

ProQinase™ 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:

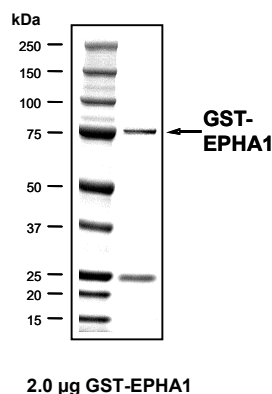
• Assay conditions:

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

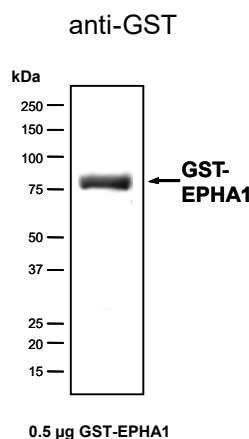
• Filter binding assay
MAFC membrane (Millipore)

Specific activity: 4 pmol/µg×min

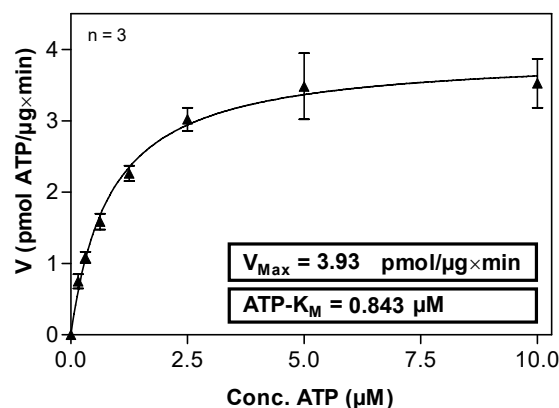
Coomassie stain:



Western blot analysis:



Determination of K_m value for ATP:



Recombinant Proteins