

CUTANA™ CUT&Tag Core Reagents Bundle



EpiCypher®

Catalog No 14-1101
Lot No 23026004-02
Secondary Ab Anti-Rabbit
Pack Size 50 Reactions

Product Description:

The CUTANA™ CUT&Tag Core Reagents Bundle brings together the essential components of a CUT&Tag kit, offering a significant cost savings over purchasing each individual reagent. Detailed instructions for using the bundled reagents can be found in the EpiCypher CUT&Tag protocol (epicypher.com/protocols). Included are ConA magnetic beads for immobilizing nuclei or cells, control antibodies, anti-rabbit secondary antibody, spike-in controls, pAG-Tn5 for antibody-targeted chromatin tagmentation, and PCR master mix for library amplification. The positive and negative control antibodies (H3K4me3, IgG) can be paired with the SNAP-CUTANA™ K-MetStat Panel of spike-in controls for assay optimization and reaction monitoring.

Some necessary reagents including CUT&Tag assay buffers, primers, and DNA purification beads are not included in the bundle (see Required Materials). For detailed buffer recipes, multiplexing primer sequences and recommended vendors, see the CUT&Tag protocol.

The EpiCypher CUT&Tag protocol is compatible with a variety of inputs, though nuclei are preferred to reduce mitochondrial background. While it is recommended to start with 100,000 cells, comparable data can be generated using as few as 1,000 cells. CUT&Tag provides robust results when profiling histone post-translational modifications (PTMs), but is not recommended for chromatin associated proteins (CAPs). The inclusion of controls, low cell tolerance, and robust PTM profiling potential make the CUT&Tag bundle the ideal choice for experienced users looking for high throughput chromatin mapping. For new users, projects mapping CAPs, or projects mapping histone PTMs where >5,000 cells / sample can be attained, check out the CUTANA™ ChIC/CUT&RUN Kit (EpiCypher 14-1048) for an all-inclusive and robust epigenomic profiling solution.

Bundle Contents:

<u>Item</u>	<u>Pack Size</u>	<u>CAT</u>
H3K4me3 Ctrl Ab	10 Reactions	13-0041t
Rabbit IgG Ctrl Ab	10 Reactions	13-0042t
Anti-rabbit secondary	50 Reactions	13-0047
pAG-Tn5	50 Reactions	15-1017
PCR master mix	50 Reactions	15-1018
SNAP-CUTANA K-Met	20 Reactions	19-1002t
ConA Beads	50 Reactions	21-1401

Required Materials Not Supplied:

The following reagents are not included. See the CUT&Tag Protocol (epicypher.com/protocols) for more information:

- Assay buffers
- Multiplexing primers (requires custom oligo synthesis)
- SPRI Beads (e.g. AMPure XP, SPRIselect)
- 8-strip tubes & magnets
- Qubit™ 4 Fluorometer & 1X dsDNA HS Kit
- Agilent BioAnalyzer® or TapeStation®
- **If using a mouse primary antibody**, the anti-mouse secondary antibody (EpiCypher 13-0048) must be purchased separately.

Storage and Stability:

DO NOT FREEZE ConA BEADS. Upon receipt, remove Concanavalin A (ConA) Beads and store at 4°C. Store the remainder of the components at -20°C. Stability from date of receipt as indicated on each reagent.

Instructions for use:

Download the CUT&Tag protocol for additional required buffers, materials, and equipment, please see epicypher.com/protocols. Use 0.5 µg per CUT&Tag reaction for each control antibody.

This product is for *in vitro* research use only and is not intended for use in humans or animals.