



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX[®] SARS-CoV-2 Panel 22 Kit
Kit Catalog Numbers: K15559-Series; K15560-Series; K15561-Series; K15562-Series; K15563-Series; K15564-Series; K15565-Series
Kit Lot Number: K0082027
Expiration Date: 31 JUL 2023

Kit Components:

Description	Lot Number	Storage Temperature	Expiration Date
SARS-CoV-2 Plate 22	Z0057107	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	D0081708	2-8°C	31 JUL 2023
ACE2 Calibration Reagent 2	A0080327	2-8°C	31 OCT 2025
SULFO-TAG [™] Anti-Human IgG Antibody	D00V0019	2-8°C	30 APR 2025
SULFO-TAG [™] Anti-Human IgM Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG [™] Anti-Human IgA Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG [™] Anti-Mouse IgG Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG [™] Anti-Mouse IgM Antibody	Not Kit Specific	2-8°C	See Label
SULFO-TAG [™] 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		
RBD (Omicron) (B.1.1.529; BA.1)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (Beta) (B.1.351)	≤ 25%	≤ 15%	≤ 18%	Pass	1,500 – 1,000,000
RBD (Gamma) (P.1)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (Alpha) (B.1.1.7)	≤ 18%	≤ 10%	≤ 13%	Pass	1,500 – 1,000,000
RBD (Delta) (B.1.617.2)	≤ 18%	≤ 10%	≤ 13%	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 (Omicron) (B.1.1.529; BA.1)	Pass
2	RBD (Beta) (B.1.351)	Pass
4	RBD (Gamma) (P.1)	Pass
6	RBD (Alpha) (B.1.1.7)	Pass
9	RBD (Delta) (B.1.617.2)	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 (Omicron) (B.1.1.529; BA.1)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (Beta) (B.1.351)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (Gamma) (P.1)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (Alpha) (B.1.1.7)	50 – 200%	70 – 130%	< 20%	± 25%	0.80 – 1.2
RBD (Delta) (B.1.617.2)	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 22 (Z0057107), SULFO-TAG Anti-Human IgG Antibody (D00V0019), 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		08 JUN 2022