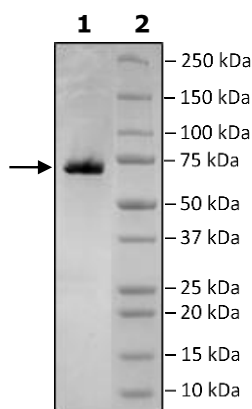


Product Information

Construct:	TDP2 (His-GST-2-362(end))
Concentration:	0.22 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 250 mM imidazole
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_016614
MW:	69 kDa
Purity:	≥90%
Specific Activity:	≥186 pmol/min/μg
Assay Conditions:	Enzyme reaction was performed in a reaction buffer (200 μl) containing 50 mM Tris (pH 7.4), 1 mM MgCl ₂ , 50 mM KCl, 1 mM DTT, 0.1 mg/ml BSA, 2 mM T5PNP (Sigma #T4510) and 600 ng TDP2. Reaction was done at 415 nm at room temperature for 60 min. Specific activity was calculated based on the molar extinction coefficient of the product (p-nitrophenol), $e = 15,000 \text{ M}^{-1}\text{cm}^{-1}$
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



TDP2 Activity

