

CUTANA™ DNA Purification Kit

Catalog No. 14-0050
Lot No. 21131003-02
Pack Size 50 samples



EpiCypher®

Product Description:

Cleavage Under Targets & Release Using Nuclease (CUT&RUN) utilizes an immunotethering approach to specifically release genomic fragments of interest into solution, enabling next-generation sequencing and genomic mapping with unprecedented sensitivity. High yield purification and concentration of partitioned chromatin is an essential step; one that requires a workflow specifically compatible with low starting material. Utilizing tapered cleanup columns designed for low elution volumes (as low as 6 µL), this kit produces purified, concentrated DNA ready for library preparation and next-generation sequencing (**Figure 1**). The user-friendly workflow can be completed in just 30 minutes. The columns recover > 50 bp DNA fragments, making the DNA Purification Kit ideal for a variety of CUT&RUN targets, including histone PTMs and transcription factors.

Kit Contents:

50 DNA Cleanup Columns + Collection Tubes
DNA Binding Buffer
DNA Wash Buffer - Add 20 mL ≥ 95% EtOH prior to use
DNA Elution Buffer

Storage and Stability:

Store at room temperature. Stable for 6 months from date of shipment.

Instructions for use:

See kit manual for complete instructions.

References:

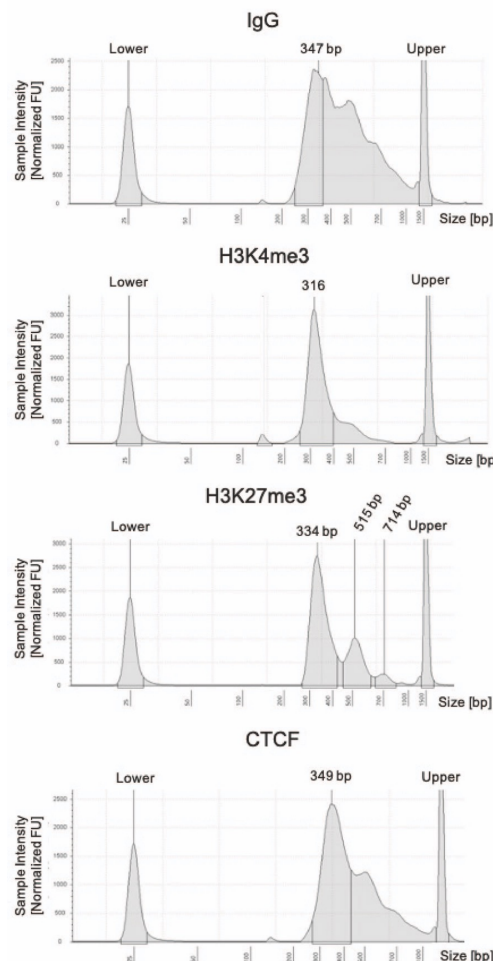


Figure 1: CUT&RUN DNA Fragment Size Distribution Analysis. 500,000 K562 cells per sample were processed with the CUTANA ChIC/CUT&RUN Kit following treatment with IgG negative control (EpiCypher Catalog No. 13-0042), H3K4me3 (EpiCypher Catalog No. 13-0041), H3K27me3 (EpiCypher Catalog No. 13-0030) and CTCF (EMD Millipore Catalog No. 07-729) antibodies. Isolated DNA was used to prepare paired-end Illumina® sequencing adaptors, and libraries were analyzed by Agilent TapeStation®, which confirmed enrichment of mononucleosomes (~300 bp peak represents 150 bp nucleosomes + sequencing adaptors).

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

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

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com