

Product Information

TGF β R1(TGFBR1)

Product Number : 09-141

Product description

Human TGF β R1, catalytic domain [200-503 amino acids of accession number NP_004603.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-TGF β R1 was purified by using glutathione sepharose chromatography.

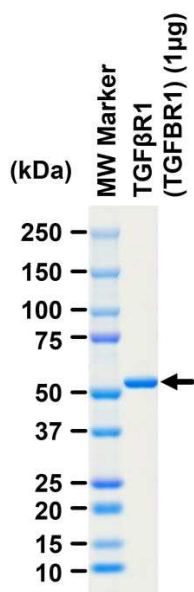
Storage buffer:

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.

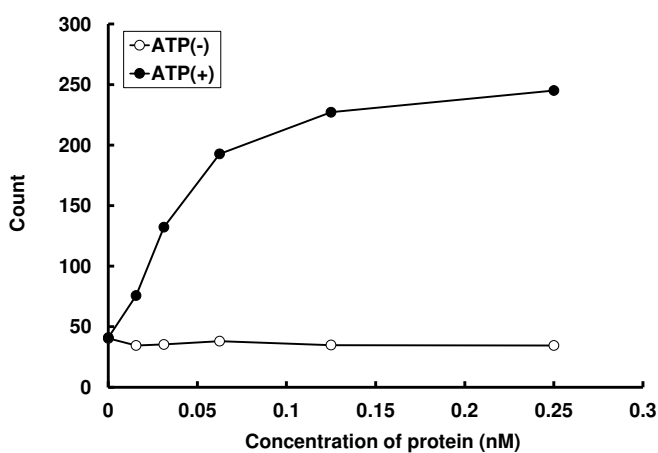
SDS-PAGE



Purity: 99 %

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 μ M