



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX Respiratory Panel 7 Kit
Lot Number: K0082844
Expiration Date: 31 MAR 2027
Catalog Numbers: K15735U-Series; K15736U-Series

Kit Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Respiratory Plate 7	Z0057518	2-8°C	31 MAR 2027
MSD GOLD SULFO-TAG™ Anti-Human IgG Antibody	D00V0043	2-8°C	31 JUL 2027
Reference Standard 2	A0080403	≤ -70°C	30 JUN 2034
Serology Control 2.1	A00C0959	≤ -70°C	30 JUN 2034
Serology Control 2.2	A00C0960	≤ -70°C	30 JUN 2034
Serology Control 2.3	A00C0961	≤ -70°C	30 JUN 2034
SULFO-TAG™ Anti-Human IgA Antibody	Not Kit Specific	2-8°C	See Label
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

Plate Uniformity Testing Results:

Parameter	Precision			Uniformity	Signal
Metric	CV of Intra-plate Averages	Average Intra-plate CV	Max Intra-plate CV	Average Uniformity Metric	Average Signal
SARS-CoV-2 Spike	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu A/Wisconsin/67/2022 H1	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 N	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (XBB.1.5 (v2))	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu A/Ghana/39/2021 H5	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RSV Pre-Fusion F	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu A/Massachusetts/18/2022 H3	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu B/Austria/2021 HA	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (KP.2)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000



CERTIFICATE OF ANALYSIS

Page 2 of 2

Coating Confirmation Testing Results:

Spot	Description	Result
1	SARS-CoV-2 Spike	Pass
2	Flu A/Wisconsin/67/2022 H1	Pass
3	SARS-CoV-2 N	Pass
5	SARS-CoV-2 Spike (XBB.1.5 (v2))	Pass
6	Flu A/Ghana/39/2021 H5	Pass
7	RSV Pre-Fusion F	Pass
8	Flu A/Massachusetts/18/2022 H3	Pass
9	Flu B/Austria/2021 HA	Pass
10	SARS-CoV-2 Spike (KP.2)	Pass

Functional Testing Results:

Sample Type	Calibrator	Controls		Samples	
Metric	Avg Signal Ratio Test:Reference	Control Recovery	Control Conc. CV	Avg % Difference Test:Reference	Slope Test:Reference
SARS-CoV-2 Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu A/Wisconsin/67/2022 H1	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
SARS-CoV-2 N	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
SARS-CoV-2 Spike (XBB.1.5 (v2))	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu A/Ghana/39/2021 H5	50 – 200%	70 – 130%	< 20%	N/A	N/A
RSV Pre-Fusion F	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu A/Massachusetts/18/2022 H3	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu B/Austria/2021 HA	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
SARS-CoV-2 Spike (KP.2)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2


Additional Comments:

Functional testing was executed using Respiratory Plate 7 (Z0057518), SULFO-TAG Anti-Human IgG Antibody (D00V0036), Reference Standard 2 (A0080403), and Serology Controls (A00C0959, A00C0960, A00C0961).

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		16 DEC 2025