

Product Information

Construct:	Nucleocapsid (P.1) (SARS-CoV-2) (1-419(end)-Avi-His)
Mutation:	P80R
Concentration:	0.32 mg/ml
Species:	SARS-CoV-2
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% 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:	QHO62115.1
MW:	48 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

