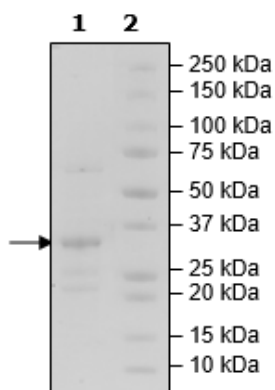


Product Information

Construct:	NMNAT2 (His-2-307(end))
Concentration:	0.19 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 200 mM Imidazole.
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_015039
MW:	35 kDa
Purity:	66%
Specific Activity:	390 pmol/min/μg
Assay Conditions:	50 mM Tris/HCl, pH 7.4, 5 mM MgCl ₂ , 0.05% Tween-20, 0.1 mM NMN, 0.1mM ATP, and NMNAT2 (3.2-200 ng) in a final volume of 50 μL. The reaction was incubated for 30 min at 37°C. The amount of ATP consumed was calculated using Kinase-Glo plus reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



NMNAT2 Activity

