

Product Information

ACVRL1(ALK1)

Product Number : **09-133**

Product description

Truncated human ACVRL1 [139-503(end) amino acids of accession number NP_000011.2] was expressed as N-terminal GST-fusion protein(68 kDa) using baculovirus expression system. GST-ACVRL1 was purified by using glutathione sepharose chromatography.

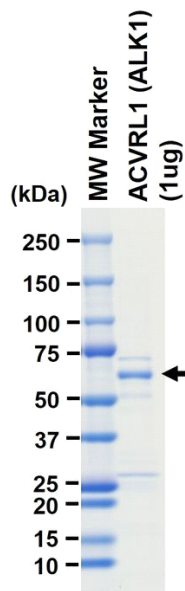
Storage buffer:

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.

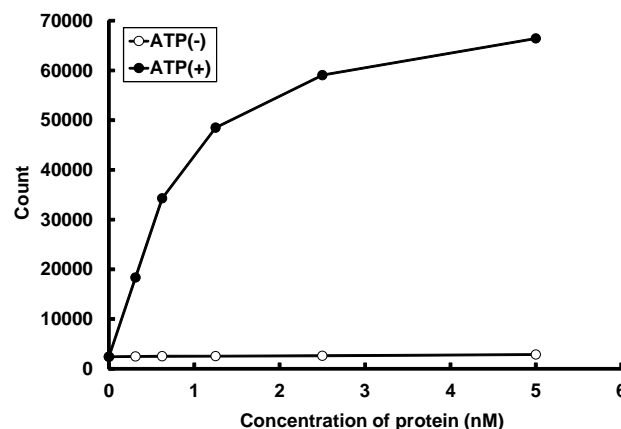
SDS-PAGE



Purity: 65 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was determined by TR-FRET assay. The enzyme was incubated with ULight-labeled peptide and Mn/ATP. Phosphorylation was detected by Eu-labeled anti-phospho antibody.

Substrate : Topo II a peptide

ATP : 100 μ M