detailed descriptions and examples of SNAP-CUTANA Spike-in applications, including antibody specificity profiling, flagging failed reactions, and next-generation sequencing (NGS) normalization. In **Section III**, we offer detailed instructions on extracting SNAP-CUTANA Spike-in data from CUT&RUN/CUT&Tag datasets using the shell script and Excel template provided on the EpiCypher product page. **Table 7** contains the DNA barcodes associated with each nucleosome in the K-MetStat Spike-in Panel ([EpiCypher 19-1002](#)).

Section I: An Introduction to SNAP-CUTANA™ Spike-in Controls

What are SNAP-CUTANA Spike-ins?

SNAP-CUTANA Spike-in Controls are panels of semi-synthetic/recombinant nucleosomes carrying defined histone post-translational modifications (PTMs), which can be distinguished in NGS by unique PTM-specific DNA barcodes. Panels are grouped by PTM class, such as that related to histone lysine methylation (K-MetStat Panel, **Figure 7**) and contain widely studied and disease-relevant modifications. This strategy enables a direct readout of assay success, *in situ* antibody validation against closely related PTMs (which have high risk for cross-reactivity) and provides quantitative tools for valuable epigenetic targets.

panels of semi-synthetic/recombinant nucleosomes carrying defined histone post-translational modifications (PTMs), which can be distinguished in NGS by unique PTM-specific DNA barcodes. Panels are grouped by PTM class, such as that related to histone lysine methylation (K-MetStat Panel, **Figure 7**) and contain widely studied and disease-relevant modifications. This strategy enables a direct readout of assay success, *in situ* antibody validation against closely related PTMs (which have high risk for cross-reactivity) and provides quantitative tools for valuable epigenetic targets.

K-MetStat Panel

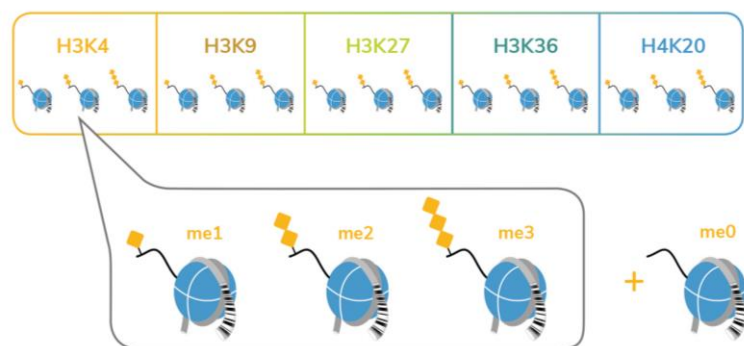


Figure 7: Histone methyl-lysine PTMs in the SNAP-CUTANA™ K-MetStat Panel.

How are the spike-ins detected? What is the “barcode”?

As in **Figure 8**, each nucleosome is wrapped with a DNA template containing the Widom 601 nucleosome positioning sequence (Lowary & Widom, *J. Mol. Biol.* 1998), and a 22 bp barcode unique to each

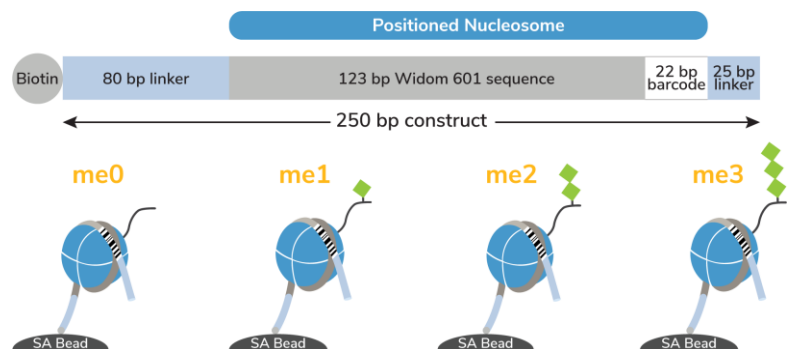


Figure 8: Schematic showing the DNA barcoding and magnetic bead immobilization used in SNAP-CUTANA™ Spike-in Controls.

PTM in the panel. This nucleosome assembly DNA is flanked by linker DNA, providing an appropriate substrate for pAG-MNase cleavage or pAG-Tn5 tagmentation in CUT&RUN or CUT&Tag, respectively. The 5' linker DNA also contains a biotin tag (**Figure 8**), allowing spike-ins to be immobilized to streptavidin (SA) magnetic beads for streamlined incorporation in CUT&RUN/CUT&Tag workflows alongside cells (or nuclei) immobilized onto ConA magnetic beads.

One important feature of SNAP-CUTANA Spike-in Panels is that each histone PTM is present in “duplicate,” allowing scientists to monitor technical variation of spike-in recovery. This means that for a given panel we prepare each modified nucleosome using two unique DNA-barcoded templates. As an example, the **SNAP-CUTANA K-MetStat Spike-in panel** ([EpiCypher 19-1002](#)) contains 16 PTM-defined states (15 methyl-lysines: me1, me2, me3 at H3K4, H3K9, H3K27, H3K36 and H4K20, and an unmodified control). However, each modified histone octamer was wrapped with two different barcoded DNAs, so the final panel contains 32 distinct DNA-barcoded nucleosomes to be monitored by NGS (**Table 7**).

SNAP-CUTANA K-MetStat Spike-in Panel : DNA Barcodes		
PTM	Barcode A (Nuc Replicate 1)	Barcode B (Nuc) Replicate 2
Unmodified	TTCGCGCGTAACGACGTACCGT	CGCGATACGACCGCGTTACGCG
H3K4me1	CGACGTTAACGCGTTTCGTACG	CGCGACTATCGCGCGTAACGCG
H3K4me2	CCGTACGTCGTGTCGAACGACG	CGATACGCGTTGGTACGCGTAA
H3K4me3	TAGTTCGCGACACCGTTCGTTCG	TCGACGCGTAAACGGTACGTCG
H3K9me1	TTATCGCGTCGCGACGGACGTA	CGATCGTACGATAGCGTACCGA
H3K9me2	CGCATATCGCGTCGTACGACCG	ACGTTTCGACCGCGGTTCGTACGA
H3K9me3	ACGATTCGACGATCGTCGACGA	CGATAGTCGCGTCGCACGATCG
H3K27me1	CGCCGATTACGTGTCGCGCGTA	ATCGTACCGCGCGTATCGGTTCG
H3K27me2	CGTTCGAACGTTTCGTTCGACGAT	TCGCGATTACGATGTTCGCGCGA
H3K27me3	ACGCGAATCGTCGACGCGTATA	CGCGATATCACTCGACGCGATA
H3K36me1	CGCGAAATTCGTATACGCGTTCG	CGCGATCGGTATCGGTACGCGC
H3K36me2	GTGATATCGCGTTAACGTTCGCG	TATCGCGCGAAACGACCGTTTCG
H3K36me3	CCGCGCGTAATGCGCGACGTTA	CCGCGATACGACTCGTTCGTTCG
H4K20me1	GTCGCGAACTATCGTCGATTTCG	CCGCGCGTATAGTCCGAGCGTA
H4K20me2	CGATACGCCGATCGATCGTTCGG	CCGCGCGATAAGACGCGTAACG
H4K20me3	CGATTCGACGGTCGCGACCGTA	TTTCGACGCGTCGATTTCGGCGA

Table 7: SNAP-CUTANA K-MetStat Spike-in DNA barcode sequences.

