



CERTIFICATE OF ANALYSIS

Product Description: Orthopoxvirus Serology Calibrator 1
Catalog Number: C0686-2
Lot Number: A0080406
Expiration Date: 31 OCT 2034
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
VACV A27L	258	AU/mL
VACV A33R	896	AU/mL
VACV B5R	209	AU/mL
VACV D8L	180	AU/mL
VACV L1R	111	AU/mL
MPXV M1R	94.7	AU/mL
MPXV E8L	117	AU/mL
MPXV B6R	156	AU/mL
MPXV A35R	604	AU/mL
MPXV A29L	195	AU/mL

Dilution to Top of Curve: 10X

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		10 DEC 2024