

Product Information

EPHB2

Product Number : **08-129**

Product description

Human EPHB2, cytoplasmic domain [581-987(end) amino acids of accession number NP_004433.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHB2 was purified by using glutathione sepharose chromatography.

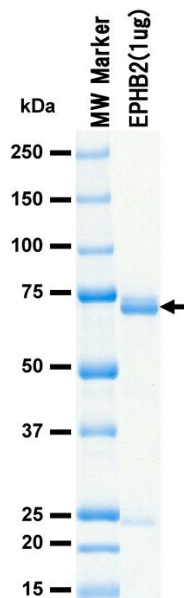
Storage buffer:

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

Storage and Handling:

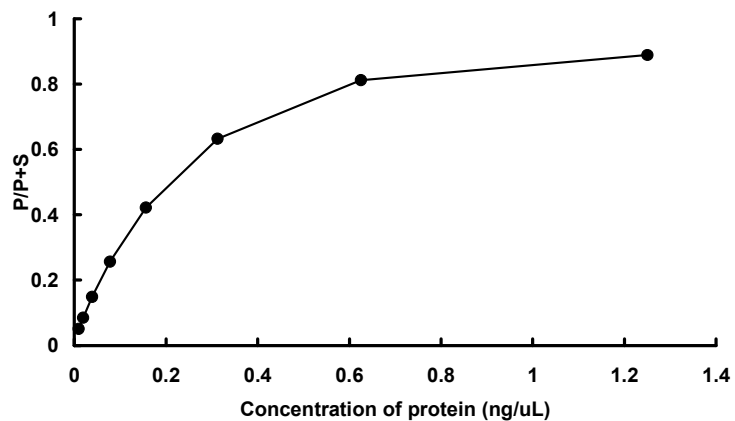
Store at -80C.
Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 83 %
The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : Blk/Lyntide

ATP : 100 μ M