Importantly, the DNA barcodes do not overlap with known sequences in the human, mouse, fly, or yeast genome and can therefore be distinguished from sample chromatin. We provide instructions on DNA barcode detection in this Appendix (see **Section III**).

How are they “spiked” into assays?

SNAP-CUTANA nucleosomes are supplied pre-immobilized to magnetic SA beads via 5' biotinylated linker DNA (**Figure 8**), allowing them to be processed alongside ConA bead-conjugated cells in CUT&RUN and CUT&Tag reactions. SNAP-CUTANA Spike-in controls are added to CUT&RUN and CUT&Tag reactions just prior to antibody addition, in one simple pipetting step (**Figure 9**). The antibody will bind its target epitope in sample cells and the spike-in panel. Addition and activation of pAG-MNase (in CUT&RUN) or pAG-Tn5 (in CUT&Tag) will cleave or tagment the associated spike-in nucleosome, respectively. In both cases, the targeted standard will be processed with sample chromatin through library preparation, PCR amplification, and NGS (**Figure 9**).

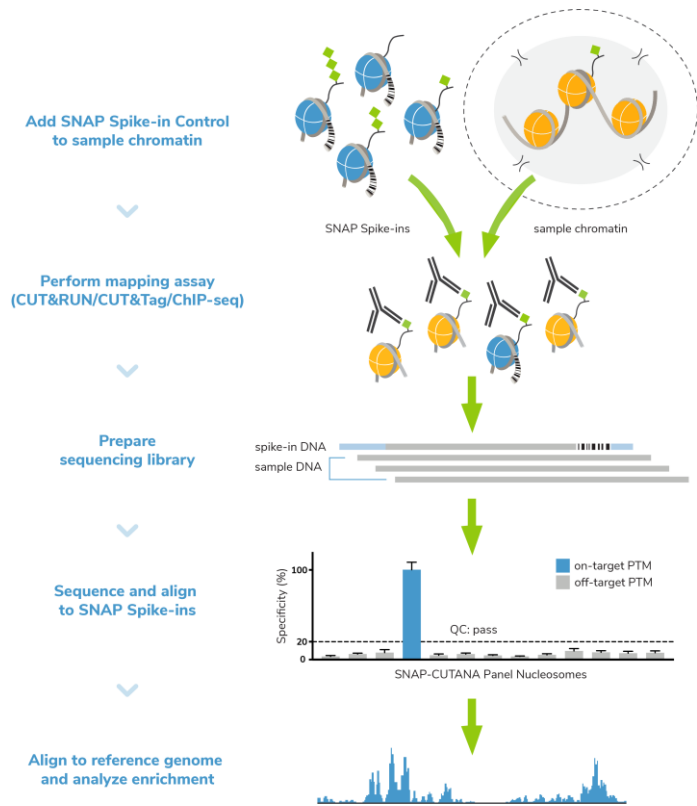


Figure 9: Schematic showing addition of SNAP-CUTANA™ Spike-in Controls during CUT&RUN workflow.

How can the SNAP-CUTANA Spike-in data be used? Why should I add these into my CUT&RUN or CUT&Tag workflow?

SNAP-CUTANA Spike-ins are the only fully defined, quantitative chromatin mapping assay control that represents the nucleosome target of histone-PTM antibodies. Thus, relative recovery of SNAP-CUTANA Spike-in Controls serve as a straightforward proxy for assay success. Because the spike-ins contain a panel of related PTM states (**Figure 7**), assessing the recovery of each nucleosome can also be used as a direct readout of antibody specificity. This is crucial, as many histone PTM antibodies display cross-reactivity, lot variation, and/or assay-specific performance (e.g. ChIP-seq vs. CUT&RUN/Tag). Finally, because the spike-in nucleosomes are carried through the protocol alongside sample chromatin, SNAP-CUTANA Spike-ins can be used to normalize NGS data and enable reliable cross-sample comparisons. See **Figure 10** for a general outline of how SNAP-CUTANA Spike-ins can be used to examine the success of CUT&RUN and CUT&Tag workflows; for more detail on each of these steps, see **Section II**.

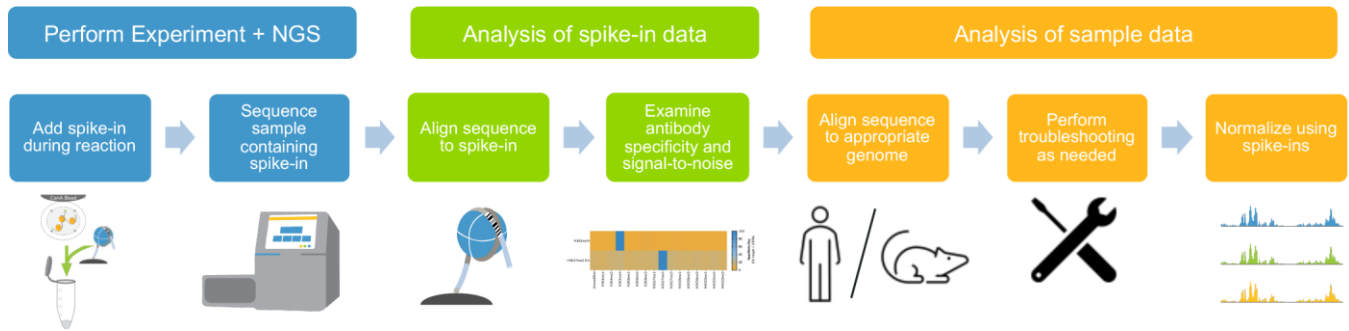


Figure 10: SNAP-CUTANA™ Spike-ins offer a quantitative control for CUT&RUN/CUT&Tag assay workflows. These defined nucleosome spike-ins can be used to determine the success of your assay, including on-target antibody specificity, as well as assay normalization and experimental comparisons.

