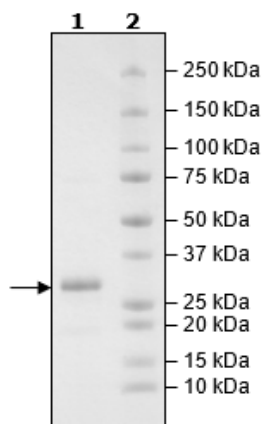


Product Information

Construct:	3CL Protease (1-306(end)-His) (SARS-CoV-1)
Concentration:	0.42 mg/ml
Species:	SARS-CoV-1
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 200 mM imidazole, and 20% glycerol
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:	1UK3_B
MW:	36 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to 3CL Protease, MBP-tagged (SARS-CoV-2) Assay Kit (BPS Bioscience #79955-1) with 3CL Protease (SARS-CoV-1) titrated at various concentrations and incubated for 1 hour at room temperature.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



3CL Protease (SARS-CoV-1) Activity

