

## H3T6pK9me2 polyclonal antibody - Classic

**Cat. No.** C15410284

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 001

**Size:** 50 µg

**Concentration:** 0.49 µg/µl

**Specificity:** Human, mouse, C. elegans, rat, chicken, Xenopus, Drosophila, plant

**Purity:** Affinity purified

**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

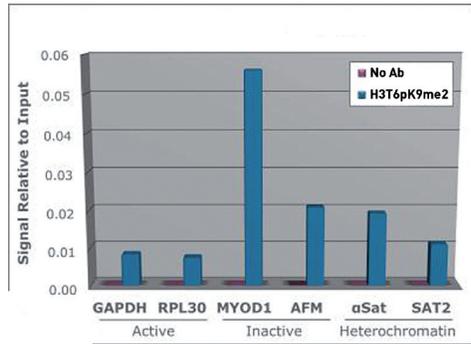
### Applications

	Suggested dilution	Results
ChIP	2-5 µg/million cells	Figure 1
Immunohistochemistry	1:100	
IF	1:100	
Western blot	1:500	

### Target description

The doubly modified dimethyl K9, phospho-T6 histone H3 is known to exist, but very little comprehensive information about the importance of this modification and its mechanism has been published. H3K9me2 is a histone post-translational modification that is enriched in transcriptionally inactive regions. It is also important in cellular differentiation and maturation. Phosphorylation of histone H3 is also associated with mitosis and meiosis, and seems to be related to developmental changes. The dual modification of H3T6pK9me2 is still under investigation and will present more insight into the complex epigenetic code.

## Results



**Figure 1. H3T6pK9me2 antibody ChIP results**

Chromatin Immunoprecipitation of H3T6pK9me2 antibody. Chromatin from one million formaldehyde cross-linked HeLa cells was used with 2 µg of H3T6pK9me2 antibody to IP DNA from fixed HeLa cells alongside a no antibody (No Ab) control. DNA was measured by qPCR and normalized to total input.

### Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

### Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

Last update: September 10, 2014