
H3F3A Protein

(Catalog No. E35003)

Description

Histone H3.1, an H3 variant that has thus far only been found in mammals, is replication dependent and is associated with gene activation and gene silencing.

Molecular Weight (kDa)

15.3 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.
H3F3A Protein	50 µg	E35003 -1

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