

## Product Information

### BMPR2

Product Number : **09-139**

#### Product description

Human BMPR2, cytoplasmic domain [174-1038(end) amino acids of accession number NP\_001195.2] was expressed as N-terminal GST fusion protein (123 kDa) using baculovirus expression system. GST-BMPR2 was purified by using glutathione sepharose chromatography.

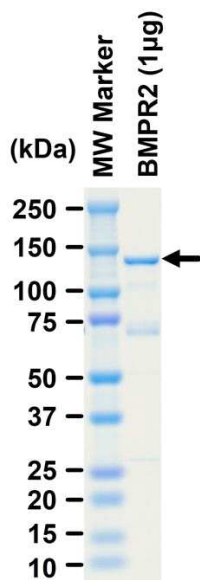
#### Storage buffer:

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.

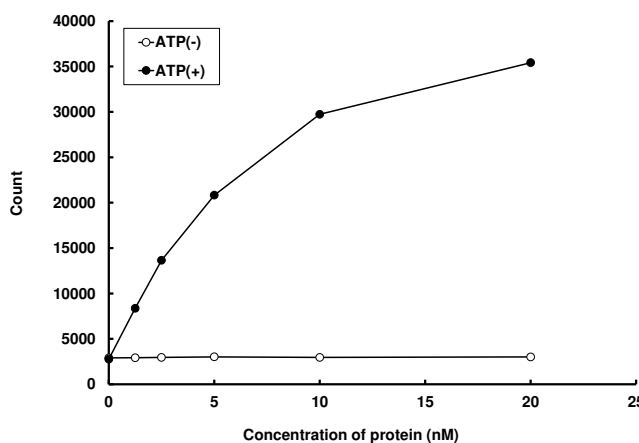
#### SDS-PAGE



Purity: 67 %

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 µM