



CERTIFICATE OF ANALYSIS

Product Description: SARS-CoV-2 N Calibrator
Catalog Number: C00ADH-2
Lot Number: A00S0104
Expiration Date: 30 SEP 2026
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SARS CoV-2 N	23200000	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		10 JUL 2024