

Product Information

ACVR2A

 Product Number : **09-131**

Product description

Human ACVR2A, cytoplasmic domain [162-513(end) amino acids of accession number NP_001607.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-ACVR2A was purified by using glutathione sepharose chromatography.

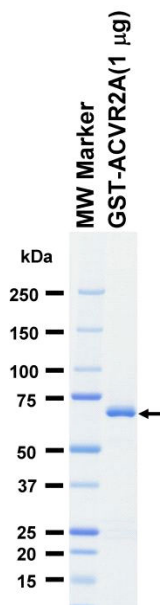
Storage buffer:

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.

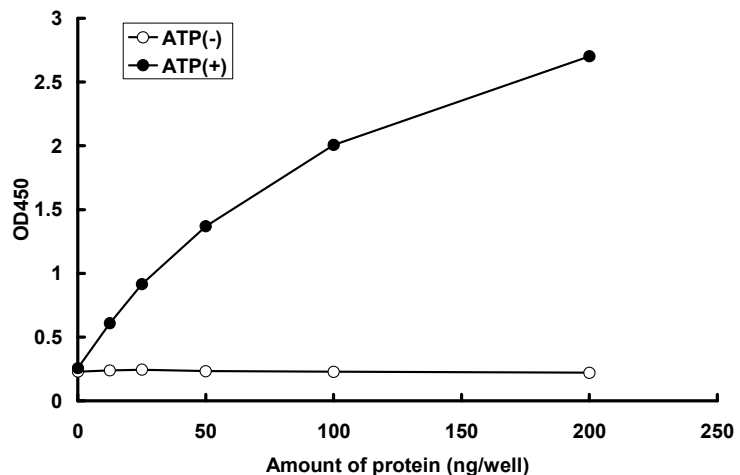
SDS-PAGE



Purity: 95 %

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

Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated substrate protein, and after stopping kinase reaction by EDTA, the reaction solution was transferred into streptavidin-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG.

Substrate : ALK4 inactive mutant

 ATP : 100 μ M