



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Test Sample 2

Catalog Number: C1ADM-1

Lot Number: A00C0735

Expiration Date: 30 JUN 2025

Storage Temperature: ≤ -70 °C

Concentration Assignment:

Analyte	Concentration of IgG (viald)	Concentration of IgA (viald)	Concentration of IgM (viald)	Unit of Measure
SARS-CoV-2 Spike	12500	1130	2450	AU/mL
HCoV-NL63 Spike	4150	3645		AU/mL
SARS-CoV-2 Nucleocapsid	11100	2000	1620	AU/mL
SARS-CoV-1 Spike	2210	695		AU/mL
SARS-CoV-2 S1 NTD	NA	328		AU/mL
HCoV-HKU1 Spike	17000	3295		AU/mL
HCoV-OC43 Spike	82500	11200		AU/mL
HCoV-229E Spike	25100	5450		AU/mL
SARS-CoV-2 S1 RBD	4880	680	1740	AU/mL
MERS-CoV Spike	4160	1600		AU/mL
Flu B/Brisbane HA	76900	21400		AU/mL
Flu A/Shanghai H7	6200	1890		AU/mL
Flu A/Michigan H1	52800	11800		AU/mL
RSV Pre-Fusion F	183000	35350		AU/mL
Flu A/HongKong H3	63100	18700		AU/mL
Flu B/Phuket HA	91500	19350		AU/mL

NA : not assigned

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		17 NOV 2020