Section II: How to use SNAP-CUTANA™ Spike-ins for CUT&RUN and CUT&Tag Workflows

1. Combine SNAP-CUTANA Spike-ins with positive and negative control antibodies for CUT&RUN/CUT&Tag assay optimization. When first starting CUT&RUN/CUT&Tag, or when trying these approaches under new experimental conditions (new cell types, changing cell number inputs, fixed vs. native preps, etc.) we strongly recommend using the SNAP-CUTANA K-MetStat spike-ins with positive (e.g. H3K4me3, [EpiCypher 13-0041](#)) and negative (e.g. IgG, [EpiCypher 13-0042](#)) control antibodies to validate your CUT&RUN workflow. These antibodies are lot-validated for superior performance and reliability in CUT&RUN/Tag. In fact, EpiCypher includes these exact positive and negative control antibodies and SNAP-CUTANA K-MetStat Spike-ins in every experiment, and have found them to be essential tools for initial workflow optimization and daily performance monitoring.

As an example, we used CUT&RUN to map H3K27me3 in K562 cells, with IgG and H3K4me3 antibodies as controls, and SNAP-CUTANA K-MetStat Spike-ins added to each reaction (**Figure 11**). The first confirmation of optimized experimental conditions was relative recovery of each spike-in DNA barcode in NGS data (see **Section III** for detailed analysis instructions). Here the IgG control showed no target preference (as expected), while the H3K4me3 control and H3K27me3 antibodies each enriched their target of interest with very low background (**Figure 11A**). This provides confidence that multiple technical aspects of the CUT&RUN experiment (*i.e.* antibodies and pAG-MNase cleavage conditions) were successful, supporting further analysis of the sample. NGS data for the sample chromatin showed the expected PTM enrichment patterns (e.g. H3K4me3 showed tight peaks at transcription start sites [TSS] with minimal IgG background; **Figure 11B-C**). These robust controls allowed us to have high confidence in the accuracy of our H3K27me3 sequencing results.

Note: If the spike-in analysis reveals off-target cleavage or low signal-to-noise this indicates problems with the assay workflow. Results from the spike-ins can guide troubleshooting experiments, as outlined in **Table 8**, below.

Use SNAP-CUTANA™ Spike-ins and control antibodies to develop new CUT&RUN assays

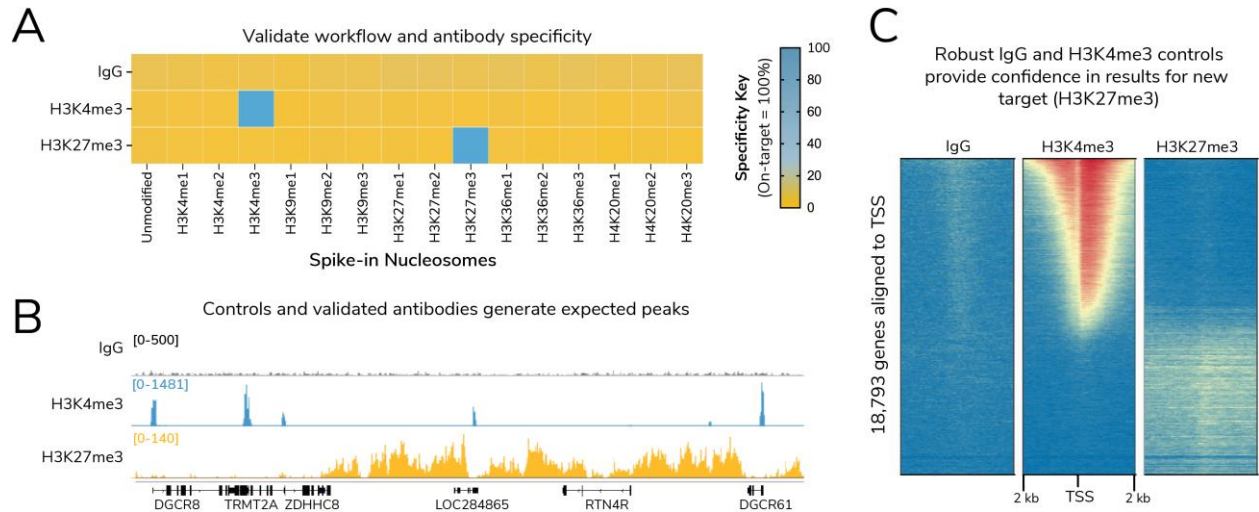


Figure 11: SNAP-CUTANA™ Spike-in Controls and control antibodies were used to validate CUT&RUN workflows for K562 cells. (A) PTM antibodies recovered on-target spike-in, while IgG antibody displayed no preference. Heatmap data normalized to DNA barcodes from on-target PTM; IgG normalized to total reads. **(B,C)** CUT&RUN generated expected H3K4me3 and H3K27me3 maps profiles from K562 sample chromatin. **(B)** RPKM normalized peaks on representative regions. **(C)** Heatmaps display signal intensity +/- 2 kb from transcription start site (TSS). Genes in each heatmap are ordered by signal intensity from H3K4me3 profiles.

2. Use SNAP-CUTANA Spike-ins to find fit-for-purpose antibodies to lysine methylation targets. Test your favorite antibody or inquire at info@epicypher.com for recommendations (**Figure 12**). EpiCypher performs extensive (and ongoing) antibody testing in epigenomic applications (to date: >1,000 commercial reagents) and has determined that the vast majority (>70%!!) of histone PTM antibodies are non-specific and/or display poor enrichment¹. From these studies we have determined that:

- Antibody capability on modified histone peptide arrays does not transfer to epigenomic assays that consider full nucleosome structure (e.g. CUT&RUN, CUT&Tag, or ChIP-seq).
- Validation of an antibody in one chromatin mapping assay does not guarantee success in another (e.g. ChIP validation does not transfer to CUT&RUN or CUT&Tag, see **FAQs**).
- Specific PTM recognition under one set of conditions does not guarantee specific PTM recognition in all conditions. Antibodies must always be validated for user-specific conditions with spike-in controls.

Thus, application-specific (*i.e. in situ*) testing with PTM-defined, target-representative controls is essential to success. As minimum criteria, EpiCypher recommends selecting antibodies that show <20% cross-reactivity to each off-target PTM in the K-MetStat panel. Below, we address some common questions when using nucleosome spike-ins for antibody specificity testing:

Question: What does antibody cross-reactivity look like?

EpiCypher recommends a cutoff of <20% binding to each off-target nucleosome in the K-MetStat panel. This is determined as described in **Section III** of this Appendix, using SNAP-CUTANA Spike-in data with **on-target** PTM barcode recovery set to 100%. Cross-reactivity <20% supports the accuracy of biological findings and interpretations¹. As an example of how antibody cross-reactivity can affect NGS data, we show specificity profiles for one H3K4me3 and three H3K27me2 antibodies (CUT&RUN with SNAP-CUTANA K-MetStat panel in K562 cells: **Figure 13**). Read counts for each DNA-barcoded nucleosome were normalized to the stated PTM target and used to generate the heatmaps in **Figure 13A**. Note the anti-H3K4me3 and anti-H3K27me2 showing high specificity for their respective PTM target, relative to two H3K27me2 antibodies that displayed substantial cross-reactivity to H3K4me3 spike-in nucleosome. This cross-reactivity identified by SNAP-CUTANA Spike-ins was reflected by CUT&RUN sequencing from the cell sample (**Figure 13B**, discussed in detail below). Without the SNAP-CUTANA Spike-in Controls to flag antibody cross-reactivity, such data could misinform biological conclusions.

