



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Test Sample 1

**Catalog Number:** C1ADL-1

**Lot Number:** A00C0734

**Expiration Date:** 30 JUN 2025

**Storage Temperature:**  $\leq -70$  °C

### Concentration Assignment:

Assay	Concentration of IgG (viald)	Concentration of IgA (viald)	Concentration of IgM (viald)	Unit of Measure
SARS-CoV-2 Spike	41300	5750	6200	AU/mL
HCoV-NL63 Spike	5230	1045		AU/mL
SARS-CoV-2 Nucleocapsid	104000	4910	7190	AU/mL
SARS-CoV-1 Spike	12200	2775		AU/mL
SARS-CoV-2 S1 NTD	817	436		AU/mL
HCoV-HKU1 Spike	28300	4470		AU/mL
HCoV-OC43 Spike	128000	11950		AU/mL
HCoV-229E Spike	32200	2665		AU/mL
SARS-CoV-2 S1 RBD	11200	5250	3510	AU/mL
MERS-CoV Spike	2780	1730		AU/mL
Flu B/Brisbane HA	90000	14250		AU/mL
Flu A/Shanghai H7	11600	2485		AU/mL
Flu A/Michigan H1	75400	9100		AU/mL
RSV Pre-Fusion F	205000	18850		AU/mL
Flu A/HongKong H3	70700	7600		AU/mL
Flu B/Phuket HA	281000	19950		AU/mL

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		11 NOV 2020