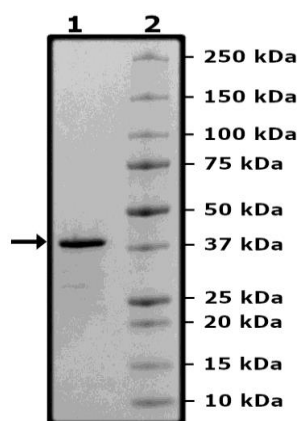


Product Information

Description:	Human Recombinant TRAF6 (TNF receptor associated factor 6), encompassing amino acids 51-159. This construct contains an N-terminal GST-tag followed by a PreScission Site to facilitate removal of the GST-tag.
Species:	Human
Construct:	TRAF6 (GST-Pre-51-159)
Concentration:	1.73 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 2 mM Glutathione
MW:	39 kDa
Genbank Accession:	NM_145803
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.
Assay Conditions:	TRAF6 activity was measured using varying amounts of TRAF6 and incubated with E1, E2, and ATP in the presence of an “intrachain Ubiquitin” mixture for 60 minutes at 30°C in a 20-μl reaction. After incubation, TR-FRET was detected (340=>620 nm, and 340=>665 nm). The TR-FRET is proportional to the auto polyubiquitination of TRAF6.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



TRAF6 Activity

