



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX<sup>®</sup> Respiratory Panel 1 Kit  
**Kit Catalog Numbers:** K15365-Series; K15366-Series; K15367-Series; K15377-Series  
**Kit Lot Number:** K0082037  
**Expiration Date:** 31 JAN 2024

### Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
Respiratory Plate 1	Z0057180	2-8°C	31 OCT 2024
Reference Standard 1	A0080286	≤ -70°C	28 FEB 2026
Serology Control 1.1	A00C0825	≤ -70°C	31 MAR 2027
Serology Control 1.2	A00C0826	≤ -70°C	31 MAR 2027
Serology Control 1.3	A00C0827	≤ -70°C	31 MAR 2027
SULFO-TAG <sup>™</sup> Human ACE2 Protein	D0081760	2-8°C	31 JAN 2024
ACE2 Calibration Reagent	A0080301	2-8°C	31 MAR 2025
SULFO-TAG <sup>™</sup> Anti-Human IgG Antibody	D00V0025	2-8°C	30 SEP 2025
SULFO-TAG <sup>™</sup> Anti-Human IgM Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG <sup>™</sup> 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

See product insert regarding which components are provided with each IgG, IgA, IgM, and ACE2 kit.

### Plate Uniformity Testing Results:

Parameter	Precision			Uniformity	Signal
	CV of Intra-plate Averages	Average Intra-plate CV	Max Intra-plate CV		
Flu B/Brisbane HA	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu A/Shanghai H7	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu A/Michigan H1	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RSV Pre-Fusion F	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu A/HongKong H3	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Flu B/Phuket HA	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000

### Coating Confirmation Testing Results:

Spot	Description	Result
1	Flu B/Brisbane HA	Pass
2	Flu A/Shanghai H7	Pass
4	Flu A/Michigan H1	Pass
7	RSV Pre-Fusion F	Pass
8	Flu A/HongKong H3	Pass
10	Flu B/Phuket HA	Pass



## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Functional Testing Results:

Sample Type	Calibrator	Controls		Samples	
Metric	Average Signal Ratio Test:Reference	Control Recovery	Control Conc. CV	Avg % Difference Test:Reference	Slope Test:Reference
Flu B/Brisbane HA	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu A/Shanghai H7	50 – 200%	70 – 130%	< 20%	± 25%	0.78
Flu A/Michigan H1	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RSV Pre-Fusion F	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu A/HongKong H3	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Flu B/Phuket HA	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2

### Additional Comments:

Functional testing was executed using Respiratory Plate 1 (Z0057180), SULFO-TAG Anti-Human IgG Antibody (D00V0025), Reference Standard 1 (A0080286), and Serology Controls (A00C0825, A00C0826, A00C0827).

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		08 JUN 2022