

## Product Information

<b>Construct:</b>	Nucleocapsid (B.1.1.529 Variant) (1-419(end)-Avi-His) (SARS-CoV-2)
<b>Substitutions:</b>	P13L, R203K, G204R
<b>Deletions:</b>	Δ31-33ERS
<b>Concentration:</b>	0.34 mg/ml
<b>Species:</b>	SARS-CoV-2
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 250 mM imidazole
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	YP_009724397.2
<b>MW:</b>	48 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining

