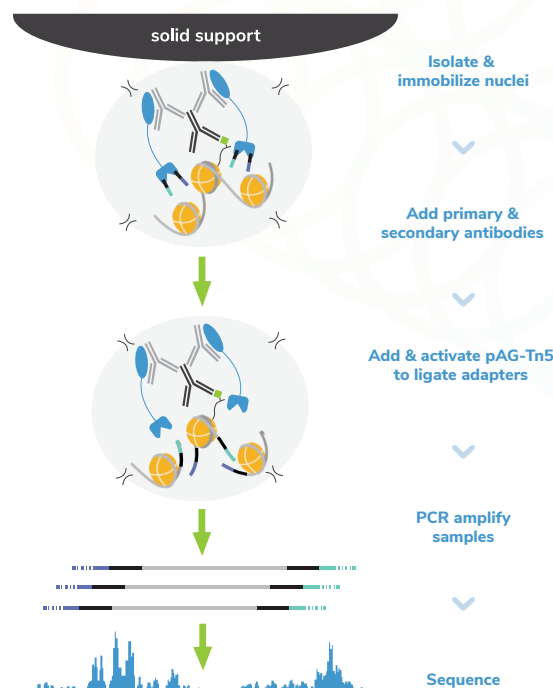


# CUTANA™ CUT&Tag Assays for chromatin mapping with low cell numbers

Cleavage Under Targets and Tagmentation (CUT&Tag) is an ultra-sensitive chromatin mapping technology that is ideal for histone post-translational modifications (PTMs) and select transcription factors (TFs).

How does CUT&Tag compare to ChIP-seq?

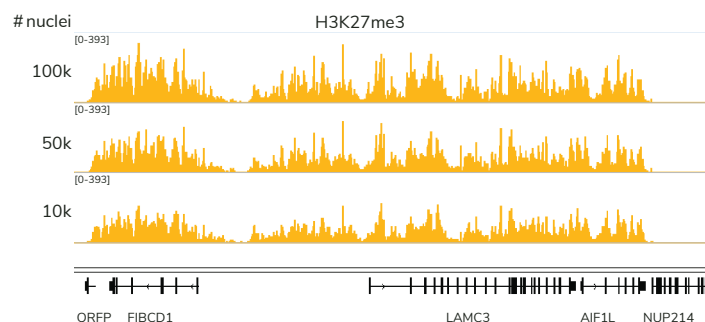
- Streamlined - no fragmentation, IP, or library prep
- Improved signal-to-noise
- Fewer cells needed
- Rapid 2-day workflow
- Reduced sequencing costs



For high-quality chromatin profiling, choose CUTANA™ Assays

FEATURES	ChIP-seq	CUT&RUN	CUT&Tag
Cells/nuclei required	>1 Million	5,000* - 500,000	10,000* - 100,000
Compatible targets	Histone PTMs, TFs	Histone PTMs, TFs & chromatin remodelers	Histone PTMs
Uniquely mapped reads	>30 Million	3-8 Million	5-8 Million
Signal-to-noise	Low	High	High

\* Success at lower inputs depends on antibody quality, cell type, and target abundance.

