

Product Information

Smad3

 Product Number : **00-004**

Product description

Full-length human Smad3 [1-425(end) amino acids of accession number NP_005893.1] was expressed as N-terminal GST-fusion protein (75 kDa) using E.coli expression system. GST-Smad3 was purified by using glutathione sepharose chromatography.

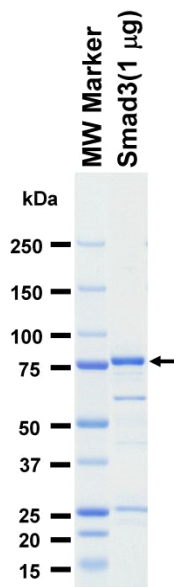
Storage buffer:

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.

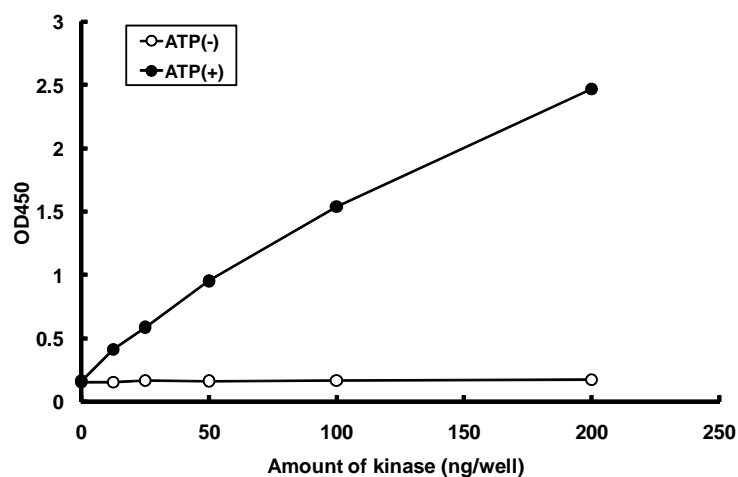
SDS-PAGE



Purity: 61 %

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

Activity data



The phosphorylational ability of the protein was determined by ELISA. The substrate protein was incubated with the indicated enzyme, 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 : TGFbR1

 ATP : 100 μ M