

EpiCypher® CUTANA™ CUT&RUN Protocol

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

1. Overview

Cleavage Under Targets & Release Using Nuclease (CUT&RUN) is a revolutionary genomic mapping strategy developed by the group of Dr. Steven Henikoff ¹. It builds on Chromatin ImmunoCleavage (ChIC) from Dr. Ulrich Laemmli ², whereby a fusion of Protein A to Micrococcal Nuclease (pA-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 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**).

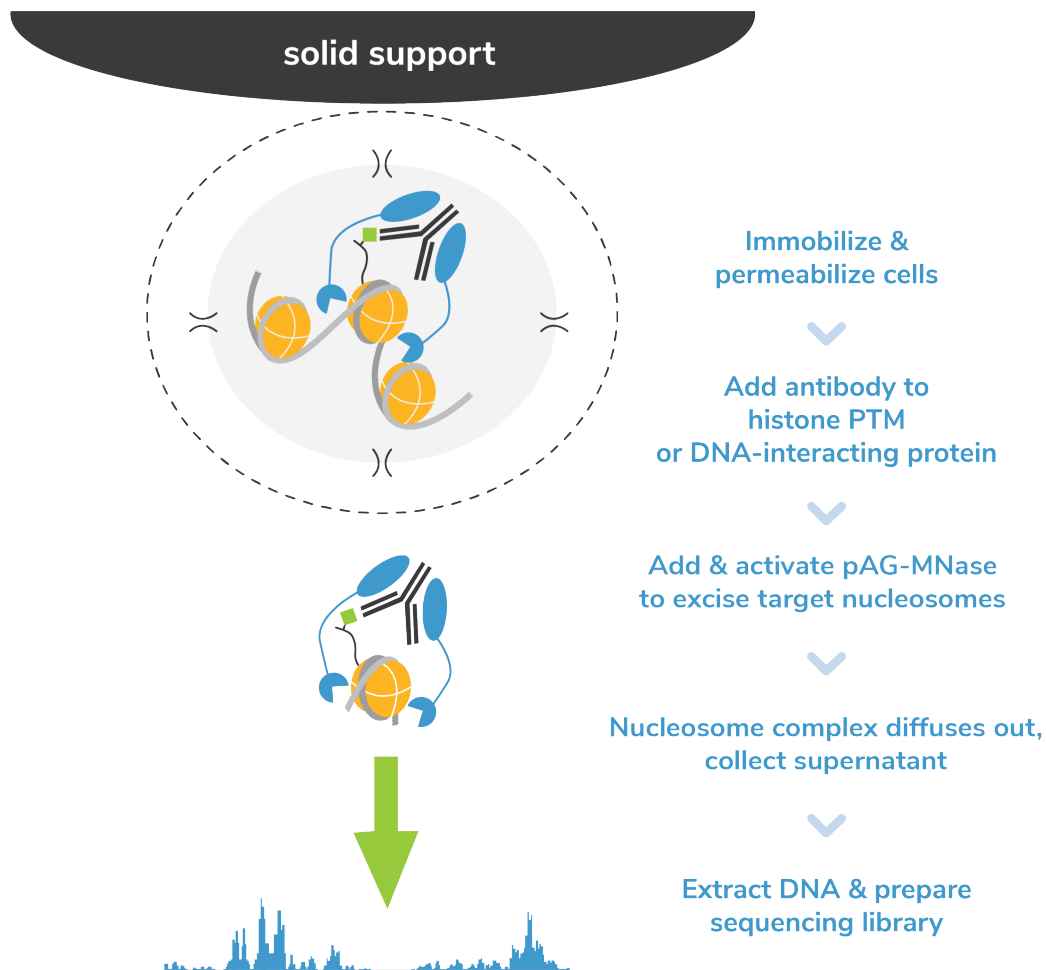


Figure 1. Overview of the CUTANA™ CUT&RUN protocol.

