

Product Information

Construct:	NSP8 (1-198(end)-His) (SARS-CoV-2)
Concentration:	1.21 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, 10% glycerol
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 009725304.1
MW:	23 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