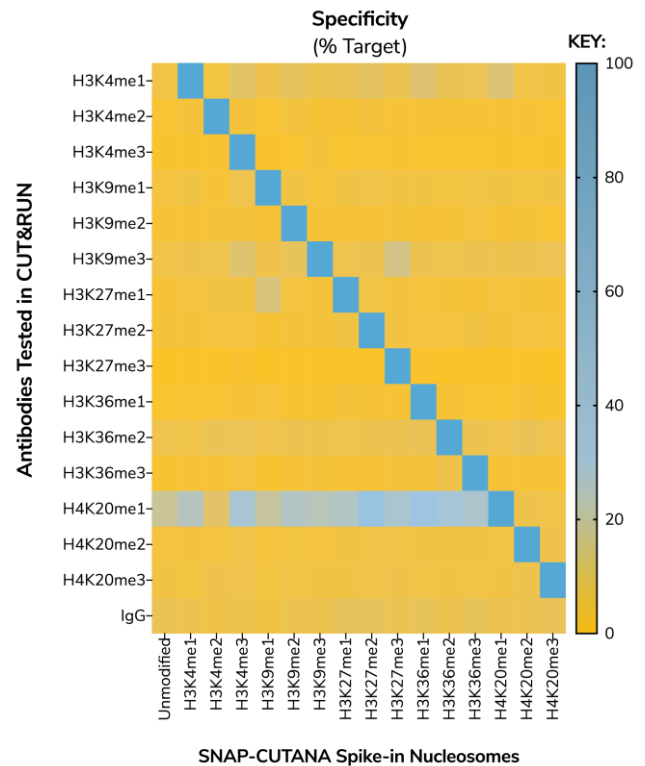


Figure 12: EpiCypher’s best-to-date CUT&RUN antibodies to lysine methylation PTMs. Each row displays SNAP-CUTANA™ K-MetStat Spike-in data for a PTM antibody validated in CUT&RUN. Heatmap data for each antibody are normalized to DNA barcodes from the on-target PTM.

Question: Is antibody cross-reactivity from spike-ins an accurate representation of antibody cross-reactivity in my sample chromatin?

Figure 13B shows heatmaps of CUT&RUN data from K562 cells aligned to gene transcription start sites (TSS, +/- 2 kb) using antibodies characterized with SNAP-CUTANA K-MetStat Spike-in Controls (**Figure 13A**). Note the “anti-correlated” enrichment pattern from highly specific H3K4me3 and H3K27me2 antibodies (**Figure 13B**). The red boxes highlight a contaminating, TSS-enriched (*i.e.* H3K4me3-like) signal using the H3K27me2 antibodies that cross-react with H3K4me3. These results demonstrate that SNAP-CUTANA Spike-ins correctly predicted antibody specificity and performance in CUT&RUN cell sample data, in agreement with our previous studies using SNAP Spike-ins for ChIP-seq¹. Thus, including SNAP-CUTANA Spike-ins allows researchers to directly assess the accuracy of NGS data, providing a powerful tool for advanced epigenomics research.

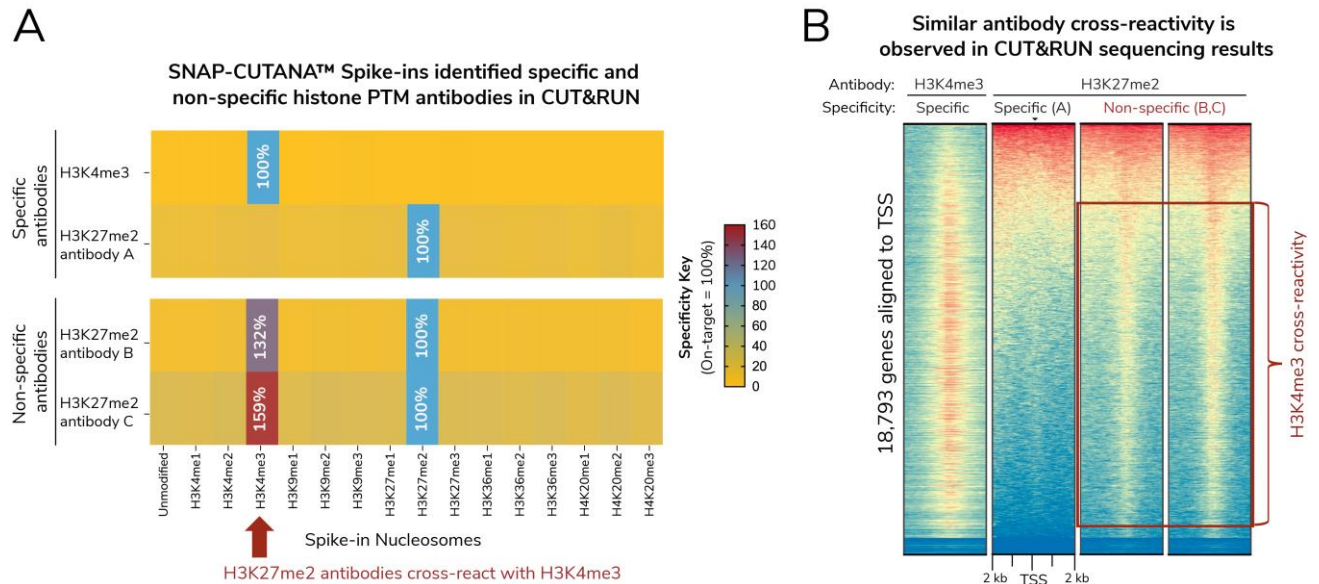


Figure 13: Antibody cross-reactivity identified by SNAP-CUTANA Spike-ins (A) is reflected in CUT&RUN sequencing results (B). (A) SNAP-CUTANA K-MetStat spike-ins were used to test H3K4me3 and H3K27me2 antibodies in CUT&RUN workflows. Data are shown normalized to on-target PTM. (B) CUT&RUN data from K562 cells was generated using the antibodies characterized in (A). Heatmaps display signal intensity +/- 2 kb from TSS. Gene rows in each heatmap are linked and ordered by signal intensity sorted by H3K4me3 specific antibody profile.

- 3. Use SNAP-CUTANA Spike-ins to monitor experimental success in every possible reaction.** While EpiCypher CUTANA protocols are extensively optimized and robust, CUT&RUN/CUT&Tag reactions can occasionally go awry. Including a SNAP-CUTANA Spike-in Control Panel in every reaction identifies problematic experiments and guides troubleshooting. For now, we **minimally** suggest adding SNAP-CUTANA K-MetStat Spike-ins to control reactions containing the positive (e.g. H3K4me3) and negative (IgG) control antibodies. We strongly recommend including these controls in every experiment.

