



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® SARS-CoV-2 Panel 5 Kit  
**Kit Catalog Numbers:** K15429-Series; K15432-Series  
**Kit Lot Number:** K0081688  
**Expiration Date:** 31 DEC 2022

### Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
SARS-CoV-2 Plate 5	Z0056894	2-8°C	31 AUG 2023
SULFO-TAG™ Anti-Human IgG Antibody	D00V0005	2-8°C	29 FEB 2024
Reference Standard 1	A0080270	≤ -70°C	30 SEP 2025
Serology Control 1.1	A00C0731	≤ -70°C	30 JUN 2025
Serology Control 1.2	A00C0732	≤ -70°C	30 JUN 2025
Serology Control 1.3	A00C0733	≤ -70°C	30 JUN 2025
SULFO-TAG™ Human ACE2 Protein	D0081619	2-8°C	31 DEC 2022
ACE2 Calibration Reagent	A00V0001	2-8°C	31 DEC 2023
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, and ACE2 kit.

### Plate Uniformity Testing Results:

Parameter	Precision			Uniformity	Signal
	CV of Intra-plate Averages	Average Intra-plate CV	Max Intra-plate CV		
CoV-2 Spike	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
CoV-2 N	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Spike P.1	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Spike B.1.1.7	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
Spike B.1.351	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000

### Coating Confirmation Testing Results:

Spot	Description	Result
1	CoV-2 Spike	Pass
3	CoV-2 N	Pass
7	Spike P.1	Pass
8	Spike B.1.1.7	Pass
9	Spike B.1.351	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	Average % Difference Test:Reference	Slope Test:Reference
CoV-2 Spike	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
CoV-2 N	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Spike P.1	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Spike B.1.1.7	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
Spike B.1.351	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2

### Additional Comments:

Functional testing was executed using SARS-CoV-2 Plate 5 (Z0056894), SULFO-TAG Anti-Human IgG Antibody (D00V0005), Reference Standard 1 (A0080270), 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
Review/Approval	Karen Hamilla	Quality		25 MAR 2021