

CUTANA™ CUT&RUN Protocol for Histone PTMs

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), whereby a fusion of Protein A to Micrococcal Nuclease (pA-MNase) selectively cleaves antibody-bound chromatin *in situ* ². In CUT&RUN, cells or nuclei are immobilized to a solid support, with pA-MNase cleaved DNA fragments isolated from solution. The workflow is compatible with next-generation sequencing (NGS) to generate high resolution genome-wide profiles of histone post-translational modifications (PTMs) and chromatin-associated proteins (**Figure 1**).

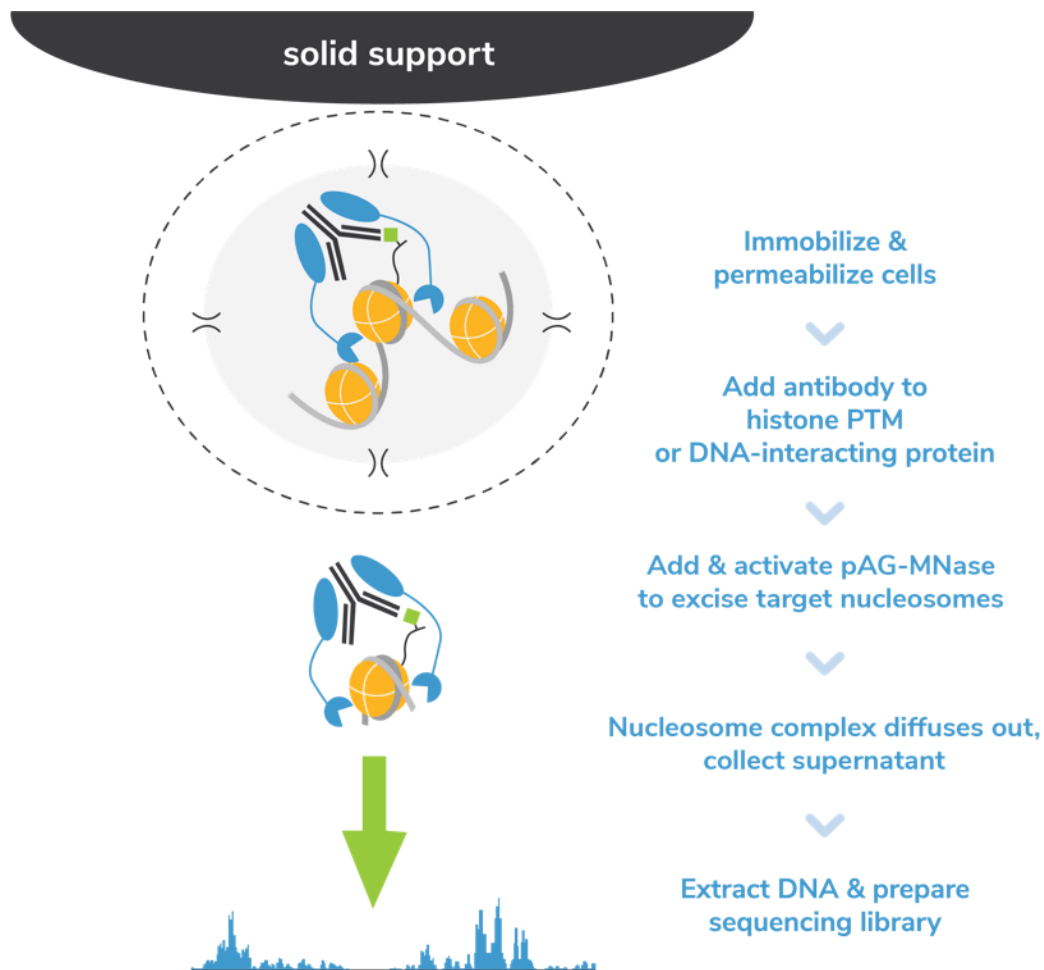


Figure 1. Overview of the CUTANA™ EpiCypher CUT&RUN protocol.

