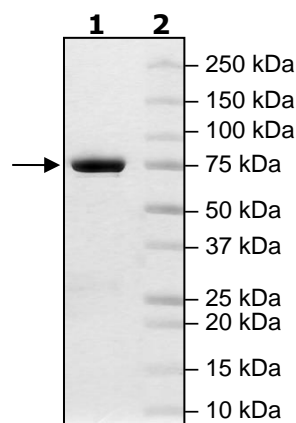


Product Information

| | |
|---------------------------|---|
| Construct: | NAMPT (GST-2-491) |
| Concentration: | 1.70 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris pH8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween, 10 mM GSH, 20% glycerol, 3 mM DTT. |
| 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_005746 |
| MW: | 82 kDa |
| Purity: | ≥90% |
| Specific Activity: | ≥1.0 pmol/min/μg |
| Assay Conditions: | The reaction mixture contained 50 mM Tris-HCl, pH 8.0, 12.5 mM MgCl ₂ , 0.1 mM nicotinamide, 0.4 mM PRPP, 2 mM ATP, 30 μg/mL of alcohol dehydrogenase, 10 μg/mL of NMNAT (lot 130419), 1.5% alcohol, 1 mM DTT, 0.02% BSA, and various amount of NAMPT in a final volume of 50 μL. Reactions were incubated at room temperature for 30 min, and the increase in NADH fluorescence (Ex 340 nm/Em 460 nm) was measured. |
| Applications: | Useful for the study of protein binding assays, enzyme kinetics, screening inhibitors, and selectivity profiling. |

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

