



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX<sup>®</sup> COVID-19 Coronavirus Panel 3 Kit  
**Kit Catalog Numbers:** K15400-Series; K15401-Series; K15402-Series; K15399-Series  
**Kit Lot Number:** K0081889  
**Expiration Date:** 31 MAR 2023

### Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
Coronavirus Plate 3	Z0056799	2-8°C	31 MAR 2023
SULFO-TAG <sup>™</sup> Anti-Human IgG Antibody	D00V0005	2-8°C	29 FEB 2024
SULFO-TAG <sup>™</sup> Anti-Human IgM Antibody	D00V0011	2-8°C	31 AUG 2024
SULFO-TAG <sup>™</sup> Anti-Human IgA Antibody	D00V0010	2-8°C	31 JUL 2024
Reference Standard 1	A00V0004	≤ -70°C	30 JUN 2025
Serology Control 1.1	A00C0771	≤ -70°C	28 FEB 2026
Serology Control 1.2	A00C0772	≤ -70°C	28 FEB 2026
Serology Control 1.3	A00C0773	≤ -70°C	28 FEB 2026
SULFO-TAG <sup>™</sup> Human ACE2 Protein	D0081708	2-8°C	31 JUL 2023
ACE2 Calibration Reagent	A0080285	2-8°C	31 AUG 2024
MSD Blocker A Kit	Not Kit Specific	Room Temperature	See Label
Diluent 100	Not Kit Specific	2-8°C	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, IgM, IgA, and ACE2 kit.

### Plate Uniformity Testing Results:

Parameter	Precision			Uniformity	Signal
	CV of Intra-plate Averages	Average Intra-plate CV	Max Intra-plate CV		
SARS-CoV-2 Spike	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
HCoV-NL63 Spike	≤ 25%	≤ 15%	≤ 18%	Pass	1,500 – 1,000,000
SARS-CoV-2 Nucleocapsid	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-1 Spike	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
MERS-CoV Spike	≤ 25%	≤ 15%	≤ 18%	Pass	1,500 – 1,000,000
HCoV-HKU1 Spike	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
HCoV-OC43 Spike	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
HCoV-229E Spike	≤ 25%	≤ 15%	≤ 18%	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	SARS-CoV-2 Spike	Pass
2	HCoV-NL63 Spike	Pass
3	SARS-CoV-2 Nucleocapsid	Pass
4	SARS-CoV-1 Spike	Pass
6	MERS-CoV Spike	Pass
7	HCoV-HKU1 Spike	Pass
8	HCoV-OC43 Spike	Pass
9	HCoV-229E Spike	Pass
10	SARS-CoV-2 S1 RBD	Pass

### Functional Testing Results:

Sample Type	Calibrator	Controls		Samples	
Metric	Average Signal Ratio Test:Reference	Control Recovery	Control Conc. CV	Average % Difference Test:Reference	Slope Test:Reference
SARS-CoV-2 Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
HCoV-NL63 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-1 Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
MERS-CoV Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
HCoV-HKU1 Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
HCoV-OC43 Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
HCoV-229E Spike	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 Coronavirus Plate 3 (Z0056799), SULFO-TAG Anti-Human IgG Antibody (D00V0005), Reference Standard (A00V0004), and Serology Controls (A00C0731, A00C0732, A00C0733).

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		28 OCT 2021