



**CERTIFICATE OF ANALYSIS**

**Product Description:** ACE2 Calibration Reagent 3

**Catalog Number:** C01APE-2

**Lot Number:** A0080375

**Storage Temperature:** ≤-70°C

**Expiration Date:** 31 JAN 2027

**Concentration Assignment of Neutralizing Antibodies:**

Assay	Alternate Name	Stock Concentration	Unit of Measure
<b>SARS-CoV-2</b>			
SARS-CoV-2 S1 RBD		327	IAU/mL
SARS-CoV-2 Spike		331	IAU/mL
SARS-CoV-2 S1 RBD (AY.3)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.4)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.4.2)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.5)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.6)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.7)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.12)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (AY.14)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (B.1.1.529; BA.1)	RBD (Omicron; BA.1)	47.5	IAU/mL
SARS-CoV-2 S1 RBD (B.1.1.7)	N501Y; RBD (Alpha)	232	IAU/mL
SARS-CoV-2 S1 RBD (B.1.351)	K417N, E484K, N501Y; RBD (Beta)	169	IAU/mL
SARS-CoV-2 S1 RBD (B.1.351.1)	K417N, E484K, N501Y; RBD (Beta)	169	IAU/mL
SARS-CoV-2 S1 RBD (B.1.617.2)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (B.1.617.2+ΔY144)	L452R, T478K; RBD (Delta)	214	IAU/mL
SARS-CoV-2 S1 RBD (BA.2)	RBD (Omicron; BA.2)	65.7	IAU/mL
SARS-CoV-2 S1 RBD (BA.2.12.1)	RBD (BA.2.12.1)	75.7	IAU/mL
SARS-CoV-2 S1 RBD (BA.2.75)	RBD (BA.2.75)	52.5	IAU/mL
SARS-CoV-2 S1 RBD (BA.2.75.2)	RBD (BA.2.75.2)	60.9	IAU/mL
SARS-CoV-2 S1 RBD (BA.4; BA.5)	RBD (BA.4; BA.5)	98.5	IAU/mL
SARS-CoV-2 S1 RBD (BA.4.6; BF.7)	RBD (BA.4.6;BF.7)	104	IAU/mL
SARS-CoV-2 S1 RBD (BQ.1)	RBD (BQ.1)	84.9	IAU/mL
SARS-CoV-2 S1 RBD (BQ.1.1)	RBD (BQ.1.1)	83.6	IAU/mL
SARS-CoV-2 S1 RBD (XBB.1)	RBD (XBB.1)	82.4	IAU/mL
SARS-CoV-2 Spike (B.1.1.529; BA.1)		65.0	IAU/mL
SARS-CoV-2 Spike (B.1.1.7)		259	IAU/mL
SARS-CoV-2 Spike (B.1.351)		168	IAU/mL
SARS-CoV-2 Spike (B.1.617.2)		241	IAU/mL
SARS-CoV-2 Spike (BA.2)		92.0	IAU/mL
SARS-CoV-2 Spike (BA.2.12.1)		103	IAU/mL
SARS-CoV-2 Spike (BA.2.75)		80.0	IAU/mL
SARS-CoV-2 Spike (BA.2.75.2)		86.2	IAU/mL



### CERTIFICATE OF ANALYSIS

SARS-CoV-2 Spike (BA.2.86)		59.6	IAU/mL
SARS-CoV-2 Spike (BA.4.6)		131	IAU/mL
SARS-CoV-2 Spike (BA.5)		124	IAU/mL
SARS-CoV-2 Spike (BF.7)		116	IAU/mL
SARS-CoV-2 Spike (BN.1)		86.1	IAU/mL
SARS-CoV-2 Spike (BQ.1)		87.1	IAU/mL
SARS-CoV-2 Spike (BQ.1.1)		81.8	IAU/mL
SARS-CoV-2 Spike (EG.5.1)		51.3	IAU/mL
SARS-CoV-2 Spike (FL.1.5.1)		54.2	IAU/mL
SARS-CoV-2 Spike (HV.1)		29.7	IAU/mL
SARS-CoV-2 Spike (JD.1.1)		24.5	IAU/mL
SARS-CoV-2 Spike (JN.1)		24.9	IAU/mL
SARS-CoV-2 Spike (XBB.1)		87.7	IAU/mL
SARS-CoV-2 Spike (XBB.1.16.1)		90.7	IAU/mL
SARS-CoV-2 Spike (XBB.1.16)		81.7	IAU/mL
SARS-CoV-2 Spike (XBB.1.16.6)		52.1	IAU/mL
SARS-CoV-2 Spike (XBB.2.3)		90.4	IAU/mL
SARS-CoV-2 Spike (XBB.1.5)		70.6	IAU/mL

**Note:** IAU/mL refers to MSD Inhibitory Antibody Units/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	Signature	Date
<b>Review/Approval</b>	Karen Hamilla		06 May 2024