

## Product Information

### EPHA4

Product Number : **08-123**

#### Product description

Human EPHA4, cytoplasmic domain [586-986(end) amino acids of accession number NP\_004429.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-EPHA4 was purified by using glutathione sepharose chromatography.

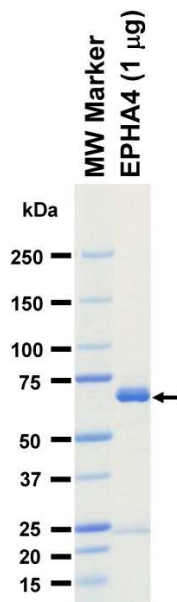
#### Storage buffer:

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.

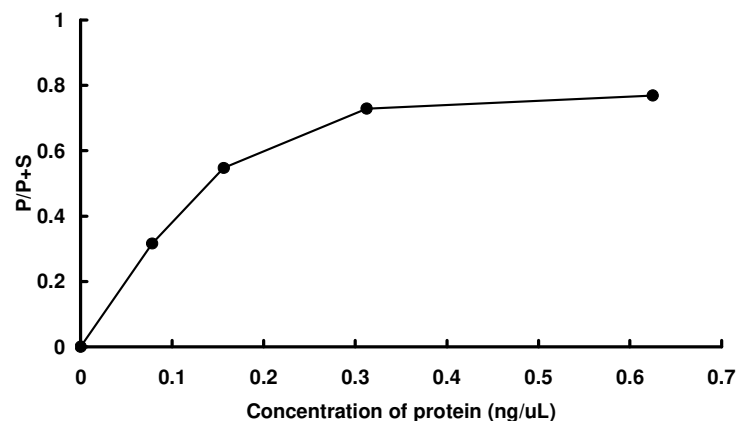
#### SDS-PAGE



Purity: 87 %

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  $\mu$ M