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 < 1 week
- Empowers benchtop sequencers (e.g. Illumina® MiniSeq or MiSeq)
- Dramatically increased experimental throughput (changing 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) enables direct compatibility with a broad range of antibody isotypes
- 50 and 250 reaction pack sizes, enabling greater experimental throughput
- Check back at www.epicypher.com for regular updates of optimized protocols
- Additional CUTANA™ products coming soon:
 - CUT&RUN Kits
 - CUT&RUN spike-in controls
 - pAG-Tn5 (for CUT&Tag applications)
 - CUT&Tag Kits
 - CUT&Tag spike-in controls
 - *Inquire for more information or to connect to EpiCypher scientists:*
info@epicypher.com

3. Specific Protocol Notes

1. Protocol is optimized using 0.5 million human K-562 cells per sample.
2. Although protocols with shortened antibody and/or CaCl₂ incubation times have been published, in our hands such changes adversely impact reproducibility.
3. Take caution throughout to avoid ConA beads sticking to the sides / caps of tubes, as they dry out easily and can result in sample loss.
4. We recommend a 5% digitonin stock solution be prepared in DMSO (as opposed to dissolution in heated H₂O).
5. To avoid digitonin precipitation use the minimal digitonin concentration in the **Digitonin Buffer** and **Antibody Buffer** required for efficient cell permeabilization (see Protocol Part III, Step 6). This should be empirically determined for different cell types before proceeding with the full CUT&RUN experiment.

4. Reagents & Materials Needed

Buffer Stocks

Item	Vendor	Catalog No.	Notes
HEPES	Amresco	7365	
KCl	EMD Millipore	PX1405	
CaCl ₂	Amresco	1B1110	
MnCl ₂	Sigma-Aldrich	M1787	
NaCl	Amresco	0241	
Spermidine	Sigma-Aldrich	S2501	Prepare 1 M stock in ddH ₂ O and make single use aliquots stored at -20°C. The stock is good for 1 month.
Protease Inhibitor	Sigma-Aldrich	11873580001	
Digitonin	EMD Millipore	300410	Prepare as 5% (wt/vol) stock in DMSO at room temperature (RT). The stock is good for 3 months when stored at -20°C.
EDTA	Amresco	M101	
EGTA	Amresco	0732	
Trypan Blue	Thermo Fisher Scientific	T10282	
RNase A	Thermo Fisher Scientific	5056489001	
Glycogen	Sigma-Aldrich	10901393001	

Reagents

Item	Vendor	Catalog No.	Notes
Concanavalin A (ConA) Magnetic Beads	Bangs Laboratories	BP531	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
Histone PTM Antibody	User-dependent		EpiCypher continues to conduct extensive studies of histone PTM antibodies ³ . Contact us for recommendations: info@epicypher.com
MinElute PCR Purification Kit	Qiagen	28004	
Qubit 1x dsDNA HS Assay Kit	Thermo Fisher Scientific	Q33230	
Library Preparation Kit	User-dependent		

Equipment

Item	Vendor	Catalog No.	Notes
Magnetic Stand	Thermo Fisher Scientific	12321D	
Qubit 4 Fluorometer	Thermo Fisher Scientific	Q33226	
Agilent 2100 Bioanalyzer	Agilent	G2939A	Or other capillary electrophoresis instruments (e.g. Agilent TapeStation)
ThermoMixer	Eppendorf	5384000020	
Tube Rotator	Grant-Bio	PTR-35	
Tube Nutator	Labnet	S0500	

