

CUTANA™ Antibodies for CUT&RUN and CUT&Tag assays

As the leading authority in antibody validation for chromatin mapping assays, EpiCypher knows good antibodies are hard to find. That's why each CUTANA antibody is meticulously validated in CUT&RUN and/or CUT&Tag, ensuring successful on-target profiling.

Histone PTM antibodies

Many commercial PTM antibodies cross-react, show poor target recovery, and generate unreliable data. ONLY EpiCypher can accurately confirm antibody performance in CUT&RUN and/or CUT&Tag.

Metrics include:

- Proven target specificity using exclusive SNAP Spike-in nucleosomes
- Robust target enrichment at low cell numbers
- Consistent results in lot-to-lot testing

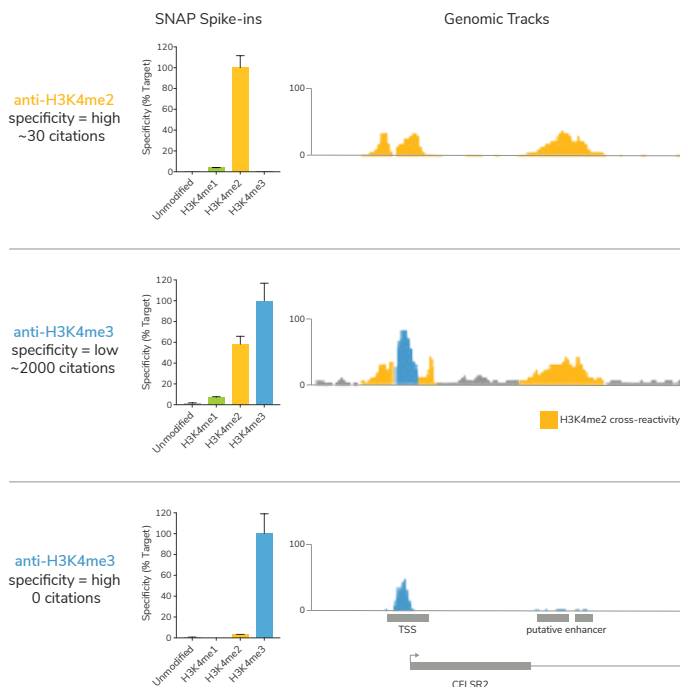
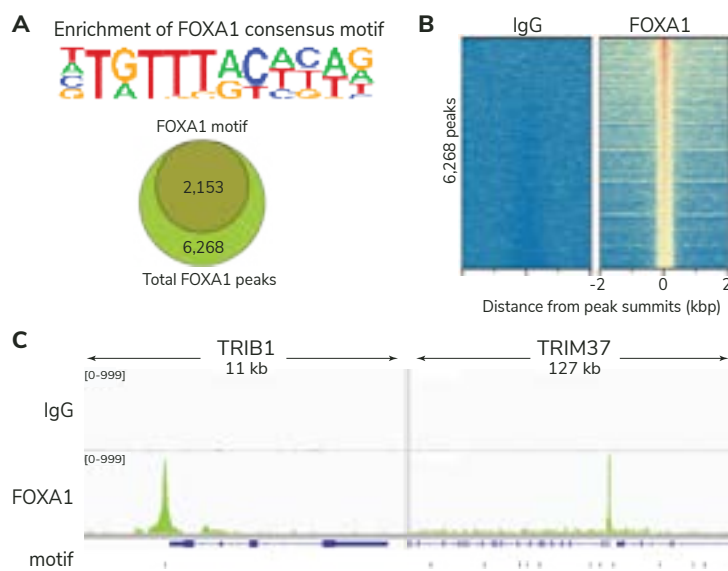


FIGURE 1

A highly cited H3K4me3 antibody (middle panel) fails SNAP Spike-in testing and shows extensive cross-reactivity to H3K4me2. Highly specific H3K4me2 (top panel; orange) and H3K4me3 (bottom panel; blue) antibodies are shown for comparison. See Shah et al. 2018.

