

Product Information

Construct:	PLP2(His-1565-1894)(HCoV-NL63)
Concentration:	5 μ M
Species:	NL63 Coronavirus
Formulated In:	50 mM Tris-HCl pH 8.5, 150 mM NaCl, 1 mM DTT, 10% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	P0C6U6.1
MW:	37 kDa