



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® SARS-CoV-2 Panel 28 Kit  
**Kit Catalog Numbers:** K15614-Series; K15615-Series; K15616-Series; K15617-Series; K15618-Series; K15619-Series; K15620-Series  
**Kit Lot Number:** K0082053  
**Expiration Date:** 31 JAN 2024

### Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
SARS-CoV-2 Plate 28	Z0057188	2-8°C	30 NOV 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™ Human ACE2 Protein	D0081760	2-8°C	31 JAN 2024
ACE2 Calibration Reagent 2	A0080327	2-8°C	31 OCT 2025
SULFO-TAG™ Anti-Human IgG Antibody	D00V0024	2-8°C	31 AUG 2025
SULFO-TAG™ Anti-Human IgM Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG™ Anti-Human IgA Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG™ Anti-Mouse IgG Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG™ Anti-Mouse IgM Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG™ Anti-Mouse 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, or ACE2 kit.

### Plate Uniformity Testing Results:

Parameter	Precision			Uniformity	Signal
	CV of Intra-plate Averages	Average Intra-plate CV	Max Intra-plate CV	Average Uniformity Metric	Average Signal
RBD (BA.2.12.1)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (Beta) (B.1.351)	≤ 25%	≤ 15%	≤ 18%	Pass	1,500 – 1,000,000
RBD (BA.2)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (BA.2+L452M)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (BA.2+L452R)	≤ 18%	≤ 10%	15.4%	Pass	1,500 – 1,000,000
RBD (Alpha) (B.1.1.7)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (BA.4; BA.5)	≤ 18%	≤ 10%	17.0%	Pass	1,500 – 1,000,000
RBD (BA.3)	≤ 18%	11.5 %	17.8%	Flag	1,500 – 1,000,000
RBD (Delta) (L452R,T478K)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 S1 RBD	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000



## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Coating Confirmation Testing Results:

Spot	Description	Result
1	RBD (BA.2.12.1)	Pass
2	RBD (Beta) (B.1.351)	Pass
3	RBD (BA.2)	Pass
4	RBD (BA.2+L452M)	Pass
5	RBD (BA.2+L452R)	Pass
6	RBD (Alpha) (B.1.1.7)	Pass
7	RBD (BA.4; BA.5)	Pass
8	RBD (BA.3)	Pass
9	RBD (Delta) (L452R,T478K)	Pass
10	SARS-CoV-2 S1 RBD	Pass

**RBD** indicates SARS-CoV-2 Receptor Binding Domain with mutation(s) listed.

### 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
RBD (BA.2.12.1)	N/A	N/A	N/A	N/A	N/A
RBD (Beta) (B.1.351)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (BA.2)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (BA.2+L452M)	N/A	N/A	N/A	N/A	N/A
RBD (BA.2+L452R)	N/A	N/A	N/A	N/A	N/A
RBD (Alpha) (B.1.1.7)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (BA.4; BA.5)	N/A	N/A	N/A	N/A	N/A
RBD (BA.3)	N/A	N/A	N/A	N/A	N/A
RBD (Delta) (L452R,T478K)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
SARS-CoV-2 S1 RBD	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2

### Additional Comments:

Functional testing was executed using SARS-CoV-2 Plate 28 (Z0057188), SULFO-TAG Anti-Human IgG Antibody (D00V0024), 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
<b>Review/Approval</b>	Karen Hamilla	Quality		17 JUN 2022