

# NSP8\_NSP7/NSP12 (SARS-CoV-2) Complex Recombinant

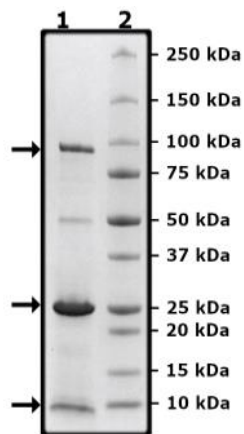
Catalog: 101466  
Lot: 230227-G2

## Product Information

<b>Description:</b>	Recombinant complex of three SARS-CoV-2 proteins NSP7 (nonstructural protein 7), NSP8 (nonstructural protein 8) and NSP12 (nonstructural protein 12). The NSP8_NSP7 construct contains an N-terminal His-tag (14xHis) followed by a TEV protease target sequence that allows cleavage of His-tag (TEV: tobacco etch virus cysteine protease BPS Bioscience #50308). The recombinant NSP12 construct is full length encompassing amino acids 1-932 (end). These recombinant proteins are co-expressed in an E. coli expression system and affinity purified.
<b>Species</b>	SARS-CoV-2
<b>Construct:</b>	NSP8_NSP7 (His-TEV-1-242(end)) / NSP12 (1-932(end)) (SARS-CoV-2)
<b>Concentration:</b>	0.53 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	20 mM HEPES, pH 8.0, 300 mM NaCl, 10% glycerol, and 3 mM DTT
<b>MW:</b>	NSP12: 108 kDa; NSP8: 23 kDa; NSP7: 10 kDa
<b>Genbank Accession:</b>	NSP7: YP_009742614; NSP8: YP_009725304.1; NSP12: YP_009725307
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	Activity was measured using the RdRp (SARS-CoV-2) TR-FRET Assay Kit (BPS Bioscience #78394) with Nsp8_Nsp7/Nsp 12 (Sars-CoV-2) complex titrated at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



RdRp Activity

