
H1 Protein

(Catalog No. E35000)

Description

Histone H1 acts on the linker region of polynucleosome DNA to condense the chromatin into structures of ~30 nm. It is not necessary for octamer or nucleosome core particle formation. Eight different Histone H1 proteins have been identified in the human genome. Histone H10 is a non replication-dependent histone that is highly expressed in cells that have terminally differentiated.

Molecular Weight (kDa)

20.7 kDa

Formulation

20 mM sodium phosphate, 300 mM NaCl, 1 mM EDTA

Source

Expressed in E.coli expression system

Size

50 µg

Storage Temp & Conditions

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

Purification Notes

SDS-PAGE purified

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array

Ordering Information:

Products	Size	Cat. No.
H1 Protein	50 µg	E35000 -1

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