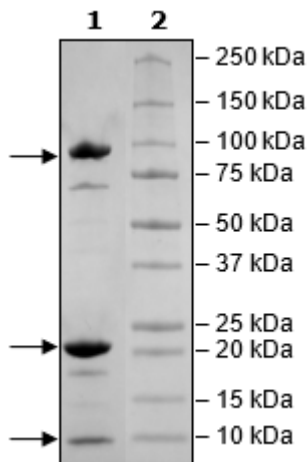


Product Information

Construct:	RdRp (FLAG-1-932(end)) / NSP7 (1-84(end)-His) / NSP8 (1-198(end)-His)
Concentration:	1.14 mg/ml
Species:	SARS-CoV-2
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 4 mM MgCl ₂ , 1 mM TCEP, and 10% glycerol
Expression System:	RdRp: HEK293; NSP7: <i>E. coli</i> ; NSP8: <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:	RdRp: YP_009725307; NSP7: YP_009725303.1; NSP8: YP_009725304.1
MW:	RdRp: 108 kDa; NSP7: 10 kDa; NSP8: 23 kDa
Purity:	≥90%
Assay Conditions:	RNA-dependent RNA polymerase was incubated with FRET Dig-RNA Duplex in the presence of RNaseOUT and biotinylated ATP for 80 min at 37°C in a 10- μ l reaction. Eu-labeled α -Dig antibody and Streptavidin-acceptor dye was then added followed by TR-FRET detection (317 => 620 nm, and 317 => 665 nm). Increase in TR-FRET is proportional to the amount of ATP incorporated in the RNA.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



RdRp Activity

