

## Product Information

### BMPR1A

Product Number : **09-137**

#### Product description

Catalytic domain of constitutive active human BMPR1A [Q233D] derived from wild type BMPR1A [187-532(end) amino acids of accession number NP\_004320.2] with N-terminal GST tag was co-expressed with BMPR2 [174-1038(end) amino acids of accession number NP\_001195.2] using baculovirus expression system. GST-BMPR1A (66 kDa) was purified by using glutathione sepharose chromatography.

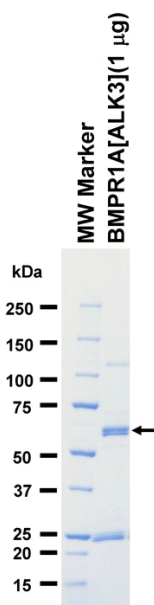
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

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

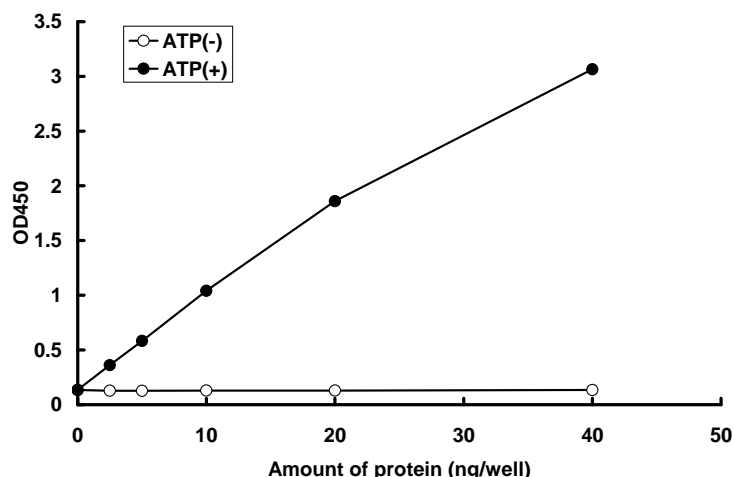
#### SDS-PAGE



Purity: 52 %

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

#### Activity data



The activity was determined by ELISA. The enzyme was incubated with GST-fused substrate protein, and after stopping kinase reaction by EDTA, the reaction solution was transferred into glutathione-coated plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG(or HRP-labeled anti-mouse IgG) .

Substrate : Smad1

ATP : 20 µM