

EPHA8 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # E20-35G

Lot # S261-3

Product Description

Recombinant human EPHA8 (565-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_020526](#).

Gene Aliases

EEK, HEK3, KIAA1459

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

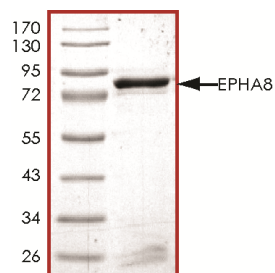
Scientific Background

EPHA8 is a member of the ephrin receptor subfamily of the protein-tyrosine kinase family in which EPH and EPH-related receptors have been implicated in mediating developmental events, particularly in the nervous system. Receptors in the EPH subfamily typically have a single kinase domain and an extracellular region containing a Cys-rich domain and 2 fibronectin type III repeats (1). EPHA8 receptors play a role in axonal pathfinding during development of the mammalian nervous system (2).

References

- Chan, J., et.al: Eek and erk, new members of the eph subclass of receptor protein-tyrosine kinases. *Oncogene* 6: 1057-1061, 1991.
- Park, S. et.al: Aberrant axonal projections in mice lacking EphA8 (Eek) tyrosine protein kinase receptors. *EMBO J.* 16: 3106-3114, 1997.

Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **82 kDa**.

EPHA8 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number **E20-35G**

Specific Lot Number **S261-3**

Purity	>90%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.