

Phospho-Histone H3 (Thr11) Polyclonal Antibody

(Catalog # A-4052)

Background

Histone H3 along with H2A, H2B, and H4 is involved in the structure of chromatin in eukaryotic cells. Histone H3 can undergo several different types of epigenetic modifications that influence cellular processes. These modifications, including acetylation, phosphorylation, methylation, ubiquitination, and ADP-ribosylation, occur on the N-terminal tail domains of histone H3, which results in remodeling of the nucleosome structure into an open conformation more accessible to transcription complexes. In most species, histone H3 is primarily phosphorylated at serine 10, serine 28, threonine 3, and threonine 11.

Description

Phospho-Histone H3 (Thr11) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. PBS, pH 7.4, containing 0.02% sodium azide as Preservative and 50% Glycerol.

Specificity

Mouse, Rat, Human

Isotype

IgG

Uniprot ID

P68431/Q71DI3/P84243

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Storage

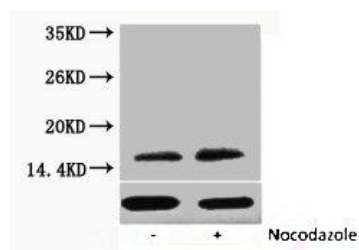
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

HIST1H3A; H3FA; HIST1H3B; H3FL; HIST1H3C; H3FC; HIST1H3D; H3FB; HIST1H3E; H3FD; HIST1H3F; H3FI; HIST1H3G; H3FH; HIST1H3H; H3FK; HIST1H3I; H3FF; HIST1H3J; H3FJ; Histone H3.1; Histone H3/a; Histone H3/b; Histone H3/c; Histone H3/d; Histone H3/f; Histone H3

Application

WB, ELISA; Recommended dilution: WB:1:500-1:1000, ELISA:1:20000



Western blot analysis of extracts from Hela cells, untreated (-) or treated, 1:2000.