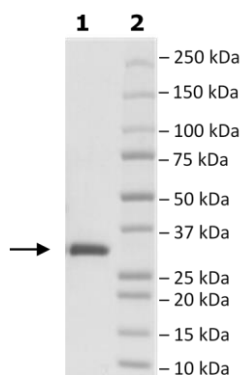


Product Information

Construct:	3CL-Protease (SARS-CoV-2) (1-306(end))
Concentration:	9.0 mg/ml
Species:	SARS-CoV-2
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 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:	YP_009725301
MW:	34 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to the 3CL Protease, MBP-tagged (SARS-CoV-2) Assay Kit (BPS Bioscience #79955-1) with 3CL Protease titrated from 60 ng/rxn to 0.9 ng/rxn.
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-2) Activity

