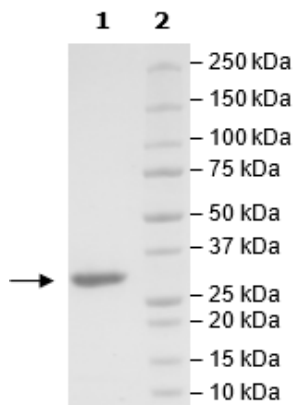


Product Information

Construct:	3CL Protease (B.1.1.529, Omicron Variant) (1-306(end)) (SARS-CoV-2)
Mutation:	P132H
Concentration:	5.99 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 (B.1.1.529, Omicron Variant) titrated from 90 ng/rxn to 0 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 (Omicron Variant) Activity

