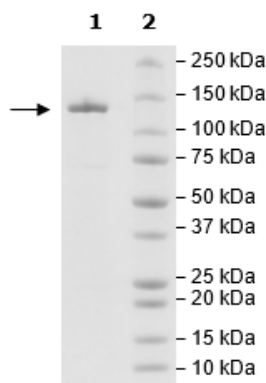


Product Information

Construct:	USP7 (His-FLAG-2-1102(end))
Concentration:	7.80 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_003470
MW:	130 kDa
Purity:	≥90%
Specific Activity:	545 pmol/min/μg
Assay Conditions:	This Assay was performed according to USP7 Inhibitor Screening Assay Kit (BPS Bioscience #79256) with USP7 titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



USP7 Activity

