

Product Information

Construct:	NSP10 (FLAG-1-139(end)) / NSP16 (His-1-298(end)) (SARS-CoV-2)
Concentration:	0.13 mg/ml
Species:	SARS-CoV-2
Formulated In:	20 mM HEPES, pH 7.4, 150 mM NaCl, 4 mM MgCl ₂ , and 10% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NSP10: YP_009742617; NSP16: YP_009725311
MW:	NSP10: 16 kDa; NSP16: 34 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

