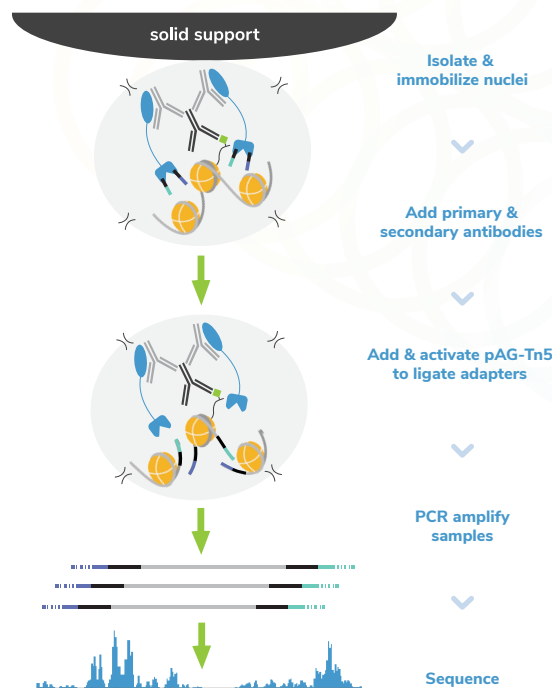


CUTANA™ CUT&Tag Assays for chromatin mapping with low cell numbers

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
- High signal-to-noise and low background
- Fewer cells needed
- Rapid 2-day workflow
- Reduced sequencing costs



Reliable data at low cell numbers

FEATURES	ChIP-seq	CUT&RUN	CUT&Tag
Cells Required	>1 million	5,000 - 500,000	1,000 - 100,000
Compatible Targets	Histone PTMs, TFs	Histone PTMs, TFs & Chromatin Remodelers	Histone PTMs & Select TFs
Sequencing Depth (Reads)	>30 Million	3-8 Million	3-8 Million
Signal-to-Noise	Low	High	High

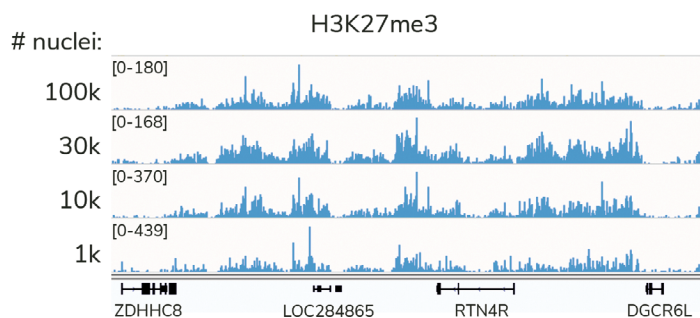
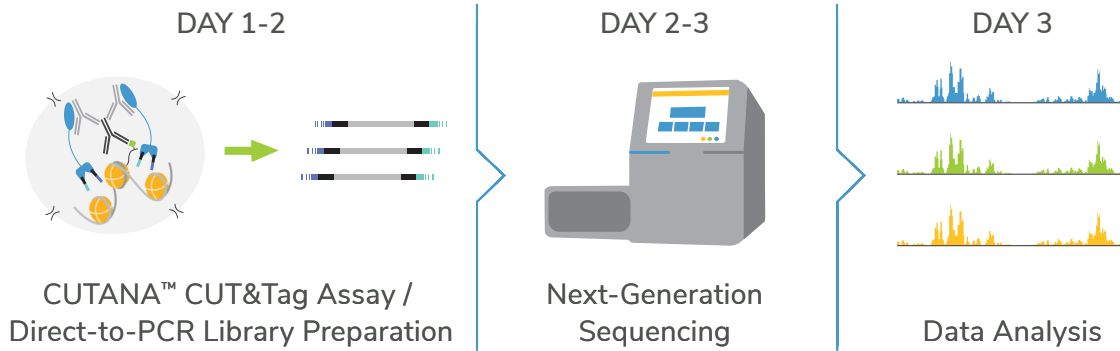


FIGURE 1 CUT&Tag was performed using an H3K27me3 antibody and decreasing amounts of K562 nuclei as input.

Advantages of CUTANA™ CUT&Tag

- Reliable data down to 1,000 cells
- Exclusive single-tube workflow
- User-friendly protocol with FAQs and troubleshooting tips
- Spike-in controls available to maximize experimental success

Go from cells to sequencing in just a few days



PROTOCOL

Learn more about EpiCypher's optimized CUTANA™ Direct-to-PCR CUT&Tag Protocol.

[CUTANA CUT&Tag Protocol](#)

ADVANTAGES

- Cells to DNA in one tube
- Use as few as 1,000 cells
- Includes modifications for cells, nuclei, cross-linking, fresh and frozen samples
- Perform assay in <2 days
- Recommended for histone PTMs
- For chromatin-associated protein targets, contact info@epicypher.com for recommendations



PRODUCTS

Get started with our CUT&Tag Core Reagents Bundle

50 reactions / Cat. No. 14-1101

Includes the essential reagents and controls you need for CUT&Tag assays!

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

ANTIBODIES

Rabbit IgG Negative Control
Cat. No. 13-0042

H3K4me3
Cat. No. 13-0041

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

SPIKE-IN CONTROLS

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

TOOLS

8-strip 0.2 mL PCR Tubes
120 strips
Cat. No. 10-0009

0.2 mL Magnetic Separation Rack
Cat. No. 10-0008

1.5 mL Magnetic Separation Rack
Cat. No. 10-0012

RESOURCES

PROTOCOLS

Visit epicypher.com/protocols for:

- CUT&Tag protocol
- SNAP-CUTANA™ Spike-in User Guide
- Cross-linking protocol
- Protocols are routinely updated!

LITERATURE

- Single cell analysis - Kaya-Okur et al. Nat. Comm. 2019 (PMID : 31036827)
- Automated CUT&Tag - Janssens et al. Nat. Genet. 2021 (PMID : 34663924)
- Chromatin accessibility using CUT&Tag - Henikoff et al. eLife. 2020 (PMID : 33191916)

BLOGS

Check out our recent CUT&Tag blogs at epicypher.com/resources/blog

CONTACT US - Let's talk about your project

info@epicypher.com

855.374.2461

epicypher.com



EpiCypher
Bringing Epigenetics to Life