

**CERTIFICATE OF ANALYSIS****Control Pack Description:** Orthopoxvirus Serology Control Pack 1**Catalog Number:** C4686-1**Lot Number:** CP000188**Expiration Date:** 31 DEC 2035**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Orthopoxvirus Serology Control 1	A00C1019	≤ -70°C	31 DEC 2035
Orthopoxvirus Serology Control 2	A00C1020	≤ -70°C	31 DEC 2035
Orthopoxvirus Serology Control 3	A00C1046	≤ -70°C	31 DEC 2035

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3		Unit of Measure
			Sector	QuickPlex Ultra	
VACV A27L	6.93	1.58	<0.1	<0.3	AU/mL
VACV A33R	35.8	9.41	<0.1	<0.2	AU/mL
VACV B5R	11.9	7.37	<0.1	<0.4	AU/mL
VACV D8L	10.7	2.09	<0.1	<0.2	AU/mL
VACV L1R	6.94	1.46	<0.2	<0.4	AU/mL
MPXV M1R	5.75	1.17	<0.2	<0.4	AU/mL
MPXV E8L	6.71	1.30	<0.1	<0.2	AU/mL
MPXV B6R	9.65	7.07	<0.1	<0.4	AU/mL
MPXV A35R	19.8	8.12	<0.1	<0.2	AU/mL
MPXV A29L	5.22	1.16	<0.1	<0.3	AU/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 MAR 2026