Chromatin-associated protein antibodies



CUT&RUN for transcription factors and other proteins rely on high-quality antibodies. Our unique validation strategy includes multiple genome-wide analyses, so you can be confident in your data.

We confirm:

- High signal over background
- Peak structures and motif enrichment consistent with biological function
- Reliable performance in lot-to-lot testing

FIGURE 2

EpiCypher antibody to FOXA1 (Cat. No. 13-2001) enriches for the FOXA1 consensus motif in CUT&RUN assays using K562 cells (A,C) with high signal-to-noise as shown by heatmap analysis (B).

Find the best antibody for your experiment



CUTANA™ CUT&RUN Antibodies

Histone PTMs (SNAP-Certified)

| | |
|----------|---------|
| H3K4me1 | 13-0057 |
| H3K4me2 | 13-0027 |
| H3K4me3 | 13-0060 |
| H3K27ac | 13-0059 |
| H3K27me3 | 13-0055 |
| H3K36me3 | 13-0058 |
| H4K20me3 | 13-0054 |

Chromatin Modifying Enzymes & Regulators

| | |
|------------|---------|
| BRD4 | 13-2003 |
| EZH2 | 13-2026 |
| Menin | 13-2021 |
| MLL1/KMT2A | 13-2004 |

Signaling Receptors

| | |
|------------------|---------|
| AR | 13-2020 |
| EGFR | 13-2018 |
| ERα (C-terminal) | 13-2012 |
| ERα (N-terminal) | 13-2011 |
| NCOA3/SRC3 | 13-2013 |

Chromatin Remodeling Enzymes

| | |
|---------------|---------|
| BRG1/SMARCA4 | 13-2002 |
| BRM/SMARCA2 | 13-2006 |
| CHD1 | 13-2008 |
| CHD3 | 13-2009 |
| CHD4 | 13-2016 |
| SNF2H/SMARCA5 | 13-2007 |
| SNF2L/SMARCA1 | 13-2005 |

Transcription Factors & Regulators

| | |
|-------------|---------|
| CTCF | 13-2014 |
| ELF1 | 13-2023 |
| FOXA1/HNF3A | 13-2001 |
| JUN/c-Jun | 13-2019 |
| SP1 | 13-2024 |
| TP53/p53 | 13-2015 |

Controls & Epitope Tags

| | |
|----------------------|---------|
| IgG Negative Control | 13-0042 |
| DYKDDDDK Tag | 13-2031 |
| HA Tag | 13-2010 |



CUTANA™ CUT&Tag Antibodies

Histone PTMs (SNAP-Certified)

| | |
|----------|---------|
| H3K4me1 | 13-0057 |
| H3K4me3 | 13-0060 |
| H3K27ac | 13-0059 |
| H3K27me3 | 13-0055 |

Controls & Secondary Antibodies

| | |
|--------------------------------|---------|
| IgG Negative Control | 13-0042 |
| Anti-Mouse Secondary Antibody | 13-0048 |
| Anti-Rabbit Secondary Antibody | 13-0047 |



ChIP Antibodies

Histone PTMs (SNAP-Certified)

| | | | |
|-----------|---------|---------|---------|
| H3K4ac | 13-0034 | H3K36ac | 13-0035 |
| H3K18acyl | 13-0050 | H4K12ac | 13-0037 |



EpiCypher®

Don't see your target?
Let us know!

info@epicypher.com

855.374.2461

epicypher.com

[CUTANA CUT&RUN Antibodies](http://epicypher.com/cut-and-run-antibodies)
epicypher.com/cut-and-run-antibodies

[CUTANA CUT&Tag Antibodies](http://epicypher.com/cut-and-tag-antibodies)
epicypher.com/cut-and-tag-antibodies

[SNAP-CUTANA Spike-in Controls](http://epicypher.com/snap-spike-ins)
epicypher.com/snap-spike-ins

[CUTANA CUT&RUN and CUT&Tag Assays](http://epicypher.com/cutana-assays)
epicypher.com/cutana-assays