

Product Information

EPHA1

Product Number : **08-119**

Product description

Human EPHA1, cytoplasmic domain [586-976(end) amino acids and M900V of accession number NP_005223.4] was expressed as N-terminal GST-fusion protein (71 kDa) using baculovirus expression system. GST-EPHA1 was purified by using glutathione sepharose chromatography.

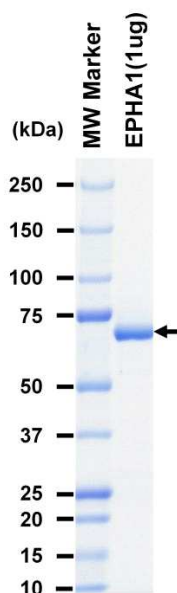
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

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

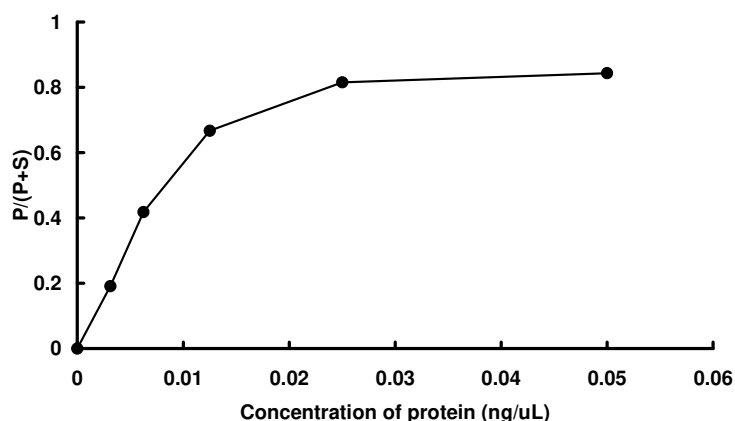
SDS-PAGE



Purity: 97 %

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

Product Information

EPHA1

Product Number : **08-119**

Product description

Human EPHA1, cytoplasmic domain [586-976(end) amino acids and M900V of accession number NP_005223.4] was expressed as N-terminal GST-fusion protein (71 kDa) using baculovirus expression system. GST-EPHA1 was purified by using glutathione sepharose chromatography.

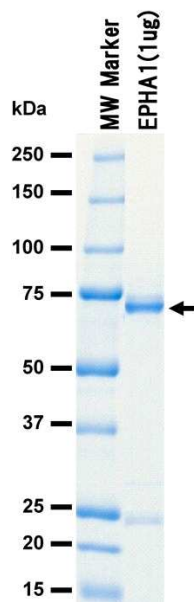
Storage buffer:

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.

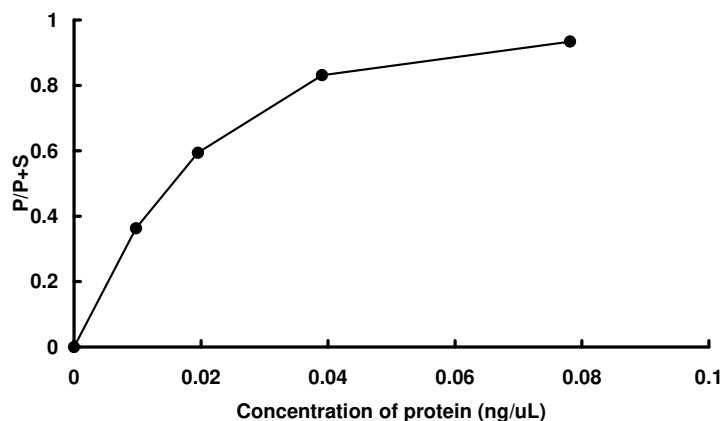
SDS-PAGE



Purity: 82 %

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 µM