

## Histone H3 (pSer10) Polyclonal Antibody

(Cat. No. A-7402)

### Background

The primary building block of chromatin is the nucleosome which is made up of DNA wound around eight core histone proteins (two each of H2A, H2B, H3, and H4). Histones play a key role in transcription regulation, DNA repair, DNA replication and chromosomal stability. DNA accessibility is regulated through an intricate series of post-translational modifications of histones, known as the histone code, and nucleosome remodeling. The N-terminal tail of histone H3 protrudes from the globular nucleosome core and can undergo several different types of epigenetic modifications that influence cellular processes. There are numerous well-characterized residues for phosphorylation within the N-terminal tail of histone H3; these include Thr-3, Ser-10, Thr-11, and Ser-28. Phosphorylation of histone H3 on Ser-10 in the N-terminal tail is considered essential during mitosis and meiosis for cell cycle progression and chromosome condensation.

### Description

Rabbit polyclonal antibody to Histone H3 (pSer10)

### Formulation

Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

### Immunogen

Synthetic peptide corresponding to human Histone H3 phosphorylated at Ser10.

### Specificity

Human, Mouse, Rat, Chicken, Drosophila, Monkey, Pig, Rabbit

### Purification

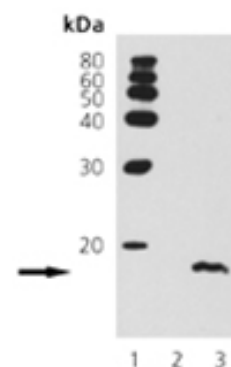
Protein A-affinity purified

### Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

### Application

WB



**Fig. 1.** Western blot analysis of MW marker (1), untreated Jurkat (2), and calyculin A-treated Jurkat (3) cell lysates probed with Histone H3 (pSer10) Polyclonal Antibody.

### Ordering Information

#### Product

Histone H3 (pSer10) Polyclonal Antibody

#### Size

100 µg

#### Cat. No.

A-7402-100

*This product is for research purposes only. Not intended for use in diagnostic procedures.*

Epigentek Group Inc. • USA Tel: 1-877-374-4368 • Int. Tel: 1-718-484-3990 • Fax: 1-718-484-3990 • [info@epigentek.com](mailto:info@epigentek.com) • [www.epigentek.com](http://www.epigentek.com)

2015-10-29-A-7402