5. CUT&RUN Protocol for Histone PTMs

Part I: ConA Bead Activation (<i>estimated time = 10 - 20 minutes</i>)	
<ol style="list-style-type: none"> 1. Gently resuspend the ConA magnetic beads and withdraw calculated total volume of the bead suspension (based on 10 μL per sample). 2. Mix into 1.5 mL of Binding Buffer in a 2 mL tube^a. 3. Place the tube on a magnet stand and remove the liquid. 4. Add 1.5 mL of Binding Buffer and mix by gentle pipetting. 5. Place the tube on a magnet stand and remove the liquid. 6. Resuspend the beads in (N + 1) x 10 μL Binding Buffer (where N = number of samples)^b. 	<ol style="list-style-type: none"> a. After each wash, remove the liquid from the cap by a pulse spin on a microcentrifuge. b. The activated beads may be held on ice for a day before use.
Part II: Binding Cells to Activated Beads^c (<i>estimated time = 0.5 – 1 hours</i>)	
<ol style="list-style-type: none"> 1. Count and harvest calculated number of fresh cells by centrifuging for 3 min at 600 x g^d. 2. Decant supernatant and resuspend cells in 1.5 mL room temperature (RT) Wash Buffer and transfer to a 2 mL screw-cap tube. 3. Centrifuge for 3 min at 600 x g. 4. Repeat steps 2 & 3 for two total washes. 5. Resuspend cells in 1 mL in RT Wash Buffer by gently pipetting. 6. While gently vortexing the cells, add activated bead suspension from Part I. 7. Place the cell bead mixture on a tube rotator and rotate at RT for 10 min. 	<ol style="list-style-type: none"> c. Keep buffers and steps in this section at RT to avoid cell stress. d. Tested at EpiCypher: 0.5 million human K-562 cells per sample.
Part III: Binding of Antibodies (<i>estimated time = 10 – 30 min + overnight</i>)	
<ol style="list-style-type: none"> 1. Quick spin, then place the tube on a magnet stand and remove the liquid. 2. Place the tube at a low angle on the vortex mixer, squirt (N + 1) x 50 μL Antibody Buffer along the side while gentle vortexing^e. 3. Gently tap to dislodge all beads. 4. Aliquot 50 μL bead suspension to N x Eppendorf tubes and add 0.5 - 1 μL antibody to each tube. 5. Place the tubes on the tube nutator and nutate overnight at 4°C^f. 6. (Optional) <i>Test cell permeabilization efficiency by digitonin^g:</i> After nutating the cell bead suspension in Antibody Buffer for 10 min at RT, mix 10 μL with 10 μL Trypan blue dye. Apply 10 μL to a glass slide and observe under a microscope. >95% cells should be stained blue if successfully permeabilized. 	<ol style="list-style-type: none"> e. Tested at EpiCypher: 0.01% digitonin and 1:100 dilution of antibodies (~ 0.5-1 mg/mL stocks) for 0.5 million K-562 cells. f. CRITICAL STEP: In contrast to shorter incubation times, overnight antibody incubation increases protocol reproducibility. g. Empirically determine the minimal [digitonin] in Digitonin buffer and Antibody buffer for efficient cell permeabilization.

<ol style="list-style-type: none"> 7. Quick spin, then place the tube on a magnet stand and remove the liquid. 8. Add 1 mL of Digitonin Buffer, thoroughly mix with the beads by inversion^h. 9. Quick spin, then place the tube on a magnet stand and remove the liquid. 10. Repeat steps 8 & 9 for a total of two washes. 	<p>h. Avoid pipetting as much as possible to avoid bead loss on pipet tips.</p>
--	--

Part IV: Binding of pAG-MNase (estimated time = 30 min)

<ol style="list-style-type: none"> 1. Place the tube at a low angle on the vortex mixer, squirt 50 µL of Digitonin Buffer along the side while gentle vortexing. 2. Gently tap to dislodge all beadsⁱ. 3. Add 2.5 µL of EpiCypher CUTANA™ pAG-MNase (20X stock). 4. Place the tubes on a tube nutator and nutate for 10 min at room temperature. 5. Quick spin, then place the tube on a magnet stand and remove the liquid. 6. Add 1 mL of Digitonin Buffer, thoroughly mix with the beads by inversionⁱ. 7. Quick spin, then place the tube on a magnet stand and remove the liquid. 8. Repeat steps 6 & 7 for a total of two washes. 	<p>i. Avoid pipetting as much as possible to avoid bead loss on pipet tips.</p>
---	--

Part V: Targeted Chromatin Digestion and Release (estimated time = 3 - 4 hours)