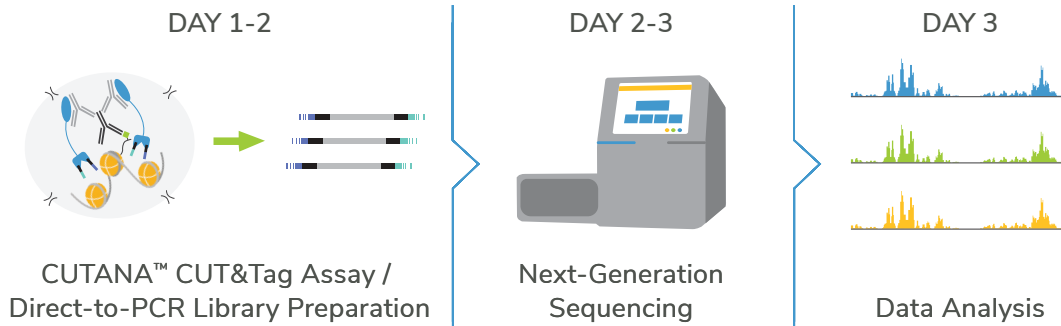


**FIGURE 1** CUT&Tag was performed using an H3K27me3 antibody and decreasing amounts of K562 nuclei as input.

## Advantages of CUTANA CUT&Tag

- Reliable data with ultra-low cell inputs
- Exclusive single-tube workflow
- User-friendly protocol with FAQs and troubleshooting tips
- Defined spike-in controls ensure experimental success

# Go from cells to sequencing in just a few days



## KITS

### Get started with our CUTANA™ CUT&Tag Kit

#### ADVANTAGES:

- Streamlined, single-tube protocol
- Lowest price per reaction vs. competitors
- Includes all the reagents and controls you need for successful CUT&Tag

#### ORDERING INFO:

##### CUT&Tag Kit 48 reactions

Cat. No. 14-1102 - Primer Set 1  
Cat. No. 14-1103 - Primer Set 2

#### CUTANA CUT&Tag Kit



## RESOURCES

### PROTOCOLS & RESOURCES

EpiCypher offers a detailed CUT&Tag protocol and quantitative spike-in controls to support robust histone PTM profiling.

CUT&Tag Protocol: [epicypher.com/protocols](https://epicypher.com/protocols)

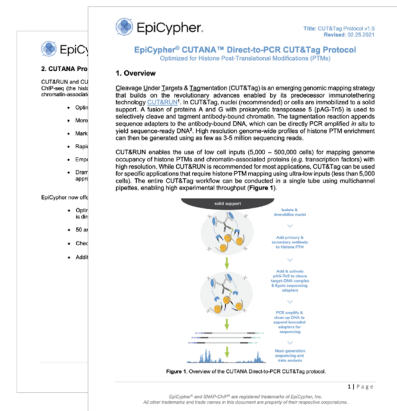
SNAP-CUTANA™ Spike-in User Guide: [epicypher.com/protocols](https://epicypher.com/protocols)

CUT&Tag vs. CUT&RUN Video: <https://youtu.be/90hD69eQ41g>

#### BLOGS

Visit [epicypher.com/blog](https://epicypher.com/blog) for information and guidance:

- The Complete Guide to CUT&Tag Experiments
- ChIP-seq vs. CUT&RUN vs. CUT&Tag: Which should you use?
- Starting CUT&RUN or CUT&Tag for a new target: What you need to know



## PRODUCTS

### ENZYMES & REAGENTS

pAG-Tn5  
50 / 250 reactions  
Cat. No. 15-1017  
Cat. No. 15-1117

ConA Conjugated Paramagnetic Beads  
50 / 250 reactions  
Cat. No. 21-1401  
Cat. No. 21-1411

Non-Hot Start 2X PCR Master Mix  
50 reactions  
Cat. No. 15-1018

### PRIMARY ANTIBODIES

H3K27me3 Antibody  
Cat. No. 13-0055

H3K4me1 Antibody  
Cat. No. 13-0057

Rabbit IgG Negative Control  
Cat. No. 13-0042

### SPIKE-IN CONTROLS

SNAP-CUTANA™ K-MetStat Panel  
Cat. No. 19-1002

### SECONDARY ANTIBODIES

Anti-Rabbit Secondary Antibody  
50 / 250 reactions  
Cat. No. 13-0047  
Cat. No. 13-1047

Anti-Mouse Secondary Antibody  
50 / 250 reactions  
Cat. No. 13-0048  
Cat. No. 13-1048

### TOOLS

Magnetic Separation Racks  
Cat. No. 10-0008 (0.2 mL)  
Cat. No. 10-0012 (1.5 mL)



### Let's discuss your project

[info@epicypher.com](mailto:info@epicypher.com)

855.374.2461

[epicypher.com](https://epicypher.com)