



CERTIFICATE OF ANALYSIS

Product Description: SARS-CoV-2 Spike Calibrator
Catalog Number: C00ADJ-2
Lot Number: A00S0116
Expiration Date: 31 MAR 2027
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SARS-CoV-2 Spike	7520000	fg/ml

Dilution to Top of Curve: 20 X

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	<i>Karen Hamilla</i>	17 OCT 2024