

Product Information

ALK2(ACVR1)[R206H]

Product Number : **09-148**

Product description

Truncated human ALK2 [145-509(end) amino acids and R206H of accession number NP_001096.1] was expressed as N-terminal GST-fusion protein(69 kDa) using baculovirus expression system. GST-ALK2[R206H] was purified by using glutathione sepharose chromatography.

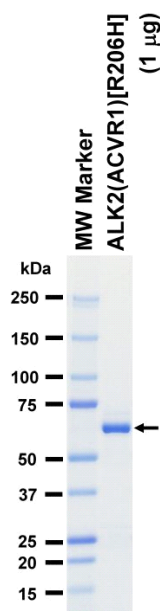
Storage buffer:

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.

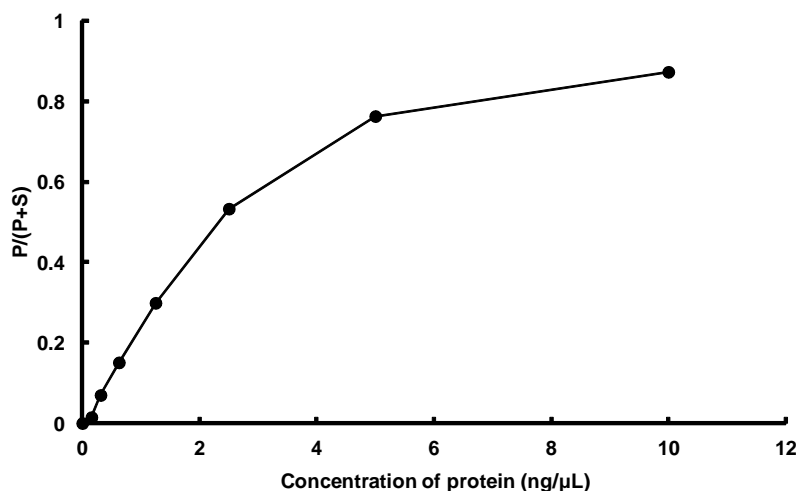
SDS-PAGE



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 µM