
H4 Protein

(Catalog No. E35005)

Description

Histone H3 combines with Histone H4 to form the H3/H4 tetramer. Two H2A/H2B heterodimers interact with an H3/H4 tetramer to form the histone octamer (1,2). Histone H4 is also modified by various enzymes and can act as a substrate for them. These modifications have been shown to be important in gene regulation.

Molecular Weight (kDa)

11.2 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 at -20°C. Aliquot to avoid repeated freezing and thawing.

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.
H4 Protein	50 µg	E35005 -1

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