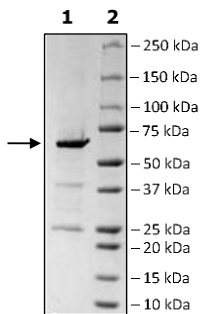


## Product Information

<b>Construct:</b>	TNKS1 (GST-1001-1327(end))
<b>Concentration:</b>	0.13 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, and 20% glycerol.
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_003747
<b>MW:</b>	63 kDa
<b>Purity:</b>	80%
<b>Assay Conditions:</b>	Enzyme reaction is conducted using the TNKS1 Histone Ribosylation Assay Kit (Biotin-labeled NAD <sup>+</sup> ) (BPS Bioscience #80573) with TNKS1 titrated from 100 ng/rxn – 1.6 ng/rxn.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



### TNKS1 Activity