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:

- Requires substantially fewer cells
- Compatible with low sequencing depths
- Markedly improved [signal : noise]
- Rapid workflow: cells → data in < 2 days
- Empowers benchtop sequencers (e.g. Illumina® MiniSeq or MiSeq)
- Dramatically increases experimental throughput (transforming the way scientists approach genomic mapping studies)

EpiCypher now offers CUTANA™ pAG-MNase, the essential reagent for ChIC/CUT&RUN workflows:

- First-in-class commercial product for ChIC/CUT&RUN assays
- Optimized fusion of Proteins A and G with Micrococcal Nuclease (pAG-MNase) is directly compatible with a broad range of antibody isotypes
- 50 and 250 reaction pack sizes, enabling greater experimental throughput
- Check www.epicypher.com for regular updates of optimized protocols
- Additional CUTANA™ products coming soon:
 - CUT&RUN Compatible Antibodies
 - CUT&RUN Kits
 - CUT&RUN spike-in controls
 - pAG-Tn5 (for CUT&Tag applications)
 - *Inquire for more information or to connect to EpiCypher scientists:*
info@epicypher.com

3. Protocol Notes

This protocol has been validated for genomic profiling of:

- Histone PTMs (e.g. lysine methylation, lysine acetylation, lysine ubiquitylation, etc.)
 - Transcription factors (e.g. CTCF)
 - Chromatin remodelers (e.g. SMARCA4-SWI/SNF, ISWI, INO80, CHD)
 - Chromatin writers & readers (e.g. MLL, BRD4)
1. Always include control conditions (e.g. positive control cells & antibody; see FAQ section). This is especially critical when evaluating CUT&RUN using previously untested targets and/or antibodies.
 2. Protocol optimized using 0.5 million human K562 cells per sample. We have also validated variations to the protocol for working with nuclei, adherent cells, cryopreserved cells/nuclei, and crosslinked material (available upon request).
 3. Protocol adapted from 1.5 ml tubes to 300 µl 8-strip PCR tubes to: **a)** minimize beads sticking to tubes; **b)** enable more rapid workflow with multichannel pipets; **c)** provide more consistent sample handling; and **d)** enable high-throughput sample preparation.
 4. Take caution throughout to avoid ConA beads sticking to the sides / caps of tubes: they dry out easily, which can result in sample loss.
 5. Although protocols with shortened antibody and/or CaCl₂ incubation times have been published³, in our hands such changes adversely impact yield and reproducibility.
 6. We recommend a 5% digitonin stock solution be prepared in DMSO (as opposed to heated H₂O) because this dramatically increases detergent solubility and standardizes cell permeabilization / protocol reproducibility.
 7. To avoid digitonin precipitation and cell lysis, use the minimal concentration of digitonin in **Digitonin Buffer / Antibody Buffer**, which are required for efficient cell permeabilization. This should be empirically determined for different samples before proceeding with the full CUT&RUN experiment.

4. Buffers, Reagents & Materials Needed

Buffer recipes

Bead Activation Buffer

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

Wash Buffer

20 mM HEPES, pH 7.5
150 mM NaCl
0.5 mM Spermidine
1x Roche Complete Protease Inhibitor-mini (CPI-mini), EDTA-free (Roche catalog # 11836170001), 1tab/10ml
Filter sterilize
Store @ 4°C for up to 1 week

Digitonin Buffer

Wash Buffer + 0.01% Digitonin *
* Optimal [Digitonin] for each cell type should be empirically determined. Starting concentration validated for K562, MCF7, and A549 cells = 0.01% digitonin.
Prepare fresh each day and store @ 4°C

Antibody Buffer

Digitonin Buffer* + 2 mM EDTA
* See note above for empirically determining optimal digitonin concentration for a given cell type
Prepare fresh each day and store @ 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 @ 4°C for up to 6 months

