

Product Information

ALK4(ACVR1B)

Product Number : **09-135**

Product description

Human ALK4, cytoplasmic domain [150-505(end) amino acids of accession number NP_004293.1] was expressed as N-terminal GST-fusion protein(68 kDa) using baculovirus expression system. GST-ALK4 was purified by using glutathione sepharose chromatography.

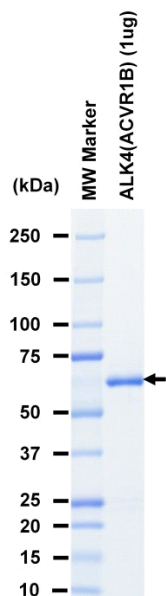
Storage buffer:

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.

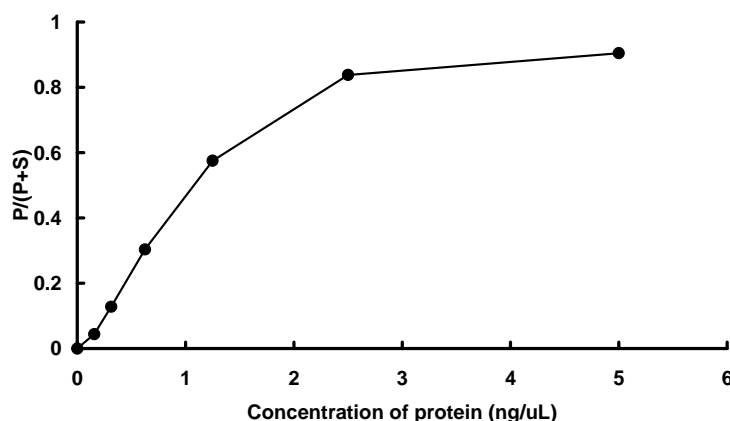
SDS-PAGE



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

Substrate : Topo II a peptide

ATP : 200 μ M