

EPHB4

aa563-987

Product No.: 0178-0000-3

Lot: 007

Description: Human EPHB4

Amino acids L₅₆₃-Y₉₈₇ (as in GenBank entry NM_004444.2)*, N-terminally fused to GST-HIS₆-Thrombin cleavage site

*Sequence may contain documented polymorphisms
Detailed sequence on request

Product identity: EPHB4, Lot 007, was confirmed as human EPHB4 by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

Theoretical MW_{Fusion Protein}: 76,794 Da

Expression: Baculovirus infected Sf9 cells

Purification: One-step affinity purification using GSH-agarose

Activation: pre-incubation with 0.1 mM ATP, followed by (NH₄)₂SO₄ precipitation

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

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

Protein concentration: 0.235 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Method for determination of K_m value & Specific activity:

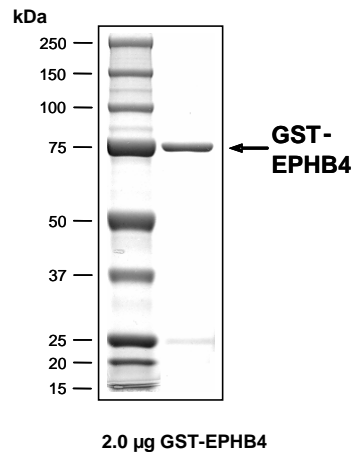
• Assay conditions:

60 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
2.5 µg / 50 µl PEG_{20,000}
ATP (variable)
Substrate: Poly(Glu,Tyr)_{4:1}
(Sigma P-0275), 0,5 µg / 50 µl
Recombinant EPHB4: 25 ng / 50 µl

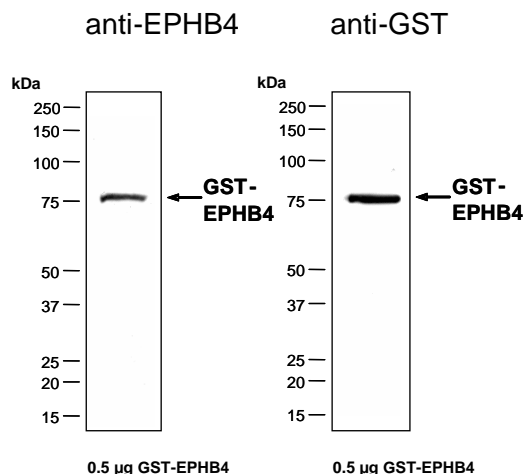
• Filter binding assay
MAFC membrane (Millipore)

Specific activity: 78 pmol/µg×min

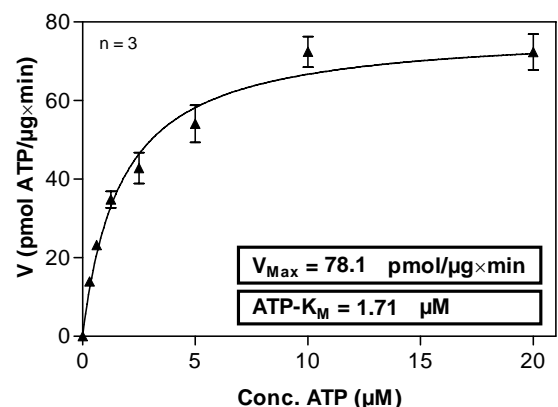
Coomassie stain:



Western blot analysis:



Determination of K_m value for ATP:



This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.