However, the K-MetStat panel has obvious application to every covered lysine methylation target (me0-1-2-3 at H3K4, H3K9, H3K27, H3K36 and H4K20). Similarly comprehensive panels or targeted reagents are in development for histone lysine acylation (K-AcylStat), ubiquitylation (K-UbStat), arginine methylation (R-MetStat), histone oncomutations, histone variants, and even chromatin associated proteins ([inquire](#) for specific progress on each). Here we will discuss how SNAP-CUTANA Spike-ins can be used to identify aberrant CUT&RUN/Tag reactions. See **Table 8** for specific troubleshooting approaches.

Question: What do successful assays look like using SNAP-CUTANA Spike-ins?

In a successful assay to a given histone PTM, spike-ins reveal specific recovery of the target PTM with minimal background and/or cross-reactivity, and genomic enrichment patterns

appear as expected. This is shown in **Figure 11** (explained above), and in **Figure 14** (below). In **Figure 14**, we used SNAP-CUTANA K-MetStat Spike-ins in CUT&RUN to map H3K4me3 and H3K27me3 from four independently prepared mouse primary B cell samples (10K cells each; protocol optimization pilot experiments with a multi-lab consortium). **Figure 14A** shows NGS results from the spike-in nucleosomes for each CUT&RUN reaction using the same antibodies. Samples 1-3 displayed expected antibody performance, with low cross-reactivity to off-target nucleosomes in the panel. These results were reflected in the sample sequencing data (**Figure 14B**), as Samples 1-3 generated expected peak structures and NGS results were consistent across samples. In contrast, Sample 4 had low signal-to-noise (**S:N**) in both the SNAP-CUTANA Spike-in data and the sample data (see next question).

Question: What do failed assays look like using SNAP-CUTANA Spike-ins?

In a failed experiment, an antibody with previously characterized specificity exhibits unusual behavior, such as cross-reactivity to an off-target PTM and/or high background. To demonstrate this in an experimental context, reference Sample 4 in **Figure 14**. In this experiment we used well-characterized H3K4me3 and H3K27me3 antibodies and added the

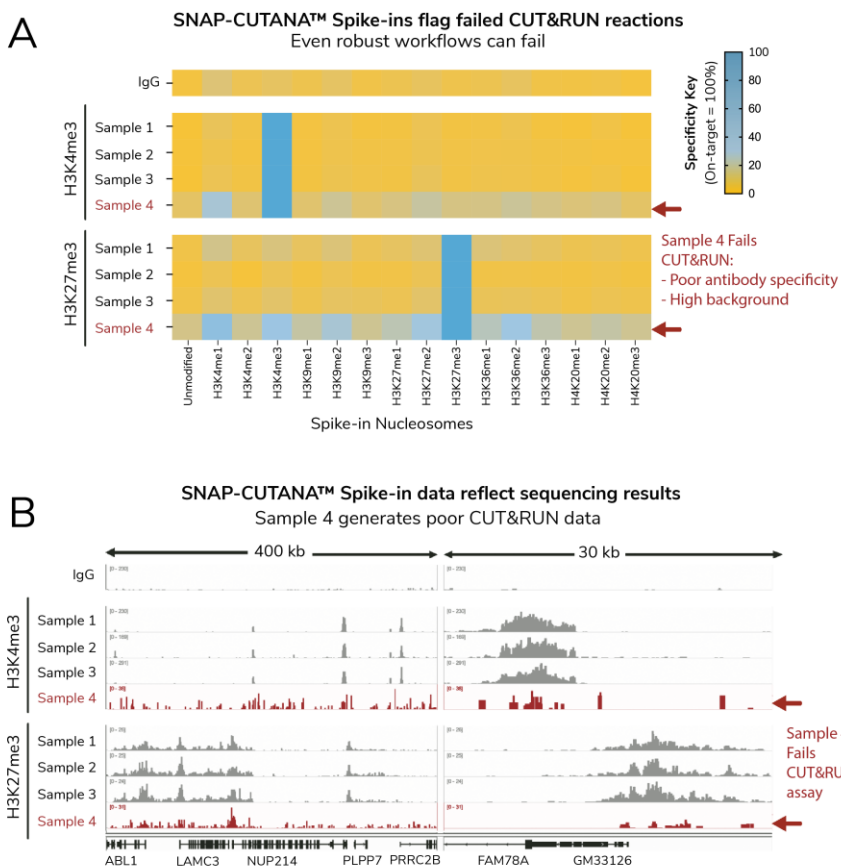


Figure 14: SNAP-CUTANA™ K-MetStat Spike-in Controls identify failed CUT&RUN reactions. CUT&RUN was used to map H3K4me3 and H3K27me3 in mouse B cells (four independently processed samples). **(A)** SNAP-CUTANA Spike-ins were included in each reaction. Heatmaps show antibody binding relative to on-target PTM; Sample 4 displays increased cross-reactivity (red arrows). **(B)** RPKM-normalized tracks from sample chromatin shows consistent peaks for Samples 1-3, while Sample 4 displayed low S:N and excessive background (red arrows).

K-MetStat Spike-in Panel to each reaction. Samples 1-3 show high antibody specificity. In contrast, Sample 4 displayed uncharacteristic high recovery of multiple off-target PTMs in the panel for both H3K4me3 and H3K27me3 reactions (**Figure 14A**, red arrows), suggesting high background and low S:N in these CUT&RUN reactions. When analyzing sample NGS

data, we observed similarly low S:N for Sample 4, for both H3K4me3 and H3K27me3 (**Figure 14B**, red arrows).

Of note, all eight CUT&RUN reactions were performed in parallel using the same antibodies, but only Sample 4 displayed problems with S:N. In addition, both H3K4me3 and H3K27me3 reactions using Sample 4 cells displayed cross-reactivity against SNAP-CUTANA Spike-ins. Combined, these results suggest overall problems with the sample material vs. a simple failed reaction, and we focused our troubleshooting approaches on this aspect (in this case the consortium lab had delivered less than the expected 10,000 cells).

Importantly, these results illustrate the utility of including SNAP Spike-in Controls in all experiments. It may be unclear from the genome tracks alone that a reaction issue had occurred. **SNAP-CUTANA Spike-ins can flag failed reactions and indicate the cause.** By flagging failed samples and using the SNAP Spike-in results to guide troubleshooting, researchers can be confident in their experimental results.

Question: *Why are SNAP-CUTANA Spike-ins better than other methods to identify failed reactions?*

Although outliers can be identified using other methods (e.g. separation in Principal Component Analysis [PCA], poor Pearson correlation), SNAP-CUTANA Spike-in Controls provide an added layer of experimental context. PCA/Pearson are only a proxy for assay stability and cannot determine which part of the experiment is at fault (e.g. antibody specificity, cell preparation, workflow). In contrast, SNAP-CUTANA Spike-in Controls provide a direct and quantitative readout of experimental success, guiding troubleshooting or providing confidence for researchers to proceed with data analysis and interpretation.

