

## ProQinase™ EPHB1

aa565-984

**Product No.:** 0317-0000-1

**Lot:** 001

**Description:** Human EPHB1  
Amino acids R<sub>565</sub>-A<sub>984</sub> (as in GenBank entry NM\_004441.2)\*, N-terminally fused to GST-HIS<sub>6</sub>-Thrombin cleavage site

\*Sequence may contain documented polymorphisms  
Detailed sequence on request

**Product identity:** EPHB1, Lot 001, was confirmed as human EPHB1 by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

**Theoretical MW<sub>Fusion Protein</sub>:** 76,906 Da

**Expression:** Baculovirus infected Sf9 cells

**Purification:** One-step affinity purification using GSH-agarose

**Storage buffer:** 50 mM Tris-HCl, pH 8.0;  
100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol

**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.500 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Method for determination of K<sub>m</sub> value & Specific activity:**

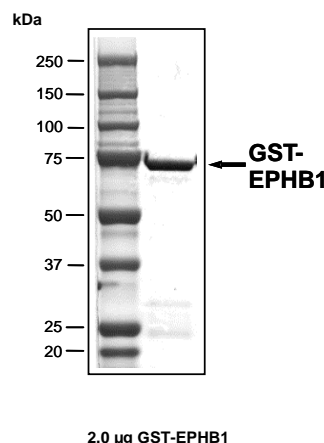
• Assay conditions:

- 60 mM HEPES-NaOH, pH 7.5
- 3 mM MgCl<sub>2</sub>
- 3 mM MnCl<sub>2</sub>
- 3 µM Na-orthovanadate
- 1.2 mM DTT
- 2.5 µg / 50 µl PEG<sub>20.000</sub>
- ATP (variable)
- Substrate: Poly(Glu, Tyr)<sub>4:1</sub>  
(Sigma P-0275), 1 µg / 50 µl
- Recombinant EPHB1: 50 ng / 50 µl

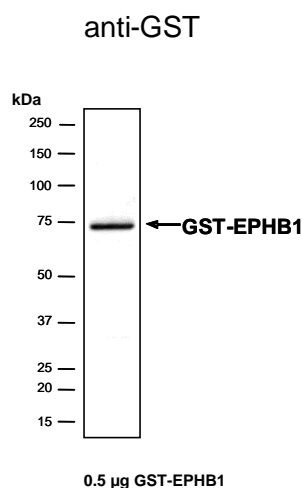
- Filter binding assay  
MAFC membrane (Millipore)

**Specific activity:** 72 pmol/µg×min

**Coomassie stain:**



**Western blot analysis:**



**Determination of K<sub>m</sub> value for ATP:**

