

Product Information

Construct:	Nucleocapsid (B.1.1.529 Variant) (1-419(end)-Avi-His) (SARS-CoV-2)
Substitutions:	P13L, R203K, G204R
Deletions:	Δ31-33ERS
Concentration:	0.58 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, 250 mM imidazole
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	YP_009724397.2
MW:	48 kDa
Purity:	74%

Quality Control Data

4-20% SDS-Page Coomassie Staining

