

**CERTIFICATE OF ANALYSIS**

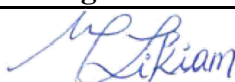
Product Description: SARS-CoV-2 Spike Calibrator
Catalog Number: C00ADJ-2
Lot Number: A00S0145
Expiration Date: 30 APR 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
SARS-CoV-2 Spike	7680000	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	Michelle Likiam	Quality		26 DEC 2025