

Product Information

BMPR1B

Product Number : **09-138**

Product description

Cytoplasmic domain of BMPR1B [149-502(end) amino acids of accession number NP_001194.1] as N-terminal GST-fusion protein(68 kDa) was co-expressed with BMPR2 [174-1038(end) amino acids of accession number NP_001195.2] using baculovirus expression system. GST-BMPR1B was purified by using glutathione sepharose chromatography.

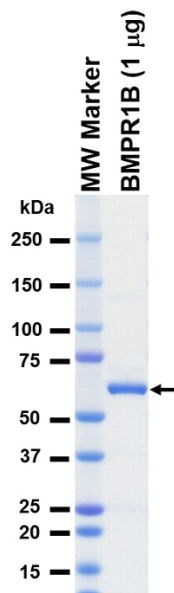
Storage buffer:

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.

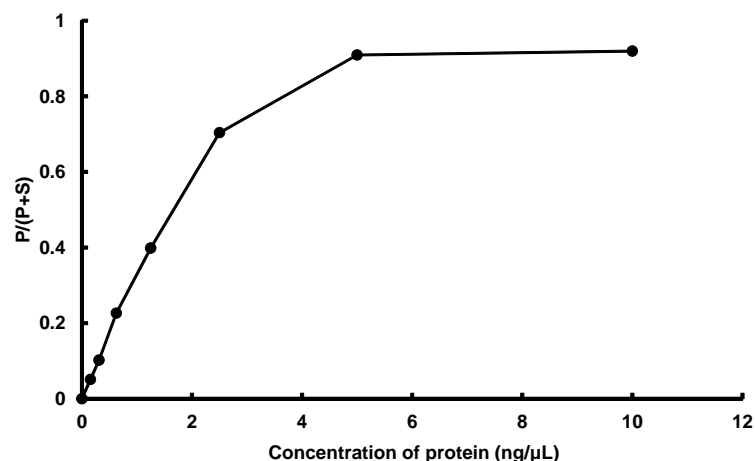
SDS-PAGE



Purity: 95 %

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 Mn/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Topo II a peptide

ATP : 200 µM