

Product Information

Description:	Recombinant human 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 the removal of the GST-tag. The construct also contains mutation of interest (D57K). The recombinant protein was affinity purified.
Species	Human
Construct:	TRAF6 (D57K) (GST-Pre-51-159)
Mutation:	D57K
Concentration:	0.66 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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

