

## Product Information

<b>Construct:</b>	RNA Polymerase, (SARS-CoV-2) (FLAG-1-932(end))
<b>Concentration:</b>	0.36 mg/ml
<b>Species:</b>	SARS-CoV-2
<b>Formulated In:</b>	45 mM Tris-HCl pH 8.0, 124 mM NaCl, 2.4 mM KCl, 4 mM MgCl <sub>2</sub> , 10% 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>	YP_009725307
<b>MW:</b>	108 kDa
<b>Purity:</b>	85%

## Quality Control Data

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