Question: *How can I use SNAP-CUTANA Spike-ins to guide troubleshooting?*

When the spike-in controls work as expected, users can trust that their antibody was specific, and most CUT&RUN/CUT&Tag experimental conditions are optimal. If the spike-in results **and/or** genomic enrichment patterns are not as expected, the specific form of the data can provide valuable insights for diagnosing and troubleshooting the problem (**Table 8**).

Spike-in results	Genomic results	Result confirms	Troubleshooting approach
Specific, high S:N	Poor S:N	<ul style="list-style-type: none"> ✓ pAG-MNase cleavage confirmed ✓ Optimal wash conditions ⚠ Cell integrity may be in question 	<ul style="list-style-type: none"> ➤ Confirm cell viability ➤ Try higher cell numbers ➤ Ensure target is present & localized to chromatin ➤ Use native or lightly cross-linked cells ➤ Optimize permeabilization conditions
Recover off-target PTMs	Good S:N, but unexpected peaks	⚠ Antibody cross-reactivity	➤ Test additional antibodies
Poor S:N	Poor S:N	⚠ Fundamental workflow failure	<ul style="list-style-type: none"> ➤ Restart with recommended control cells & antibodies ➤ Assess bead clumping ➤ Prepare fresh buffers ➤ Try using our kit (EpiCypher 14-1048)
Specific, high S:N	High S:N, expected peaks	➤ Happy data analysis!	✗ None needed!

Table 8: Interpreting SNAP-CUTANA™ Spike-in and NGS results to guide troubleshooting. S:N, signal-to-noise.

4. Use SNAP-CUTANA Spike-ins to normalize data and quantitatively compare samples.

In addition to profiling antibody specificity, guiding assay development, and monitoring assay success, SNAP-CUTANA Spike-ins can also be used to normalize NGS data for reliable, quantitative cross-sample comparisons. EpiCypher is currently investigating multiple normalization methodologies and will share recommendations when available. Below, we describe an emerging application of this exciting technology.

Question: How does normalization to spike-ins compare with other normalization methods?

To demonstrate the application of SNAP-CUTANA Spike-ins for NGS normalization and drug response studies, we show a direct comparison of CUT&Tag data normalization using the K-MetStat Spike-ins vs. standard RPKM methods (**Figure 15**). Briefly, CAL27 squamous cell carcinoma cells were treated with either vehicle (0hr) or 3 μM of the EZH2 methyltransferase inhibitor tazemetostat (24, 48hr). CUT&Tag was performed with an antibody to H3K27me3 (CST #9733) using the EpiCypher CUTANA™ Direct-to-PCR CUT&Tag Protocol (epicypher.com/protocols). The SNAP-CUTANA K-MetStat Panel was added to each sample prior to antibody addition. Sequencing data were RPKM normalized (top panel) or normalized to the K-MetStat Spike-in Controls (similar to reported methods²; bottom panel).

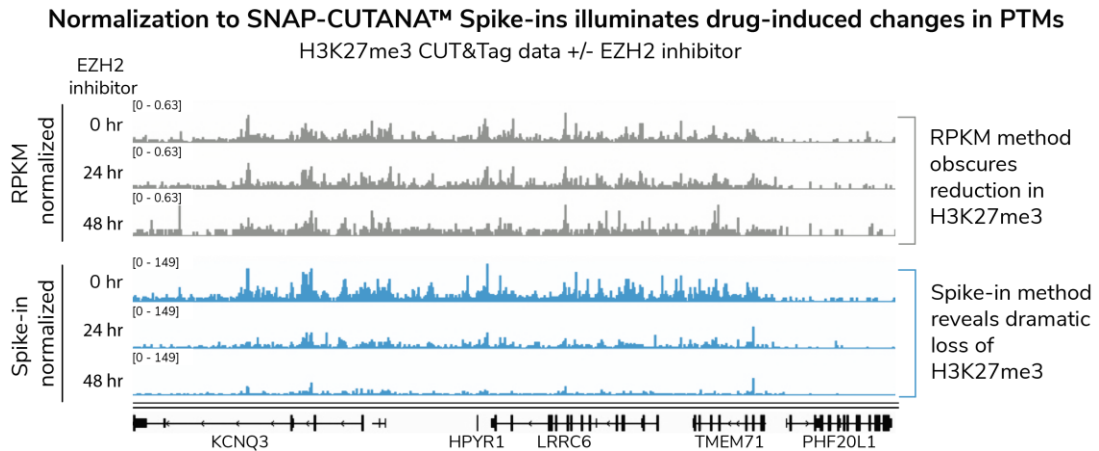


Figure 15: CUT&Tag data normalization using SNAP-CUTANA™ K-MetStat Spike-in Controls illuminates drug-induced differences in histone PTM enrichment. Cancer cells treated with the EZH2 inhibitor tazemetostat for the indicated times show a modest difference in H3K27me3 enrichment compared to vehicle (0hr) when using RPKM normalization. However, spike-in normalization reveals a pronounced and time-dependent reduction in H3K27me3 enrichment following treatment. *Special thanks to Drs. Yinglu Li, Xiao Chen and Chao Lu (Columbia University) for performing the experiment and sharing the data.*

For spike-in normalization, a scale factor was calculated for each sample by dividing the percent of total reads aligned to human genome by the percent of total reads aligned to the spike-in barcodes (Scale Factor = % Human Reads / % Spike-in Reads) and applying this factor to adjust the total sequencing reads of each respective sample. A very modest difference in H3K27me3 enrichment after EZH2 inhibition is observed in RPKM normalized samples. In contrast, spike-in normalization reveals a pronounced and time-dependent reduction in H3K27me3 enrichment following tazemetostat treatment, consistent with immunoblot analysis of total H3K27me3 levels (not shown). Thus, normalization using SNAP-CUTANA Spike-ins can illuminate drug-induced differences in histone PTM enrichment that are not discernable by other approaches.

Q: What resources are there for developing normalization strategies using SNAP-CUTANA Spike-ins?

