

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX SARS-CoV-2 Spike Kit
Lot Number: K00S0290
Expiration Date: 31 MAR 2027
Catalog Number: K150ADJS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin SARS-CoV-2 Spike Antibody	B00T0269	2-8°C	28 FEB 2029
TURBO-BOOST SARS-CoV-2 Spike Antibody	D00T0269	2-8°C	31 MAR 2029
Blocker S1	R0140014	≤ -10°C	30 SEP 2027
Diluent 61	M0610005	≤ -10°C	30 NOV 2027
S-PLEX Enhance E1	Y0550023	≤ -10°C	31 AUG 2027
S-PLEX Enhance E2	Y0560022	≤ -10°C	31 JUL 2027
SARS-CoV-2 Spike Calibrator	A00S0145	≤ -70°C	30 APR 2028
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580032	≤ -70°C	31 AUG 2027
S-PLEX Detect D2	Y0590025	≤ -70°C	30 NOV 2027
S-PLEX Enhance E3	Y0570025	≤ -70°C	31 AUG 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot Number	Name	Calibrator	
		Recovery	CV
1	SARS-CoV-2 Spike	75% - 125%	< 25%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
SARS-CoV-2 Spike	175 - 2800	284	fg/mL	75% - 125%	< 25%	47.4	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
SARS-CoV-2 Spike	20X Calibrator	7680000	fg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 FEB 2026