

# CUTANA™ Concanavalin A-Conjugated Paramagnetic Beads



EpiCypher®

**Catalog No.** 21-1401

**Lot No.** 21131006-03

**Pack Size** 550  $\mu$ L

## Product Description:

This product contains Concanavalin A (ConA) conjugated to paramagnetic microspheres. ConA is a lectin (carbohydrate-binding protein) that binds specifically to mannosyl- and glucosyl-containing extracellular glycoproteins. The ConA magnetic beads are therefore useful to immobilize cells or nuclei presenting these glycans in their extracellular matrices.

## Formulation:

Concanavalin-A conjugated to 1  $\mu$ m paramagnetic microspheres in 10 mM PBS with 0.1% sodium azide.

## Storage and Stability:

DO NOT FREEZE!! Stable for six months at 4°C from date of receipt.

## Application Notes:

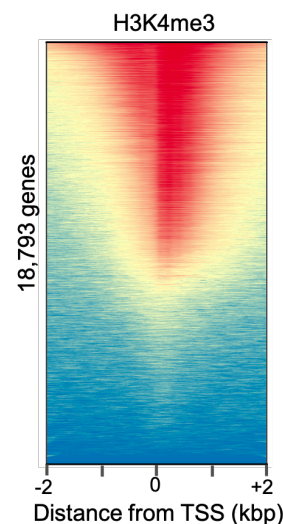
For a detailed CUT&RUN protocol using this product, see the EpiCypher CUT&RUN Protocol at:

[www.epicypher.com/resources/protocols](http://www.epicypher.com/resources/protocols).

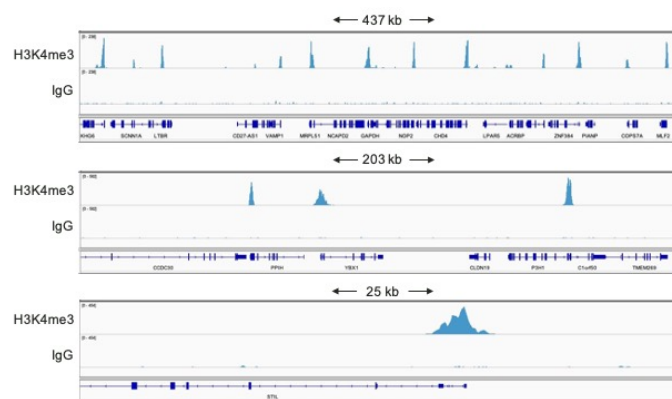
\*Note: It is recommended to use 10  $\mu$ L slurry per sample (500,000 cells or less) for ChIC/CUT&RUN. Compatible with EpiCypher Magnetic Separation Racks (Catalog Nos. 10-0008 & 10-0012).

## References:

Lai et al., *bioRxiv* (2020.06.08.140046)



**Figure 1.** CUT&RUN data was generated using EpiCypher CUTANA Concanavalin A-Conjugated Paramagnetic Beads (10  $\mu$ L) with 500,000 K562 nuclei and H3K4me3 antibody (EpiCypher 13-0041, 0.5  $\mu$ g). Shown is a heatmap for H3K4me3 CUT&RUN signal (from 7.4 million paired-end reads) aligned to the transcription start site (TSS, +/- 2kb) of 18,793 genes. High and low signal are ranked by intensity (top to bottom) and reflected by red and blue colors, respectively.



**Figure 2.** CUT&RUN data was generated as described above. Three representative loci are shown in the Integrative Genomics Viewer (IGV, Broad Institute). Window size is denoted at the top of each panel. Top tracks at each locus were generated using H3K4me3 antibody (EpiCypher Catalog No. 13-0041, 0.5  $\mu$ g), while the bottom tracks were generated using IgG Negative control antibody (EpiCypher Catalog No 13-0042, 0.5  $\mu$ g).

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