

---

## K-Ras Protein

(Catalog No. E60001)

---

### Description

KRAS Recombinant protein is a single, non-glycosylated polypeptide chain containing 205 amino acids (1-185 a.a.) and fused to 20 amino acid His-Tag at N-terminus

### Molecular Weight (kDa)

23.2 kDa

### Formulation

20mM Tris pH-8, 0.1M NaCl, 1mM DTT & 10% glycerol.

### Source

Expressed in E.coli expression system

### Size

25 µg

### Storage Temp & Conditions

KRAS Human although stable at 4°C for 1 week, should be stored desiccated -20°C. Please prevent freeze thaw cycles.

### Purification Notes

Purity :Greater than 90.0% as determined by SDS-PAGE..

### Application Notes

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

### Ordering Information:

Products	Size	Cat. No.
K-Ras Protein	25 µg	E60001 -1

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