

Product Information

Construct:	ORF9b (GST-Pre-1-97(end))(SARS-CoV-2)
Concentration:	1.01 mg/ml
Species:	SARS-CoV-2
Formulated In:	40 mM Tris-HCl, pH 8.3, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 20 mM Glutathione and 20% 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
Uniprot:	P0DTD2
MW:	38 kDa
Purity:	>90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

