

---

## HAT1 Protein

(Catalog No. E27000)

---

### Description

HAT1 Human recombinant protein containing 343 amino acids (20-341 a.a.). With N-terminal His Tag.

### Molecular Weight (kDa)

40.1 kDa

### Formulation

20mM Tris-HCl pH-8, 1mM DTT, and 10% glycerol.

### Source

E. Coli

### Size

10 µg

### Storage Temp & Conditions

6 months at -80°C. Please avoid freeze-thaw cycles.

### Purification Notes

Purity: > 90%

### Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array; Preparation Method: in vitro wheat germ expression system ;Best use within three months from the date of receipt of this protein.

### Ordering Information:

Products	Size	Cat. No.
HAT1 Protein	10 µg	E27000-1

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