

## Product Information

### BTN-ACVR2A

 Product Number : **09-431-20N**

#### Product description

Human ACVR2A, cytoplasmic domain [162-513(end) amino acids of accession number NP\_001607.1] 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.

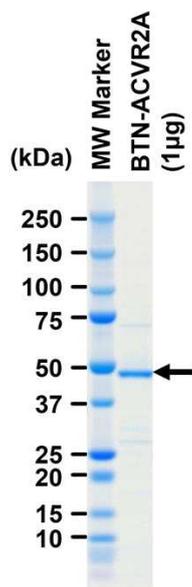
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
 1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

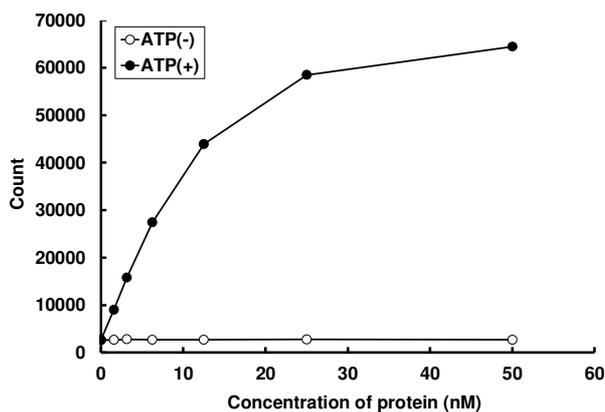
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 84 %  
 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  $\mu$ M