

Product Information

Construct:	VP39 (His-2-333(end)) (Vaccinia Virus)
Concentration:	0.95 mg/ml
Species:	Vaccinia Virus
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% 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:	P07617
MW:	40 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