<ol style="list-style-type: none"> 1. Place the tube at a low angle on the vortex mixer, squirt 150 µL of Digitonin Buffer along the side while gentle vortexing. 2. Pre-chill the tubes and 100 mM CaCl₂ on ice. 3. Add 3 µL of 100 mM CaCl₂ to each tube with gentle vortexing and immediately replace the tube on ice. 4. Move the tubes to a tube nutator and nutate for 2 hours at 4°C^{j,k}. 	<p>CRITICAL STEPS:</p> <p>j. In contrast to shorter incubation times, 2 hour digestion increases protocol reproducibility.</p> <p>k. It is critical to keep the tubes at 0-4°C to minimize background chromatin cleavage by the pAG-MNase.</p>
<ol style="list-style-type: none"> 5. Add 100 µL of STOP Buffer and mix by gentle vortexing. 6. Incubate the tubes at 37°C for 10 min on a ThermoMixer at 500 rpm to release fragments into solution. 7. Centrifuge at 4°C for 5 min at 16,000 x g. 8. Place the tube on a magnet stand and transfer the liquid to a 2 mL tube. 	
<ol style="list-style-type: none"> 9. Extract the DNA with the Qiagen MinElute PCR Purification Kit by following the manufacturer's instructions^l. 10. Quantify the CUT&RUN DNA by Qubit analysis, following the manufacturer's instructions^{m,n}. 	<p>l. Use 10 µL EB buffer for the final elution.</p> <p>m. 2 - 5 ng total DNA is desired for the library preparation.</p>

	<p>n. Tested at EpiCypher:</p> <p>For H3K27me3 CUT&RUN, ~ 30-50 ng final DNA yield is typically expected;</p> <p>For H3K4me3 CUT&RUN, ~ 4-10 ng final DNA yield is typically expected.</p>
--	---

Part VI: Sequencing Library Preparation (estimated time = 1 – 2 days)	
<ol style="list-style-type: none"> 1. Prep indexed DNA libraries following the manufacturer’s instructions^o. 2. For PCR amplification of CUT&RUN fragments, perform the following procedures to enrich for mononucleosome sized fragments: <ol style="list-style-type: none"> a. 98°C 45 s, b. 98°C 15 s, c. 60°C 10 s, d. Repeat b-c for 13x, e. 72°C 1 min f. 8°C Hold 3. Quantify the DNA library by Qubit analysis, following the manufacturer’s instructions. 4. Assess the DNA library size distribution by Agilent 2100 BioAnalyzer analysis, following the manufacturer’s instructions^p. 	<p>o. Tested at EpiCypher: Both Illumina TruSeq ChIP Sample Prep Kit and NEBNext® Ultra™ II DNA Library Prep Kit for Illumina® work well for CUT&RUN library preparation.</p> <p>p. Typical BioAnalyzer traces for H3K27me3 and H3K4me3 after library preparation can be found in the EpiCypher CUTANA™ pAG-MNase Technical Data Sheet.</p>

6. Buffer Recipes

Binding 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 Protease Inhibitor
Filter sterilize
Store @ 4°C for up to 1 week

Digitonin Buffer

Wash Buffer + x% Digitonin*
*[Digitonin] needs to be empirically determined, see Protocol Part III, Step 6
Store @ 4°C for up to 1 day

Antibody Buffer

Digitonin Buffer + 2 mM EDTA
Store @ 4°C for up to 1 hour upon antibody addition

STOP Buffer

340 mM NaCl
20 mM EDTA
4 mM EGTA
0.02% Digitonin
50 µg/mL RNase A
50 µg/mL Glycogen
Filter sterilize
Store @ 4°C for up to 1 week

7. References

- ¹ Skene PJ, Henikoff JG & Henikoff S (2018) Targeted in situ genome-wide profiling with high efficiency for low cell numbers. *Nature Protocols* **13**:1006-1019 ([doi:10.1038/nprot.2018.015](https://doi.org/10.1038/nprot.2018.015)).
- ² Schmid M, Durussel T & Laemmli UK (2004) ChIC and ChEC; genomic mapping of chromatin proteins. *Molecular Cell* **16**:147-157 ([doi:10.1016/j.molcel.2004.09.007](https://doi.org/10.1016/j.molcel.2004.09.007)).
- ³ Shah RN, Grybowski AT, Cornett EM, Johnstone AL, Dickson BM, Boone BA, Cheek MA, Cowles MW, Maryanski D, Meiners MJ, Tiedemann RL, Vaughan RM, Sun ZW, Rothbart SB, Keogh M-C & Ruthenburg AJ (2018) Examining the roles of H3K4 methylation states with systematically characterized antibodies. *Molecular Cell* **72**:162-177 ([doi:10.1016/j.molcel.2018.08.015](https://doi.org/10.1016/j.molcel.2018.08.015)).