

---

## H3F3b Protein

(Catalog No. E35004)

---

### Description

Histone H3.2, an H3 variant that is found in all eukaryotes except budding yeast, is replication dependent and is associated with 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.
H3F3b Protein	50 µg	E35004 -1

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