

## Product Information

### BTN-ACTR2B(ACVR2B)

 Product Number : **09-432-20N**

#### Product description

Human ACTR2B, cytoplasmic domain [161-512(end) amino acids of accession number NP\_001097.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (43 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

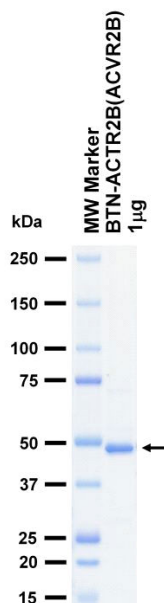
#### Storage buffer:

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.

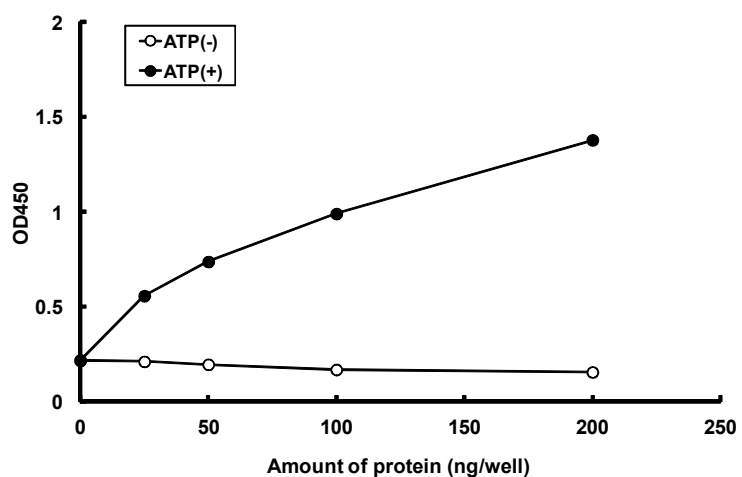
#### SDS-PAGE



Purity: 96 %

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/peptide, 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(or HRP-labeled

Substrate : ALK4 [inactive mutant]

ATP : 100  $\mu$ M