
MSK1 Protein

(Catalog No. E22001)

Description

Human RPS6KA5 full-length ORF (AAH17187.1, 1 a.a. - 549 a.a.) recombinant protein with GST-tag at N-terminal.

Molecular Weight (kDa)

86.02 kDa

Formulation

50 mM Tris-HCl, 10 mM reduced Glutathione, pH=8.0 in the elution buffer.

Source

In vitro wheat germ expression system

Size

10 µg

Storage Temp & Conditions

Store at -80°C. 50 mM Tris-HCl, 10 mM reduced Glutathione, pH=8.0 in the elution buffer. Best used within three months from the date of receipt.

Purification Notes

Glutathione Sepharose 4 Fast Flow

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
MSK1 Protein	10 µg	E22001 -1

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