



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX[®] SARS-CoV-2 Panel 15 Kit
Kit Catalog Numbers: K15499-Series; K15500-Series; K15501-Series; K15502-Series
Kit Lot Number: K0081800
Expiration Date: 31 DEC 2022

Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
SARS-CoV-2 Plate 15	Z0057003	2-8°C	31 DEC 2023
SULFO-TAG [™] Anti-Human IgG Antibody	D00V0009	2-8°C	31 JUL 2024
SULFO-TAG [™] Anti-Human IgM Antibody	D00V0011	2-8°C	31 AUG 2024
SULFO-TAG [™] Anti-Human IgA Antibody	D00V0010	2-8°C	31 JUL 2024
Reference Standard 1	A0080286	≤ -70°C	28 FEB 2026
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 [™] Human ACE2 Protein	D0081619	2-8°C	31 DEC 2022
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, IgA, IgM, and 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 Spike (AY.1)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (AY.2)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.617.2+ΔY144)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.620)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.258.17)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.466.2)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.1.7+E484K)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.351.1)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
SARS-CoV-2 Spike (B.1.618)	≤ 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	SARS-CoV-2 Spike (AY.1)	Pass
3	SARS-CoV-2 Spike (AY.2)	Pass
4	SARS-CoV-2 Spike (B.1.617.2+ ΔY144)	Pass
5	SARS-CoV-2 Spike (B.1.620)	Pass
6	SARS-CoV-2 Spike (B.1.258.17)	Pass
7	SARS-CoV-2 Spike (B.1.466.2)	Pass
8	SARS-CoV-2 Spike (B.1.1.7+E484K)	Pass
9	SARS-CoV-2 Spike (B.1.351.1)	Pass
10	SARS-CoV-2 Spike (B.1.618)	Pass

Functional Testing Results:

Sample Type	Calibrator	Controls		Samples	
Metric	Average 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 Spike (AY.1)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (AY.2)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.617.2+ ΔY144)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.620)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.258.17)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.466.2)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.1.7+E484K)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.351.1)	N/A	N/A	N/A	N/A	N/A
SARS-CoV-2 Spike (B.1.618)	N/A	N/A	N/A	N/A	N/A

Additional Comments:

Functional testing was executed using SARS-CoV-2 Plate 15 (Z0057003), SULFO-TAG Anti-Human IgG Antibody (D00V0009), Reference Standard 1 (A0080286), and Serology Controls (A00C0771, A00C0772, A00C0773).

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		28 JUL 2021