

## EPHB4 (I782S), Protein

Human recombinant protein expressed in Sf9 cells

**Catalog # E24-12G**

Lot # H2966-6

### Product Description

Recombinant human EPHB4 (I782S) (561-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_004444](#).

### Gene Aliases

HTK, MYK1, TYRO11, Mdk2

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, and 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.

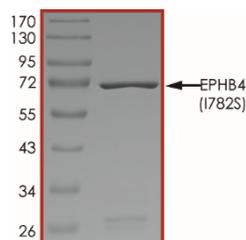
### Scientific Background

EPHB4 is a member of the Eph family of receptor tyrosine kinases which play a role in regulating cell adhesion and cell movement during embryonic development (1). Furthermore, because tissue disorganization and abnormal cell adhesion, movement, and survival characterize the more advanced stages of cancer, the inappropriate functioning of Eph receptor could play a causal role in malignancy (2). EPHB4 is uniquely expressed in the vascular endothelial and endocardial cells and animal knockout studies with EPHB4 reveal a phenotype that is similar to Ephrin B2 gene knockdown. Furthermore, EPHB4 has been shown to be the major essential interaction partner of Ephrin B2 in angiogenesis.

### References

- Andres, A.C. et al: Expression of two novel eph-related receptor protein tyrosine kinases in mammary gland development and carcinogenesis. *Oncogene*, 1994;9(8):2431.
- Kumar, S.R. et al: The receptor tyrosine kinase EphB4 is overexpressed in ovarian cancer, provides survival signals and predicts poor outcome. *Br. J. Cancer*. 2007, 96(7):1083-91.

### Purity



The purity of EPHB4 (I782S) was determined to be **>95%** by densitometry, approx. MW **75 kDa**.

## EPHB4 (I782S), Protein

Human recombinant protein expressed in Sf9 cells

Catalog #	E24-12G
Lot #	H2966-6
Purity	>95%
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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