

Histone H3 C-terminal Antibody



EpiCypher®

Catalog No 13-0001

Lot No 21264002-01

Pack Size 100 µL

Type Polyclonal

Host Rabbit

Target Size 15 kDa

Format Serum

Reactivity *C. elegans*, Chicken, *Drosophila*, Human, *S. cerevesiae*, *S. pombe*, Wide Range (predicted)

Applications WB

Product Description:

Rabbit polyclonal antibody recognizing the C-terminus of histone H3, for use in Western blot. Histone H3 is one of a core components of the nucleosome, the basic unit of chromatin.

Immunogen:

Synthetic peptide corresponding to the C-terminus of human histone H3.1, conjugated to KLH.

Formulation:

Rabbit serum with 30% glycerol and 0.035% sodium azide.

Storage and Stability:

Stable for 2 years at -20°C from date of receipt.

Recommended Dilution:

This antibody has been validated for use in Western blotting (1:2,500 - 1:5,000). This antibody serves as a good loading control for Western Blot analysis or negative control for experiments involving H3 modifications.

References:

Rizzardi LF et al (2012) *Genetics* 192:371-384.

Rothbart SB et al (2012) *Nat Struct Mol Biol* 11:1155-1160.



Western Blot Data: Western blot using Histone H3 C-terminal Antibody (1:5,000 dilution). HEK293 core histones (0.2 µg) were blotted and probed.

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