



## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Reference Standard 2

**Catalog Number:** C00AUB-2

**Lot Number:** A0080403

**Storage Temperature:** ≤ -70 °C

**Expiration Date:** 30 JUN 2034

**Values are provided in two units** – MSD Arbitrary Units/mL (AU/mL) and WHO Biological Arbitrary Units/mL (BAU/mL) (NIBSC 20/136)

WHO/NIBSC Units: MSD Units				
Antigen	Isotype	Vialed	1:10	Unit of Measure
SARS-CoV-2 Spike	IgG	11.4	1.14	BAU/mL
SARS-CoV-2 Spike	IgA	14.1		BAU/mL
SARS-CoV-2 N	IgG	2.09	0.209	BAU/mL
SARS-CoV-2 N	IgA	6.88		BAU/mL

### Concentration Assignment of IgG Antibodies:

**IgG:** Use at 1:10 for all plate types

Assay	Conc. of IgG (vialed)	Conc. of IgG (1:10)	Unit of Measure
SARS-CoV-2 Spike	1270	127	AU/mL
SARS-CoV-2 N	885	88.5	AU/mL
SARS-CoV-2 Spike (B.1.1.529; BA.1)	291	29.1	AU/mL
SARS-CoV-2 Spike (B.1.1.7)	899	89.9	AU/mL
SARS-CoV-2 Spike (B.1.351)	775	77.5	AU/mL
SARS-CoV-2 Spike (B.1.617.2; AY.4)	901	90.1	AU/mL
SARS-CoV-2 Spike (BA.2)	464	46.4	AU/mL
SARS-CoV-2 Spike (BA.2.12.1)	455	45.5	AU/mL
SARS-CoV-2 Spike (BA.2.75)	272	27.2	AU/mL
SARS-CoV-2 Spike (BA.2.86)	336	33.6	AU/mL
SARS-CoV-2 Spike (BA.5)	635	63.5	AU/mL
SARS-CoV-2 Spike (EG.5.1)	318	31.8	AU/mL
SARS-CoV-2 Spike (FL.1.5.1)	335	33.5	AU/mL
SARS-CoV-2 Spike (HV.1)	366	36.6	AU/mL
SARS-CoV-2 Spike (JD.1.1)	365	36.5	AU/mL
SARS-CoV-2 Spike (JN.1)	299	29.9	AU/mL
SARS-CoV-2 Spike (XBB.1.16)	420	42.0	AU/mL
SARS-CoV-2 Spike (XBB.1.16.6)	286	28.6	AU/mL
SARS-CoV-2 Spike (XBB.1.5)	289	28.9	AU/mL
SARS-CoV-2 Spike (XBB.2.3)	478	47.8	AU/mL



### CERTIFICATE OF ANALYSIS

#### Concentration Assignment of IgA Antibodies:

IgA: Use as vialled for all plate types

Assay	Conc. of IgA (vialled)	Unit of Measure
SARS-CoV-2 Spike	227	AU/mL
SARS-CoV-2 N	62.0	AU/mL
SARS-CoV-2 Spike (B.1.1.529; BA.1)	62.6	AU/mL
SARS-CoV-2 Spike (B.1.1.7)	158	AU/mL
SARS-CoV-2 Spike (B.1.351)	148	AU/mL
SARS-CoV-2 Spike (B.1.617.2; AY.4)	212	AU/mL
SARS-CoV-2 Spike (BA.2)	143	AU/mL
SARS-CoV-2 Spike (BA.2.12.1)	128	AU/mL
SARS-CoV-2 Spike (BA.2.75)	70.7	AU/mL
SARS-CoV-2 Spike (BA.2.86)	67.7	AU/mL
SARS-CoV-2 Spike (BA.5)	165	AU/mL
SARS-CoV-2 Spike (EG.5.1)	78.9	AU/mL
SARS-CoV-2 Spike (FL.1.5.1)	77.3	AU/mL
SARS-CoV-2 Spike (HV.1)	73.0	AU/mL
SARS-CoV-2 Spike (JD.1.1)	78.7	AU/mL
SARS-CoV-2 Spike (JN.1)	57.6	AU/mL
SARS-CoV-2 Spike (XBB.1.16)	73.0	AU/mL
SARS-CoV-2 Spike (XBB.1.16.6)	58.9	AU/mL
SARS-CoV-2 Spike (XBB.1.5)	72.5	AU/mL
SARS-CoV-2 Spike (XBB.2.3)	72.5	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
<b>Review/Approval</b>	Karen Hamilla	Quality		28 OCT 2024