

Product Information

ACTR2B(ACVR2B)

 Product Number : **09-132**

Product description

Human ACTR2B, cytoplasmic domain [161-512(end) amino acids of accession number NP_001097.2] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-ACTR2B was purified by using glutathione sepharose chromatography.

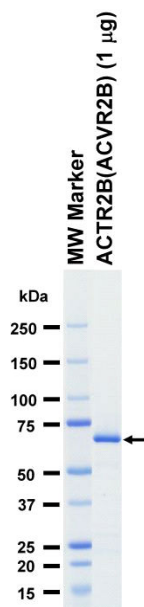
Storage buffer:

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.

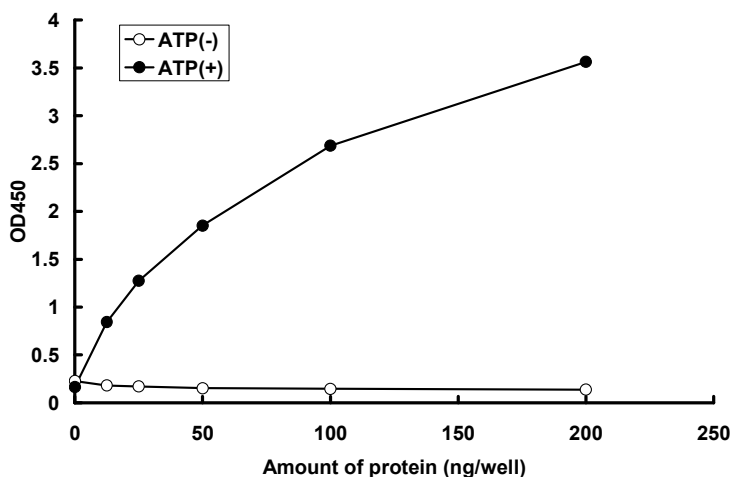
SDS-PAGE



Purity: 94 %

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