

H3R8me2(sym) polyclonal antibody - Classic

Cat. No. C15410287

Type: Polyclonal	Specificity: Human, mouse, C. elegans, rat, chicken, Xenopus, Drosophila, plant
Size: 50 µg	Isotype: NA
Concentration: 1.1 µg/µl	Source: Rabbit
Lot No.: 001	Purity: Affinity purified
Storage buffer: NA	Storage conditions: NA
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Description

Polyclonal antibody raised in rabbit against H3 (sym-dimethyl Arg8), using a KLH-conjugated synthetic peptide.

Applications

Applications	Suggested dilution	References
ChIP	2-5 µg/million cells	Figure 1
Western Blotting	2 µg/mL	Figure 2, 3

Target Description

Chromatin is the arrangement of DNA and proteins in which chromosomes are formed. Correspondingly, chromatin is formed from nucleosomes, which are comprised of a set of four histone proteins (H2A, H2B, H3, H4) wrapped with DNA. Chromatin is a very dynamic structure in which numerous post-translational modifications work together to activate or repress the availability of DNA to be copied, transcribed, or repaired. These marks decide which DNA will be open and commonly active (euchromatin) or tightly wound to prevent access and activation (heterochromatin). Common histone modifications include methylation of lysine and arginine, acetylation of lysine, phosphorylation of threonine and serine, and sumoylation, biotinylation, and ubiquitylation of lysine. Specifically, methylation of arginine 8 on histone H3 (H3 R8me2s) is associated with transcriptional repression, and modified by PRMT5, but not CARM1.

Validation Data

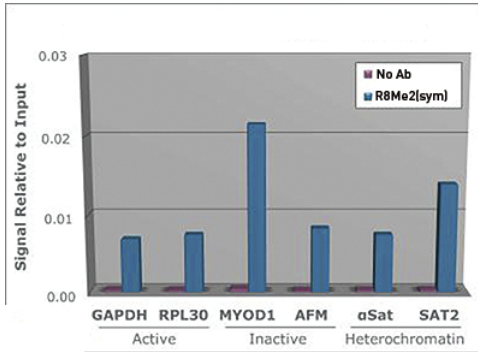


Figure 1. H3R8me2(sym) antibody ChIP results

Chromatin Immunoprecipitation using the H3R8me2(sym) antibody. Chromatin from one million formaldehyde cross-linked HeLa cells was used with 2 µg of H3R8me2(sym) to IP DNA from fixed HeLa cells alongside a no antibody (No Ab) control. DNA was measured by PCR and normalized to total input.

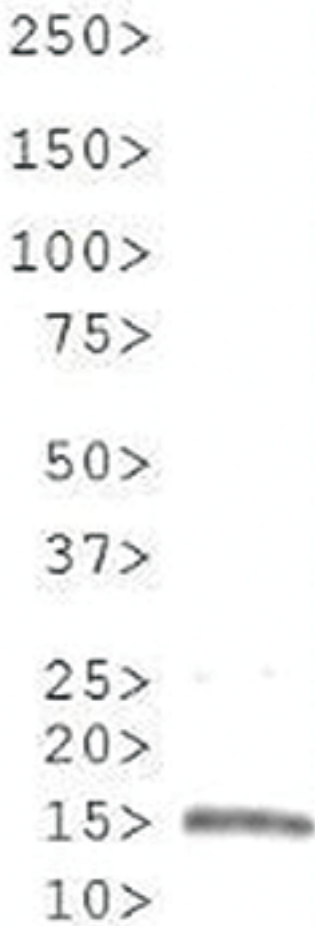


Figure 2 H3R8me2(sym) antibody western blot results

Western Blot of H3R8me2(sym) antibody. 30 µg C. elegans embryo lysate. Primary antibody: diluted 1:500 overnight at 4°C. Secondary antibody: IRDye800TM rabbit secondary antibody at 1:10,000 for 45 min at RT. Predicted/Observed size: ~15 kDa. Other band(s): None.

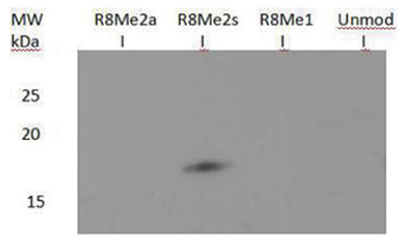


Figure 3. H3R8me2(sym) antibody Western blot results

Western Blot of H3R8me2(sym) antibody. Lane 1: R8Me2a. Lane 2: R8Me2s. Lane 3: R8Me1. Lane 4: Histone H3 R8. Load: 30 µg per lane. Primary antibody incubated at 2 µg/mL overnight at 4°C. Secondary antibody IRDye800TM rabbit secondary antibody at 1:10,000 for 45 min at RT. Predicted/Observed size: ~15 kDa. Other band(s): None.