

Data Sheet

EPHB4

Human, recombinant, N-terminal HIS tag

Catalog #: 40201

Lot #: 120716 **Conc.:** 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5,
 150 mM NaCl, 0.1 mM EGTA, 0.1 mM
 EDTA, 0.1 mM PMSF, 50% glycerol and 3
 mM DTT.

Stability: >12 months at -80°C

References:

1. Bruhl, T., *et al.*, *Circ. Res.* **94(6)**, 743-751 (2004.)
2. Fuller, T., *et al.*, *J. Cell Sci.* **116(Pt12)**, 2461-2470 (2003).

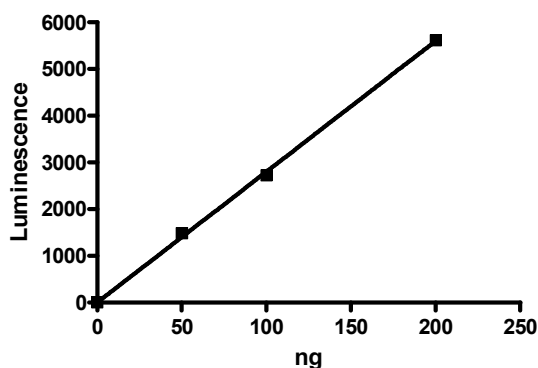
Description: Human EPHB4 (GenBank Accession No. NM_004444), (a.a. 563-987) with N-terminal His tag, MW=52 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 13.2 pmole/min/μg
 Assay conditions: ADP-Glo assay with 0.2 mg/ml poly-(Glu4:Tyr) as a substrate and 100 μM ATP at final concentration. Reaction was done for 50 min at 30°C. ADP-Glo for 45 min at RT, and Kinase detection for 55 min at RT.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

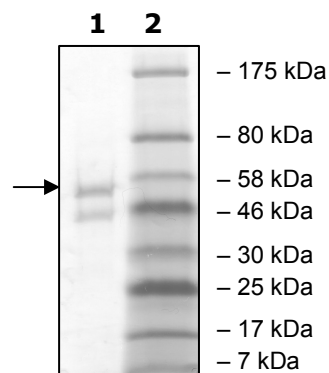
Specific Activity



4-20% SDS-PAGE
Coomassie staining

Lane 1:
 2 μg EPHB4
Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 52 kDa
Purity: ≥60%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com