



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX<sup>®</sup> SARS-CoV-2 Panel 2 Kit  
**Kit Catalog Numbers:** K15383-Series; K15384-Series; K15385-Series; K15386-Series; K15475-Series; K15476-Series; K15477-Series  
**Kit Lot Number:** K0082081  
**Expiration Date:** 31 JAN 2024

### Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
SARS-CoV-2 Plate 2	Z0057218	2-8°C	31 DEC 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	A0080309	2-8°C	30 APR 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
	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
SARS CoV-2 Nucleocapsid	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS CoV-2 S1 RBD	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000

### Coating Confirmation Testing Results:

Spot	Description	Result
1	SARS CoV-2 Spike	Pass
3	SARS CoV-2 Nucleocapsid	Pass
10	SARS CoV-2 S1 RBD	Pass



## CERTIFICATE OF ANALYSIS

Page 2 of 2

### 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
SARS CoV-2 Nucleocapsid	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 2 (Z0057218), 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		12 AUG 2022