Reagents

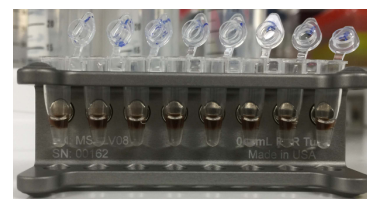
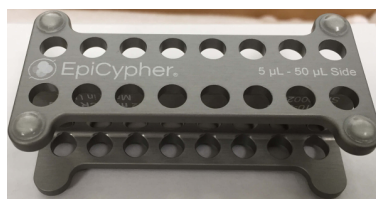
Item	Vendor	Catalog No.	Notes
Concanavalin A (ConA) Magnetic Beads	EpiCypher	21-1401	ConA is a lectin, which can cause immune activation. For technical support re. immune cell studies, contact info@epicypher.com
pAG-MNase	EpiCypher	15-1016	Supplied as 20X stock
Rabbit IgG Negative Control Antibody	EpiCypher	13-0042	Dilute 1:10 in Antibody buffer, store @ 4°C for 1 month.
H3K4me3 Positive Control Antibody: SNAP-ChIP Certified, CUTANA CUT&RUN Compatible	EpiCypher	13-0041	0.5 µg/µl → Use at 1:100 dilution in CUT&RUN.
Antibody to histone PTM, TF, or chromatin regulator	User-dependent		EpiCypher continues to conduct <i>extensive</i> antibody characterization (most particularly for those against histone PTMs ⁴). Contact us for recommendations: info@epicypher.com
NEB Monarch PCR & DNA Cleanup Kit	New England Biolabs	T1030S	Chosen for high yields of low [DNA] samples, low elution volume, and retention of >50bp fragments (thus suitable for nucleosomal (PTMs) and subnucleosomal (TFs) fragments)
Agencourt AMPure XP magnetic beads	Beckman Coulter	A63880	For PCR cleanup and size selection (i.e. 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

Equipment

Item	Vendor	Catalog No.	Notes
8-strip PCR tube magnetic stand	EpiCypher	10-0008	For all bead separation steps
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	Or comparable capillary electrophoresis instrument (e.g. Agilent TapeStation)
High Performance Multi-Channel Pipettors, 8-Channel	VWR	76169-252	For aspiration and wash steps to accelerate workflow
Tube Nutator	VWR	82007-202	For bead incubation steps (overnight antibody incubation, pAG-MNase digest reaction)
Vortex-Genie	Scientific Industries	SI-0236	For bead mixing steps.

EpiCypher's 8-strip PCR tube magnetic stand (#10-0008):

Enables streamlined sample handling for higher experimental throughput and improved reproducibility.



5.1. EpiCypher CUTANA: CUT&RUN Protocol (~5hrs)

Day 1

Section I: ConA Bead Activation (~30 min)

1. Gently resuspend the **ConA beads (Concanavalin A)** and transfer 11 μ l/sample to 1.5 ml tube for batch processing.
2. Place the tube on a magnet until slurry clears and pipet to remove supernatant (supe).
3. Add 100 μ l/sample cold **Bead Activation Buffer** and pipet to mix. Place the tube on a magnet until slurry clears and pipet to remove supe.
4. Repeat previous step for total of two washes.
5. Resuspend beads in 11 μ l/sample cold **Bead Activation Buffer**.
Split activated ConA beads into separate tubes for different cell types and/or antibodies.
6. Aliquot 10 μ l/sample of activated bead slurry into 8-strip tube. Keep beads on ice until needed.

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

7. Harvest 0.5 million cells/sample by spinning for 3 min at 600g at room temperature (RT) in 1.5ml tube, and decant supe.
8. Resuspend cells in 100 μ l/sample RT **Wash Buffer**; spin for 3 min at 600g at RT; decant supe.
9. Repeat previous step for total of two washes.
10. Resuspend cells in 100 μ l/sample in RT **Wash Buffer** and aliquot 100 μ l washed cells to each 8-strip tube containing 10 μ l of activated bead. Gently vortex (setting #7) to mix.
11. **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)

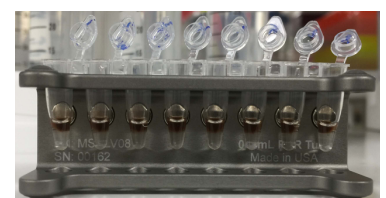
12. Place the tube on a magnet until slurry clears and pipet to remove supe.
13. Add 50 μ l cold **Antibody Buffer** to each sample and gently vortex.
14. Add 0.5 μ l **Antibody** (1:100 dilution) to each sample and gently vortex.
15. **Incubate** 8-strip tube (**caps slightly elevated**) on nutator **overnight at 4°C**.
 - a. **Critical step:** To keep beads in solution, slightly elevate cap side of 8-strip on nutator.