Several methods have been reported for experimental normalization using exogenous spike-ins; these approaches can be adopted by calculating a single scalar normalization ratio using total read counts from SNAP-CUTANA Spike-ins:

SNAP-ChIP:

Tay *et al.* Hdac3 is an epigenetic inhibitor of the cytotoxicity program in CD8 T cells. *J. Exp. Med.* 217, e20191453 (2020). (PMID: [32374402](https://pubmed.ncbi.nlm.nih.gov/32374402/))

Lam *et al.* Cell-type-specific genomics reveals histone modification dynamics in mammalian meiosis. *Nat. Commun.* 10, 3821 (2019). (PMID: [31444359](https://pubmed.ncbi.nlm.nih.gov/31444359/))

ChIP-Rx:

Orlando *et al.* Quantitative ChIP-Seq Normalization Reveals Global Modulation of the Epigenome. *Cell Rep.* 9, 1163-1170 (2014). (PMID: [25437568](#))

- 5. Use SNAP-CUTANA Spike-ins to push the boundary.** At EpiCypher, we use SNAP-CUTANA Spike-ins in all CUT&RUN and CUT&Tag reactions to monitor assay stability, define antibody specificity, and much more. We envision SNAP-CUTANA Spike-ins as an essential tool in broadening the clinical applications of CUT&RUN and CUT&Tag assays, particularly for low-input assays and the use of valuable patient samples. As these new technologies develop, quantitative controls are crucial to standardize workflows across lab environments and/or different experiments.

Use SNAP-CUTANA Spike-in Controls to address key problems and tackle advanced applications in the field:

- Optimize CUT&RUN/CUT&Tag assays for reduced cell numbers or rare sample types
- Validate your workflow for new cell types/preparation methods
- Standardize CUT&RUN/CUT&Tag assays for quantitative clinical experiments
- Identify and avoid widely cited antibodies that are not fit for CUT&RUN/CUT&Tag assays

Section III: How to Analyze Sequencing Data Using SNAP-CUTANA™ Spike-in Controls

- 1. Align paired-end sequencing reads to the SNAP-CUTANA Spike-in Controls and determine the number of reads assigned to each barcoded nucleosome in the panel.** Each spike-in nucleosome contains a unique, PTM-specific barcode (**Figure 8**) that will be detected by NGS. Importantly, each modified nucleosome is assembled using two distinct DNA barcoded templates, providing a technical replicate within each reaction. Thus, for the K-MetStat Spike-in Panel, which contains 16 PTM states, there are 32 DNA barcodes.

These barcode sequences are not contained in mouse, human, fly, or yeast genomes, and thus require an independent alignment procedure. In addition, depending on the directionality of adaptor ligation during library prep, the barcode reads may be contained with R1 reads (adaptor P5) or R2 reads (adaptor P7) from paired-end sequencing. As a result, both R1 and R2 fastq files should be aligned to DNA barcodes.

These steps are easily accomplished using the shell script available on the respective SNAP-CUTANA panel product page. A shell script is a .sh extension file that can be opened with any basic text editor program (*e.g.* TextEdit on a Mac or any text editing app). It should not be opened or saved as a PDF/Word Doc.

The instructions for aligning sequencing reads to the spike-in DNA barcodes are contained in the shell script, and expanded upon here:

- a. In Finder (on a Mac), duplicate the template shell script provided by EpiCypher to save an experiment-specific copy. It should be saved as a .sh file.
- b. Download your R1 and R2 fastq.gz files. Unzip (extract) all fastq.gz to fastq on Mac or Linux by double-clicking on the .gz files. Place these fastq files and your shell script in a specific folder on your computer.
- c. Open your experiment-specific shell script file in a text editor program.
- d. Within the provided shell script, below the instructions, there is a “template loop” for barcode alignment. Copy&Paste the lines between "# template loop begin ##" and "# template loop end ##" (from the first "echo" to the last "done") in the shell script as many times as needed: you will need 1 template loop per paired-end R1 & R2 data set.

Example: For three sequencing reactions, copy the template loop three times.

- e. Next, replace the template file names in the loops (Sample1_R1.fastq and Sample1_R2.fastq) with your R1 and R2 file names. Each loop should contain matched R1 and R2 file names for a given sequencing reaction. Add the R1 fastq. file name in the first section of the loop, and the R2 fastq. file name in the second section of the loop.

Example: I replace the dummy file names with file names for my H3K4me3 experiment. File names “H3K4me3_Rep1_R1.fastq” and “H3K4me3_Rep1_R2.fastq” go in the first loop, “H3K4me3_Rep2_R1.fastq” and “H3K4me3_Rep2_R2.fastq” go in the second loop, and so on.

- f. If using a Linux or Mac operating system, use Terminal to run your script, per the instructions in the shell file. In Terminal, be sure to change the directory of Terminal to the location of your fastq. files.

Example: My files are saved to my Desktop, in a folder called “NGS Analysis.” On a Mac, I would open Terminal and type “cd” followed by a space, and then drag the folder onto the Terminal window to copy the file path. Press Return.

Alternatively, you can also type in the path to your folder containing the fastq files (eg: cd path_to_fastq).

- g. To execute your shell script in the Terminal application, type: “sh” followed by a space, and then drag the .sh file into Terminal to copy the file path. Press return to run the script. Again, you can also type this in manually by typing “sh filepath_to_shell.sh”. Press return to run the script.
- h. Terminal will generate read counts one “loop” (*i.e.* paired R1 & R2 file set) at a time, for all 32 barcodes. For each loop, the script will first generate read counts from R1 files, and then from the R2 files. The barcodes will be in the order listed under "# Barcode identities" in the shell script instructions, and datasets will also be annotated based on the filenames.
- i. Now you can move forward with analysis in Excel.

2. Use the provided Excel file on the SNAP-CUTANA Product Page to generate a heatmap of the spike-in reads. When the barcode counts generated in Terminal are copied and pasted into the appropriate cells of the Excel file, a heatmap of the results (*e.g.* similar to that shown in **Figure 11A**) will automatically be generated. This heatmap will provide information

on the success of control reactions and antibody specificity. Instructions on using the Excel template:

- a. The template consists of two sheets: the “K-MetStat” sheet, where you paste in your R1 and R2 data, and the “Output Table” which generates the full antibody specificity heatmap.
- b. We provide space to copy in SNAP-CUTANA read count data for an IgG negative control, an H3K4me3 positive control, and 6 additional reactions (scroll down). Copy and paste to create additional sample analyses as needed.
- c. Select your target(s) in the Excel sheet from the drop-down menu in Column B. Note that the Target is pre-set for the IgG and H3K4me3 analyses.
- d. For each reaction, copy the R1 and R2 barcode read counts generated from running the script and paste into the appropriate highlighted yellow cells in Excel. Terminal generates the R1 and R2 read counts in the order we have provided on the Excel Template.
- e. Once your Target is selected and your R1 and R2 reads are pasted in, the template will auto-generate a heatmap analysis normalized to the Target PTM (*i.e.* on-target will be set as 100% binding). Note that for IgG, binding is automatically normalized to the sum total of barcode reads.
- f. Assess antibody binding specificity, starting with the control reactions. Antibody specificity data for all 8 samples are summarized in one large heatmap, found in the “Output Table” sheet in the Excel Template. Some guidance on using the controls:
 - IgG should not display specific enrichment for any SNAP-CUTANA Spike-in (all boxes yellow/orange).
 - H3K4me3 should display specific recovery of the H3K4me3 spike-in (blue) and <20% binding to all other PTMs (yellow).
 - Confirmation of these controls is considered a general readout of workflow success: *i.e.* cells were prepared properly, your pAG-MNase or pAG-Tn5 enzyme behaved as expected, library preparation and sequencing proceeded normally, etc.
- g. Once these controls have been confirmed, proceed to analysis of other reactions/antibodies. Each antibody should display 100% binding to target (blue), and <20% binding to off-target PTMs.
 - **EpiCypher considers an antibody with <20% binding to all off-target PTMs specific and suitable for downstream data analysis.**
- h. Finally, we recommend calculating the percentage of total reads attributed to SNAP-CUTANA Spike-ins for each reaction. To determine the percent of total reads assigned to the spike-ins, fill in the “Uniq align reads” with the full number of uniquely aligned reads from the NGS reaction. The total number of barcode reads will be auto-filled from R1 and R2 counts for each reaction. Percentages (% total barcode reads) will auto-populate. This is also referred to as the spike-in “bandwidth” and should be ~1% for reliable analysis of antibody specificity. However, the range may be higher or lower depending on target abundance and antibody used. For example:
 - For H3K4me3 (low abundance target in cells), panel barcode reads are typically 1-10% of total sequencing reads.



