

## Product Information

<b>Construct:</b>	Nucleocapsid (B.1.1.7) (SARS-CoV-2) (1-419(end)-Avi-His)
<b>Mutation:</b>	D3L, R203K, G204R, S235F
<b>Concentration:</b>	0.44 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
<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>	QHO62115.1
<b>MW:</b>	48 kDa
<b>Purity:</b>	>90%

## Quality Control Data

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