Bead slurry before o/n incubation @ 4C | nutator o/n | Bead slurry after o/n incubation @ 4C

Day 2

16. Place the tube on a magnet until slurry clears and pipet to remove supe.



17. While beads are on magnet, add 250 µl cold **Digitonin Buffer** directly onto beads of each sample, and then pipet to remove supe.
18. Repeat previous step for total of two washes.
19. Add 50 µl cold **Digitonin Buffer** to each 8-strip tube, and gently vortex. Beads are often clumpy at this point, but can easily be dispersed with gentle pipetting with a P200 pipet.

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

20. Add 2.5 µl **CUTANA™ pAG-MNase** (20x stock) to each sample, and gently vortex.
 - **Critical step:** *to evenly distribute pAG-MNase across cells/nuclei, ensure beads are thoroughly resuspended by gentle vortexing and/or pipetting with a P200 pipet.*
21. **Incubate** samples for **10 min at RT**, and then return 8-strip tube to magnet.
22. While beads are on magnet, add 250 µl cold **Digitonin Buffer** directly to each sample, and then pipet to remove supe.
23. Repeat previous step for total of two washes.
24. Add 50 µl cold **Digitonin Buffer** to each sample, and gently vortex. Beads are often clumpy at this point, but can easily be dispersed with gentle pipetting with a P200 pipet.

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

25. Place 8-strip on ice, add 1 µl **100 mM CaCl₂** to samples, and gently vortex.
 - **Critical step:** *For efficient pAG-MNase digestion, ensure beads are thoroughly resuspended by vortexing and/or pipetting with a P200 pipet.*
26. **Incubate** 8-strip tube on nutator for **2 hours at 4°C**. Key step wherein MNase tethered to chromatin is activated by CaCl₂ to cleave chromatin.
27. Add 33 µl **Stop Buffer** to each sample, and gently vortex to mix. Stops MNase activity by chelating Ca⁺⁺ ions.
28. **Incubate** 8-strip tube for **10 min at 37°C** in thermocycler. *Releases chromatin to supernatant and degrades RNA.*
29. Place 8-strip tube on a magnet stand until slurry clears and transfer supe to 1.5 ml tube.
30. Purify DNA using the **NEB Monarch DNA Cleanup Kit** as per manufacturer's instructions.
31. Elute DNA in 10 µl **Elution Buffer** and use 1 µl to quantify the CUT&RUN-enriched DNA using the **Qubit fluorometer** as per manufacturer's instructions.

Typical CUT&RUN DNA total yields from 500k native K562 cells, pre-PCR:

- *Note: Total yields are influenced by a number of considerations, including antibody specificity and efficiency, and epitope abundance in each cell type.*

IgG → ~ 3ng
H3K27me3 → ~200ng
H3K4me3 → ~50ng
CTCF → ~10ng
BRD4 → ~20ng
SMARCA4/BRG1 → ~15ng

5.2. Library prep (4 hrs), Bioanalyzer (1 hr), & Illumina sequencing

Section VI: NEB Ultra II Library Preparation (~4 hrs)

32. Using 5-10ng purified CUT&RUN-enriched DNA, prepare Illumina library using the **NEB Ultra II Library Prep Kit** per manufacturer's instructions (~4hrs).

- *Note: NEB Ultra II kit contains hot-start Q5 DNA polymerase.*

33. Overview of workflow, according to manufacturer's instructions.

- End repair
- Universal adapter ligation
- DNA cleanup using 1.1x Ampure beads to sample volume (retains fragments >150bp)
- PCR and primer indexing according to cycling parameters below

34. CUT&RUN-specific PCR cycling parameters:

- 45 sec @ 98°C → activation of hot-start Q5 polymerase
 - 15 sec @ 98°C → DNA melting
 - 10 sec @ 60°C → hybrid primer annealing & short extension (<700 bp)
 - Repeat "step b-c" for a total of 14x
 - 1 min @ 72°C → final extension
- **Critical step:** PCR cycling parameter are designed to enrich for DNA fragment within the sequence-able range (100bp-700bp). Thus, contaminating higher molecular weight DNA fragments are not enriched by PCR.

35. DNA cleanup using 1.1x Ampure beads to sample volume (eg: 55 µl Ampure to 50 µl PCR).

36. Elute DNA in 15 µl **0.1x TE buffer** and use 1 µl to quantify the purified PCR product using the **Qubit fluorometer** as per manufacturer's instructions.

- *Typical yield purified PCR DNA library, ~500-750ng (30-50ng/µl in 15µl)*

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

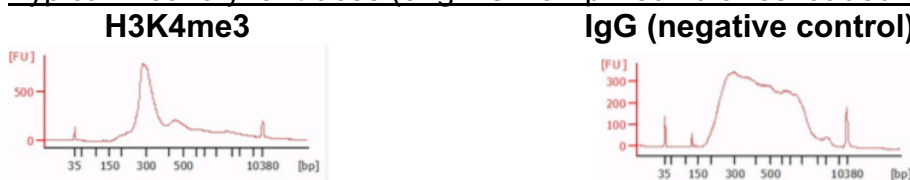
37. For each sample, prepare 5µl at 10ng/µl for loading on the Agilent Bioanalyzer.

- *Note: record dilution factor such that original sample molarity can be calculated from Bioanalyzer nM for desired DNA size range (100-1,000bp).*

38. For each purified PCR DNA library for Illumina sequencing, load 1ul of 10ng/µl sample on **Agilent High Sensitivity DNA Chip** (Cat# 5067-4626) as per manufacturer's instructions.

- *Typical molarity for 15 µl purified PCR DNA library (100-1,000bp region) = 100-200 nM*

Typical Bioanalyzer traces (5ng PCR amplified libraries loaded on Bioanalyzer)



Section VIII: Illumina sequencing

39. Based on Bioanalyzer molarity calculations, pool libraries at the desired ratios.

- Only 3-5 million paired-end (PE) reads are needed for good CUT&RUN coverage.

40. To obtain 3-5M PE reads for 8 samples, we typically load 0.8 pM (500 μ l) into a cartridge for a MiniSeq High Output Reagent Kit, 150-cycles (catalog # FC-420-1002).

- Flow cells using the MiniSeq High Output Kit typically yield 25-40M pass-filter clusters (i.e. 25-40M PE reads).

41. To obtain 3-5M PE reads for >48 samples, we typically load 0.8 pM (1,500 μ l) into a cartridge for a NextSeq 500/550 High Output Kit v2.5, 150-cycles (catalog # 20024907).

- Flow cells using the NextSeq High Output Kit v 2.5 typically yield 400M pass-filter clusters (i.e. 400M PE reads).

6. Buffer components

Components	Source	Cat #
HEPES	Sigma-Aldrich	H3375
KCl	Sigma-Aldrich	P3911
CaCl₂	Sigma-Aldrich	C1016
MnCl₂	Sigma-Aldrich	203734
Molecular biology grade H₂O (RNase, DNase free)	VWR	020201--500
NaCl	Sigma-Aldrich	S5150-1L
EDTA	Sigma-Aldrich	E5134
EGTA	Sigma-Aldrich	E3889
RNase A	Thermo Fisher Scientific	EN0531
Glycogen	Sigma-Aldrich (Roche)	Roche 10930193001
Pure Spermidine (6.4M)	Sigma-Aldrich	S0266
Digitonin	Millipore Sigma	300410
DMSO	Sigma	D8418-100ml
CaCl₂	Sigma-Aldrich	C1016
Trypan Blue	Thermo Fisher Scientific	T10282

