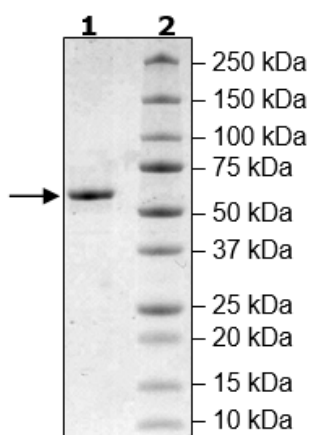


Product Information

Construct:	FYN (FLAG-2-537(end))
Concentration:	0.32 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 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_002037
MW:	61 kDa
Purity:	≥90%
Specific Activity:	240 pmol/min/g
Assay Conditions:	Activity was measured using Kinase-Glo Plus™ with 0.2 mg/ml poly(Glu-Tyr) and 20 mM ATP and increasing amounts of FYN. Assay was incubated 30°C for 50 minutes.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



FYN Activity

