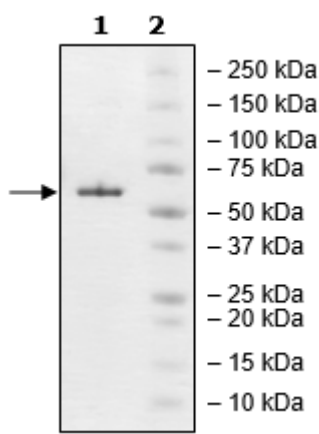


Product Information

Construct:	Tankyrase 2 (GST-849-1166(end))
Concentration:	0.24 mg/ml
Species:	Human
Formulated In:	40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT and 20% glycerol
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_025235
MW:	62 kDa
Purity:	≥90%
Assay Conditions:	Enzyme reaction was conducted following the protocol for the TNKS2 Histone Ribosylation Assay Kit (Biotin-labeled NAD+) (BPS Bioscience #80578).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



TNKS2 Activity

