



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX SARS-CoV-2 Spike Kit
Lot Number: K00S0328
Expiration Date: 30 NOV 2027
Catalog Number: K150ADJS-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Biotin SARS-CoV-2 SpikeAntibody	B00T0298	2-8°C	31 OCT 2029
TURBO-BOOST SARS-CoV-2 Spike Antibody	D00T0298	2-8°C	31 OCT 2029
Blocker S1	R0140022	≤ -10°C	31 OCT 2028
Diluent 61	M0610005	≤ -10°C	30 NOV 2027
S-PLEX Enhance E1	Y0550028	≤ -10°C	31 MAR 2028
S-PLEX Enhance E2	Y0560028	≤ -10°C	31 MAY 2028
SARS-CoV-2 Spike Calibrator	A00S0145	≤ -70°C	30 APR 2028
S-PLEX Coating Reagent C1	521331	≤ -70°C	31 OCT 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580037	≤ -70°C	30 APR 2028
S-PLEX Detect D2	Y0590029	≤ -70°C	31 MAY 2028
S-PLEX Enhance E3	Y0570030	≤ -70°C	31 MAR 2028

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

Precision, Accuracy, and Sensitivity Specifications:

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

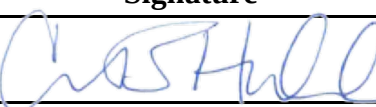
LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
SARS-CoV-2 Spike	175 - 2800	190	fg/mL	75% - 125%	< 25%	71.9	fg/mL

Stock Concentration of Calibrator:

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

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		09 JUL 2026