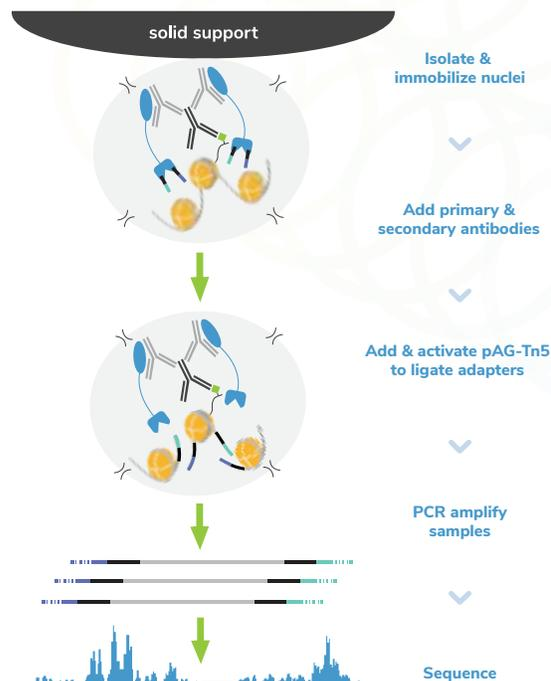


CUTANA™ CUT&Tag Assays for chromatin mapping with low cell numbers

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

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.

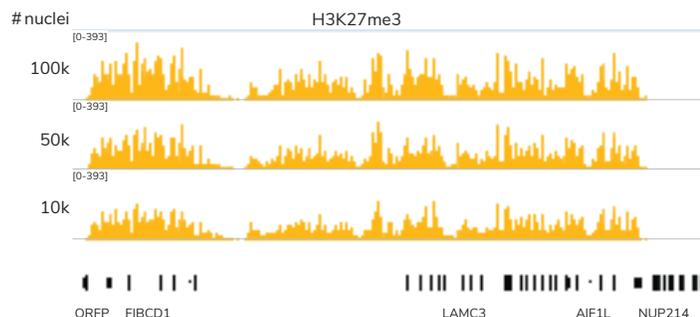
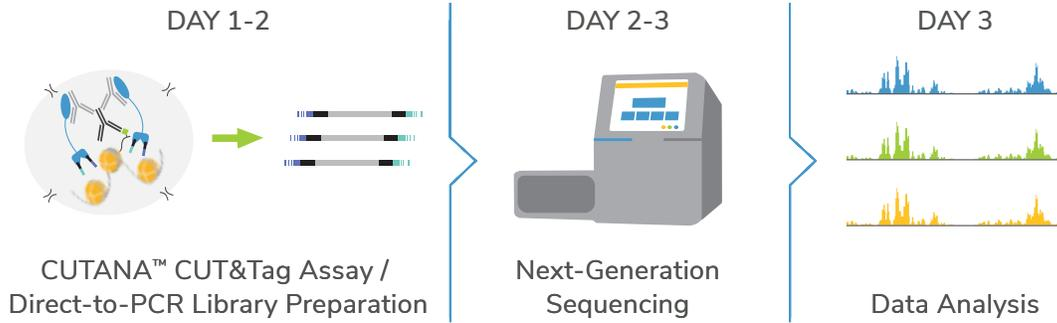


FIGURE 1 CUT&Tag generates highly reproducible H3K27me3 profiles down to 10,000 nuclei. K562 cell were used as input.

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

SNAP-CUTANA™ Spike-in User Guide: epicypher.com/protocols

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

BLOGS

Visit 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

H3K4me1 Antibody
Cat. No. 13-0057

H3K4me3 Antibody
Cat. No. 13-0060

H3K27ac Antibody
Cat. No. 13-0059

H3K27me3 Antibody
Cat. No. 13-0055

Rabbit IgG Negative Control
Cat. No. 13-0042

SECONDARY ANTIBODIES

Anti-Rabbit Secondary Antibody
Cat. No. 13-0047

Anti-Mouse Secondary Antibody
Cat. No. 13-0048

SPIKE-IN CONTROLS

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

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

855.374.2461

epicypher.com