

## Product Information

### TGFβR2(TGFBR2)

Product Number : **09-142**

#### Product description

Human TGFβR2, catalytic domain [194-567(end) amino acids of accession number NP\_003233.2] was expressed as N-terminal GST-fusion protein (70 kDa) using baculovirus expression system. GST-TGFβR2 was purified by using glutathione sepharose chromatography.

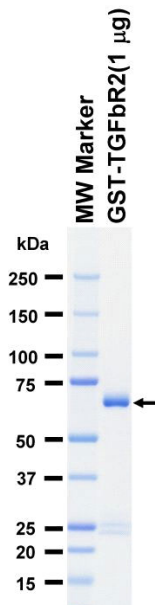
#### Storage buffer:

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.

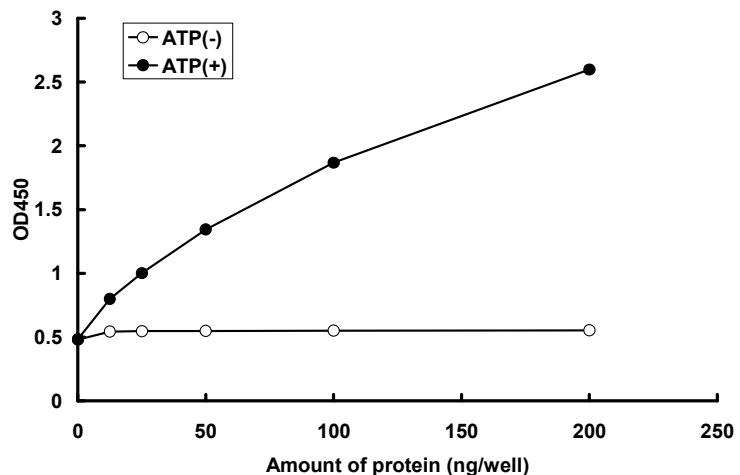
#### SDS-PAGE



Purity: 83 %

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 : Histone H1

ATP : 100 μM