

**CERTIFICATE OF ANALYSIS**

**Control Pack Description:** Orthopoxvirus Serology Control Pack 1  
**Catalog Number:** C4686-1  
**Lot Number:** CP000141  
**Expiration Date:** 31 JAN 2028

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Orthopoxvirus Serology Control 1	A00C0880	≤ -70°C	31 JAN 2028
Orthopoxvirus Serology Control 2	A00C0881	≤ -70°C	31 JAN 2028
Orthopoxvirus Serology Control 3	A00C0882	≤ -70°C	31 JAN 2028

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
VACV A27L	6.81	1.41	<0.100	AU/mL
VACV A33R	36.3	8.63	<0.100	AU/mL
VACV B5R	11.3	6.25	<0.100	AU/mL
VACV D8L	9.96	2.07	<0.100	AU/mL
VACV L1R	6.68	1.55	<0.200	AU/mL
MPXV M1R	5.88	1.33	<0.200	AU/mL
MPXV E8L	6.95	1.24	<0.100	AU/mL
MPXV B6R	9.04	6.03	<0.100	AU/mL
MPXV A35R	19.8	7.83	<0.100	AU/mL
MPXV A29L	5.85	1.28	<0.100	AU/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		30 OCT 2024