

EPHA1

aa569-976

Product No.: 0215-0000-1

Lot: 002

Description: Human EPHA1

Amino acids R₅₆₉-D₉₇₆ (as in GenBank entry XM_056982)*, N-terminally fused to GST-HIS₆-Thrombin cleavage site

*Sequence may contain documented polymorphisms
Detailed sequence on request

Product identity: EPHA1, Lot 002, was confirmed as human EPHA1 by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

Theoretical MW_{Fusion Protein}: 79,631 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: 0.071 µg/µl (Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Method for determination of K_m value & Specific activity:

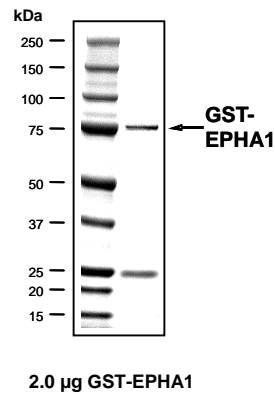
• Assay conditions:

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), 2 µg / 50 µl
Recombinant EPHA1: 50 ng / 50 µl

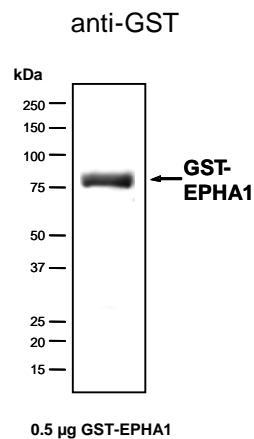
• Filter binding assay
MAFC membrane (Millipore)

Specific activity: 4 pmol/µg×min

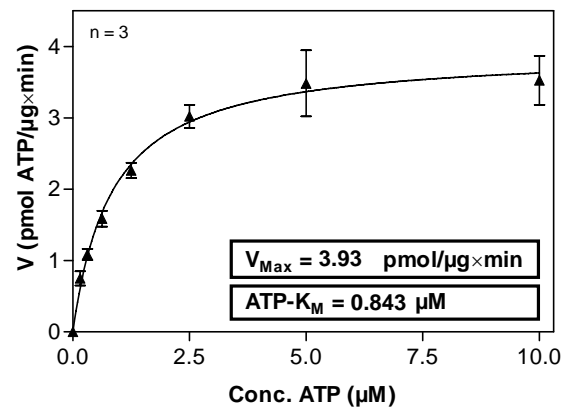
Coomassie stain:



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.