

# Histone H3 C-terminal Antibody

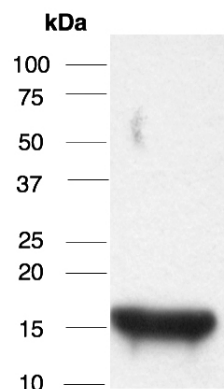
**Catalog No.** 13-0001  
**Lot No.** 12346001  
**Pack Size** 100 µl



**Type** Polyclonal      **Host** Rabbit  
**Mol. Wgt.** 15 kDa      **Reactivity** Ce, Ch, Dm, H, Sc,  
**Format** Serum      **Appl.** ChIP, WB

## Product Description:

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



**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.

## Immunogen:

Synthetic peptide corresponding the 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:

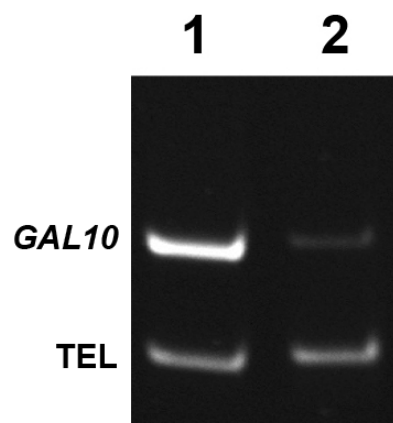
Stable for 2 years at -20°C from date of receipt.

## Application Notes:

This antibody has been validated for use in ChIP (3-5µl per reaction) and Western blotting (1:2,500 - 1:5,000). This antibody serves as a good negative control for

## References Using this Product:

Rizzardi LF et al (2012). *Genetics* 192: 371-384.  
Rothbart SB et al (2012). *Nat Struct Mol Biol* 11: 1155-1160.  
Rothbart SB et al (2013). *Genes Dev* 27: 1288-1298.



**ChIP Data:** ChIP using Histone H3 C-terminal Antibody (4.5 µl per ChIP) with *S. cerevesiae* chromatin. PCR was performed using *GAL10* gene primers or telomeric (TEL) control primers. Lane 1: Chromatin taken from cells where *GAL10* gene is not expressed. Lane 2: *GAL10* gene expressed.

**Applications Key:** ChIP: Chromatin IP; ChIP-seq: Chromatin IP sequencing; E: ELISA; FACS: Flow cytometry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; WB: Western Blotting

**Reactivity Key:** B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S.cerevesiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish