

Product Information

BTN-TGF β R1(TGFBR1)

 Product Number : **09-441-20N**

Product description

Human TGF β R1, catalytic domain [200-503(end) amino acids of accession number NP_004603.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (38 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

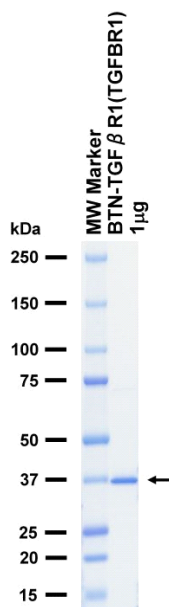
Storage buffer:

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.

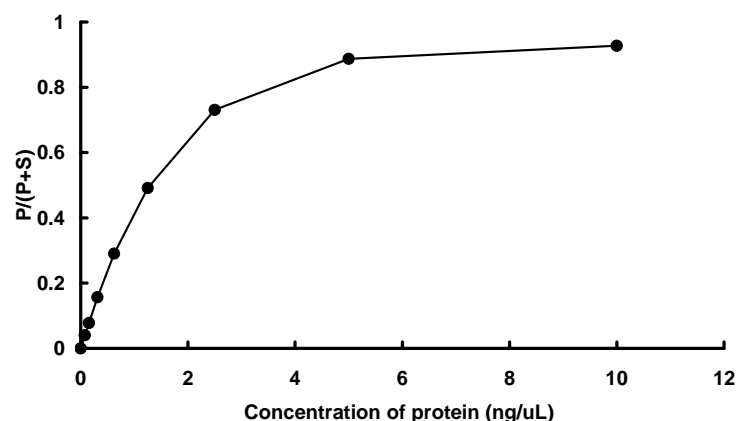
SDS-PAGE



Purity: 98 %

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

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mn/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Topo II a peptide

 ATP : 200 μ M