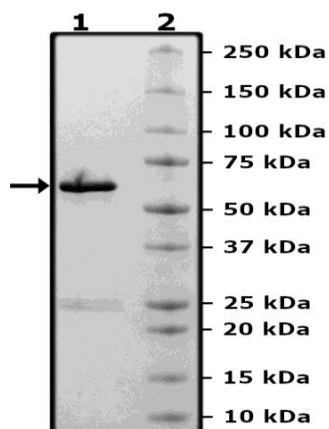


Product Information

Description:	Recombinant human TGFβ1 (transforming growth factor, beta receptor 1), also known as ALK5, encompassing amino acids 80-end. This construct contains an N-terminal GST tag. The recombinant protein was affinity purified and is kinase active.
Species	Human
Construct:	TGFbR1 (GST-80-426(end))
Concentration:	0.19 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 2 mM glutathione, 3 mM DTT, and 25% glycerol
MW:	66 kDa
Genbank Accession:	NM_001130916.2
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	1.8 pmol/min/μg
Assay Conditions:	Assay was done according to TGFβ1 (ALK5) Kinase Assay Kit (BPS Bioscience #78819) with various amounts of TGFBR1 substrate.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



TGFBR1 (ALK5) Activity

