



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX<sup>®</sup> SARS-CoV-2 Panel 30 Kit  
**Kit Catalog Numbers:** K15632-Series; K15633-Series; K15634-Series; K15635-Series; K15636-Series; K15637-Series; K15638-Series  
**Kit Lot Number:** K0082103  
**Expiration Date:** 31 JAN 2024

### Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
SARS-CoV-2 Plate 30	Z0057232	2-8°C	31 JAN 2025
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	A0080327	2-8°C	31 OCT 2025
SULFO-TAG <sup>™</sup> Anti-Human IgG Antibody	D00V0027	2-8°C	30 NOV 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
SULFO-TAG <sup>™</sup> Anti-Mouse IgG Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG <sup>™</sup> Anti-Mouse IgM Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG <sup>™</sup> 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
	Metric	CV of Intra-plate Averages	Average Intra-plate CV		
RBD (BA.2.12.1)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (B.1.351)(Beta)	≤ 25%	≤ 15%	24.2%	Flag	1,500 – 1,000,000
RBD (Omicron; BA.2)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (B.1.1.7)(Alpha)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (BA.4; BA.5)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (BA.2.75)	≤ 25%	≤ 15%	≤ 18%	Flag	1,500 – 1,000,000
RBD (L452R,T478K)(Delta)	≤ 18%	≤ 10%	14%	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 (B.1.351)(Beta)	Pass
3	RBD (Omicron; BA.2)	Pass
6	RBD (B.1.1.7)(Alpha)	Pass
7	RBD (BA.4; BA.5)	Pass
8	RBD (BA.2.75)	Pass
9	RBD (L452R,T478K)(Delta)	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)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (B.1.351)(Beta)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (Omicron; BA.2)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (B.1.1.7)(Alpha)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (BA.4; BA.5)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (BA.2.75)	N/A	N/A	N/A	N/A	N/A
RBD (L452R,T478K)(Delta)	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 30 (Z0057232), SULFO-TAG Anti-Human IgG Antibody (D00V0027), 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		02 SEP 2022