

## H3S10pT11p polyclonal antibody - Classic

**Cat. No.** C15410288

**Type:** Polyclonal

**Source:** Rabbit

**Lot #:** 001

**Size:** 50 µg

**Concentration:** 1.2 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.

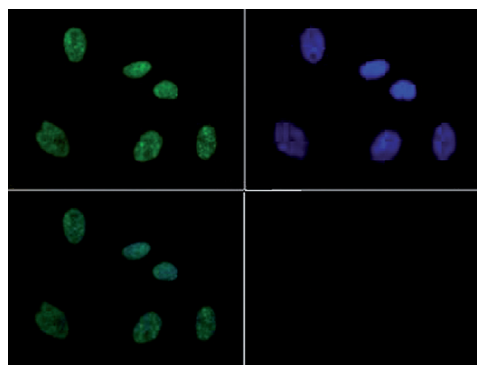
### Applications

	Suggested dilution	Results
ChIP	2-5 µg/million cells	
Immunohistochemistry	1:200	
IF	1:200	Figure 1
Western blot	1:500	Figure 2, 3

### Target description

When confronted with DNA damage or the lack of certain kinases like Chk1 and Aurora1, histone H3 is phosphorylated at T11, causing a reduction in acetylation of the same histone at K9 and ultimately leading to a reduction of active transcription in affected cells. When the T11 phosphorylation is combined with a typical S10 phosphorylation that is critical in chromosome condensation in metaphase, the effect on transcriptional regulation is amplified further. Phosphorylation of H3S10 is decreased by T11 phosphorylation. Thus, detection of the combined modification is critical in determining transcriptional effects, especially when contrasted against single modifications of H3.

## Results



**Figure 1. H3S10pT11p Immunofluorescence results**

Immunofluorescence of H3S10pT11p antibody. Tissue: HeLa cells. Fixation: 0.5% PFA. Primary antibody: H3S10pT11p antibody at a 1:200 dilution for 1 h at RT. Secondary antibody: FITC secondary antibody at 1:10,000 for 45 min at RT. Localization: H3S10pT11p is nuclear and chromosomal. Staining: H3S10pT11p is expressed in green, nuclei are counterstained with Dapi (blue).

250>  
150>  
100>  
75>  
50>  
37>  
25>  
20>  
15>  
10>

**Figure 2 H3S10pT11p Western blot results**

Western Blot of H3S10pT11p antibody. 30 µg C. elegans embryo lysate. Primary antibody: H3S10pT11p at a 1:500 dilution overnight at 4°C. Secondary antibody: IRDye800™ rabbit secondary antibody at 1:10,000 for 45 min at RT. Predicted/Observed size: ~15 kDa. Other band(s): None.

188 →  
98 →  
62 →  
49 →  
38 →  
28 →  
17 →  
14 →

← H3S10T11p

**Figure 3. H3S10pT11p Western blot results**

Western Blot of H3S10pT11p antibody. 30 µg HeLa histone extracts. Primary antibody: H3S10pT11p at a 1:500 dilution overnight at 4°C. Secondary antibody: IRDye800™ rabbit secondary antibody at 1:10,000 for 45 min at RT. Predicted/Observed size: ~15 kDa. Other band(s): None.

### 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