- For H3K27me3 (high abundance target in cells), panel barcode reads are typically 0.1-1% of total sequencing reads.
- For IgG negative control antibody (no target present in sample), panel barcode reads are typically 10-20% of total sequencing reads.

Outside of this range, consider adjusting the spike-in dilution to be optimal for future experiments. The main goal is that thousands of sequencing reads are aligned to the spike-ins for adequate sampling of the panel and reliable use in antibody specificity assessment and data normalization.

Appendix III References

¹ Shah, R. N. *et al.* (2018) Examining the Roles of H3K4 Methylation States with Systematically Characterized Antibodies.

² Orlando *et al.* (2014) Quantitative ChIP-Seq Normalization Reveals Global Modulation of the Epigenome.

Appendix IV: Use of *E. coli* Spike-in DNA for Experimental Normalization

Description: This Appendix describes the optimization and application of *Escherichia coli* (*E. coli*) Spike-in DNA ([EpiCypher 18-1401](#)) for CUT&RUN sequencing normalization, as noted in **Section V** of the CUTANA CUT&RUN Protocol. This protocol should be used in combination with **CUTANA pAG-MNase for ChIC/CUT&RUN** ([EpiCypher 15-1016](#)), which has very low levels of *E. coli* DNA. These steps were developed using 500,000 native K562 cells.

Protocol: Optimizing amount of *E. coli* Spike-in DNA for CUT&RUN

1. Add *E. coli* spike-in DNA to CUT&RUN **Stop Buffer** prior to sample addition by preparing a **Stop Buffer Master Mix** (see **Section V**). The amount of spike-in added should consider cell/nuclei input, condition (fixed, native, etc.), and target abundance. Thus, optimization may be required. Some general guidelines:
 - For reactions using 500,000 cells add 0.5 – 1 ng *E. coli* spike-in DNA/reaction. This works for most cell types and conditions.
 - The spike-in amount should be scaled linearly from the 500,000 cell input baseline; e.g. using 0.05 – 0.1 ng for 50,000 cells.
 - The major goal is to aim for the spike-in to comprise ~1% (0.2-5%) of total sequencing reads (**Table 9**).

<i>E. coli</i> Spike-in DNA	Target	Total Reads	<i>E. coli</i> Reads	% <i>E. coli</i> Reads
0.5 ng	IgG	3,644,233	155,549	4.27%
	H3K4me3	3,121,112	42,210	1.35%
	H3K27me3	5,254,299	8,511	0.16%
1.0 ng	IgG	2,569,291	241,645	9.41%
	H3K4me3	3,127,912	147,565	4.72%
	H3K27me3	9,650,258	22,419	0.23%

Table 9. EpiCypher *E. coli* Spike-in DNA (0.5 and 1.0 ng) was added to CUT&RUN samples using 500,000 K562 cells enriched for a low abundance target (H3K4me3, EpiCypher 13-0041), a high abundance target (H3K27me3, EpiCypher 13-0030) and IgG negative control (EpiCypher 13-0042). Total numbers of paired-end sequencing reads, reads aligned to *E. coli*, and percentage of total sequencing reads aligned to *E. coli* spike-in DNA are shown. Green boxes highlight the spike-in amounts recommended for each target.

2. At the appropriate step in the CUTANA CUT&RUN protocol (see **Section V**), quench MNase activity with 33 μ L **Stop Buffer Master Mix** containing *E. coli* spike-in DNA.
3. Complete the CUT&RUN protocol and prepare Illumina sequencing library.

Protocol: NGS Normalization using *E. coli* Spike-in DNA

Description: To use spike-in *E. coli* DNA for normalization and pairwise comparisons perform the following steps.

4. Align NGS data to the experimental sample reference genome (e.g. human hg19 build) and to the *E. coli* K12, MG1655 reference genome (see Illumina iGenomes Reference Sequences; support.illumina.com/sequencing/sequencing_software/igenome.html).
5. For pairwise comparisons, determine the number of *E. coli* spike-in DNA reads and normalize to total sequencing depth (i.e. total number of uniquely mapped reads). Calculate this ratio for each individual NGS dataset.
 - Example: CUT&RUN was performed to map a histone PTM target from treated and untreated cell samples.
 - Treated spike-in: 100,000 *E. coli* reads in 5,000,000 uniquely aligned reads = 2%
 - Untreated spike-in: 30,000 *E. coli* reads in 3,000,000 uniquely aligned reads = 1%
6. Calculate normalization factor^A, such that after normalization the *E. coli* spike-in signal is equal across all samples. Continuing with the example above:
 - Treated normalization factor = 1/2% spike-in bandwidth = 0.5
 - Untreated normalization factor = 1/1% spike-in bandwidth = 1.0
7. Use single scalar normalization ratio with the --scaleFactor option enabled in the [deeptools bamCoverage](https://deeptools.readthedocs.io/en/develop/content/tools/bamCoverage.html) tool to generate normalized bigwig files for visualization in IGV (<https://deeptools.readthedocs.io/en/develop/content/tools/bamCoverage.html>). Example:
 - Treatment sample --scaleFactor = 0.5
 - Untreated sample --scaleFactor = 1.0

The effect of normalization on a sample is inversely proportional to the *E. coli* spike-in bandwidth. In other words, samples with the highest bandwidth will receive the largest reduction in signal after normalization. For further information on experimental normalization using exogenous spike-in controls, see references below^{A,B}.

Appendix IV References

^A Tay *et al.* (2020) Hdac3 is an epigenetic inhibitor of the cytotoxicity program in CD8 T cells.

^B Orlando *et al.* (2014) Quantitative ChIP-Seq Normalization Reveals Global Modulation of the Epigenome.