



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® Orthopoxvirus Panel 1 Kit
Kit Catalog Numbers: K15688-Series
Kit Lot Number: K0082230
Expiration Date: 31 JUL 2025

Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
Orthopoxvirus Plate 1	Z0057335	2-8°C	31 JUL 2025
Orthopoxvirus Serology Calibrator 1	A0080357	≤ -70°C	31 JAN 2028
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
SULFO-TAG™ Anti-Human IgG Antibody	D00V0029	2-8°C	31 JUL 2026
Diluent 100	Not Kit Specific	2-8°C	See Label
MSD Blocker A Kit	Not Kit Specific	Room Temperature	See Label
MSD GOLD Read Buffer B	Not Kit Specific	Room Temperature	See Label

Coating Confirmation Testing Results:

Spot	Description	Result
1	VACV A27L	Pass
2	VACV A33R	Pass
3	VACV B5R	Pass
4	VACV D8L	Pass
5	VACV L1R	Pass
6	MPXV M1R	Pass
7	MPXV E8L	Pass
8	MPXV B6R	Pass
9	MPXV A35R	Pass
10	MPXV A29L	Pass



CERTIFICATE OF ANALYSIS

Page 2 of 2

Functional Testing Results:

Sample Type	Calibrator	Controls			Samples	
Metric	Avg Signal Ratio Test:Reference	Positive Control Recovery	Positive Control Conc. CV	Negative Control	Avg % Difference Test:Reference	Slope Test:Reference
VACV A27L	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
VACV A33R	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
VACV B5R	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
VACV D8L	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
VACV L1R	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
MPXV M1R	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
MPXV E8L	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
MPXV B6R	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
MPXV A35R	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3
MPXV A29L	50 – 200%	70 – 130%	≤20%	Pass	± 25%	0.70 – 1.3

Additional Comments:

All kit components were manufactured and tested according to MSD documents. The lots listed in the Kit Components table meet MSD's specifications.

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 APR 2023