

Data Sheet

EPHB4

Human, recombinant, N-terminal HIS tag
Catalog #: 40201
Lot #: 120719 **Conc.:** 1.34 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 200 mM
 Imidazole, 3 mM DTT, and 20% glycerol.

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).

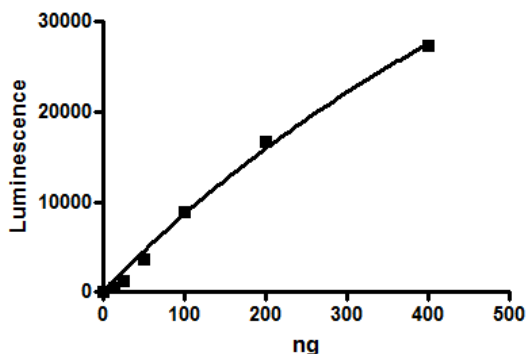
Description: Human EPHB4 (GenBank Accession No. NM_004444), (a.a. 563-987) with N-terminal His tag, MW=52 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 43 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 45 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:
5 μg EPHB4
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 52 kDa
Purity: ≥78%