7. Frequently Asked Questions (FAQs)

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 / negative controls in every experiment. If the QC checks for the positive control 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 500k K562 cells
 - ii. Positive control antibody (H3K4me3 or H3K27me3)
 - iii. Negative control antibody (IgG or no antibody)
- b. Experimental sample antibodies and cell types

Quality control checks before decision to sequence:

- a. Yield of CUT&RUN enriched DNA (~200 ng for positive control H3K27me3 antibody, ~3ng for negative control IgG antibody). See **Section V**.
- b. Yield of library prep and PCR (~750ng from 5-10ng library input). See **Section VI**.
- c. Bioanalyzer/Tapestation traces should be enriched with mono-nucleosomal size fragments (~275bp = ~150bp nucleosome+ 125bp adapters, see **Section VII**)

DNA spike-in controls

2. 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 in EpiCypher's pAG-MNase preps. However, at a typical sequencing depth of 3-5M reads, too few *E. coli* DNA fragments (~hundreds) are recovered for reliably computing sample normalization. Thus, we recommend considering other normalization strategies, such as exogenous DNA spike-in (*E. coli* / yeast / fly) or recombinant CUTANA nucleosomes (in development).
3. Does EpiCypher offer spike-in DNA controls for CUT&RUN?
EpiCypher plans to offer multiple DNA spike-in options for CUT&RUN, in addition to exogenous *E. coli* DNA spike-in. In particular, CUTANA spike-in nucleosomes (i.e. fully-defined semi-synthetic nucleosomes) for CUT&RUN are in active development and offer multiple advantages, including enabling sample input normalization and multiple control points (**1.** antibody specificity; **2.** antibody efficiency; **3.** pAG-MNase activity; and **4.** experimental success).

Cell input compatibility

4. 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 cancer lines.

EpiCypher has not yet directly tested, but a number of groups have successfully performed CUT&RUN on human and mouse primary tissue⁵⁻⁸, FACS sorted⁹ and immune cells^{10,11}.

Note that lectins (e.g. ConA-beads) play a role in the innate immune system and so immune cells types may be inadvertently stimulated via binding to ConA-beads. To circumvent this potential problem in CUT&RUN, EpiCypher recommends using nuclei⁹ or a crosslinking strategy¹².

5. Is CUT&RUN compatible with frozen or crosslinked cells?

Yes. 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. EpiCypher has also tested previously reported crosslinking conditions and recommended wash buffers¹², and found that while yields are significantly lower than native cells, the resulting data tracks display similar quality. We continue to explore crosslinking protocols and believe these will preserve labile PTMs, cell stimulation, or facilitate time-course studies.

Antibodies and targets

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

Yes. The current protocol has been used at EpiCypher to generate numerous non-PTM CUT&RUN data, including CTCF, BRD4, and SMARCA4 (BRG1). No protocol modifications were necessary to generate these data since the columns we use to purify the CUT&RUN-enriched chromatin fragments retain >50bp DNA.

However, for TFs in particular, which generate sub-nucleosomal size (<120bp) 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 sub-120bp fragments, the authors reduced the inactivation temperature after end repair from 65°C for 30m to 50°C for 1hr. Also, after the ligation reaction, Ampure bead size-selection ratio was increased from 1.1x volume to 1.75x volume. For TF CUT&RUN fragment sizes, both nucleosomal (~150bp) and sub-nucleosomal (<120bp) reflect TF chromatin occupancy locations, however, the sub-nucleosomal fragments provide higher resolution mapping for their locations.

7. What PTM antibodies do EpiCypher recommend for CUT&RUN?

We currently recommend SNAP-ChIP Certified histone PTM antibodies, since these display high specificity and efficiency in a ChIP experiment. However, we remain cautious about transferring ChIP capability directly to CUT&RUN, and our ongoing studies include the creation of *in situ* nucleosome standards to interrogate antibodies with the highest possible resolution. For non-PTM targets, we recommend testing a few different antibodies to each target and proceed with the reagent that generates the highest quality data.

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. & Fazzio, 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).