

Product Information

EPHB4

Product Number : **08-131**

Product description

Human EPHB4, cytoplasmic domain [577-987(end) amino acids of accession number NP_004435.3] was expressed as N-terminal GST-protein (73 kDa) using baculovirus expression system. GST-EPHB4 was purified by using glutathione sepharose chromatography.

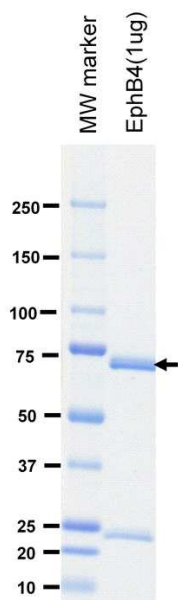
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

Store at -80C.
Avoid repeating freeze-thaws.

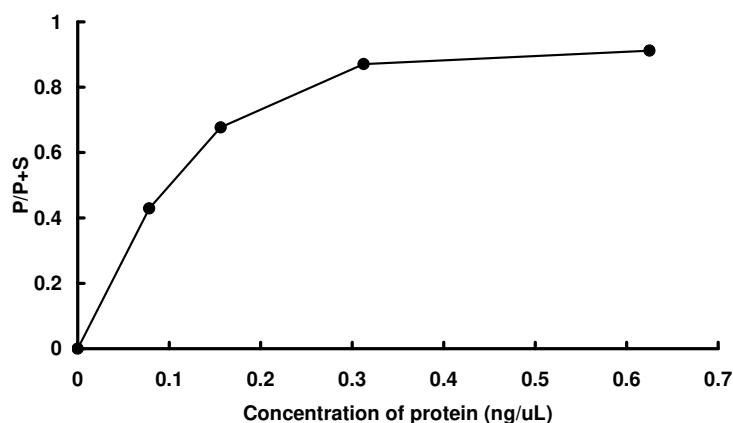
SDS-PAGE



Purity: 70 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Blk/Lyntide

ATP : 100 μ M