

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human IL-6 Kit
Lot Number: K00S0296
Expiration Date: 30 JUN 2027
Catalog Number: K151B3S-Series; K156B3S-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 Human IL-6 Antibody	B00T0274	2-8°C	30 APR 2029
Human IL-6 Calibrator	A00S0143	2-8°C	30 SEP 2030
TURBO-BOOST Human IL-6 Antibody	D00T0274	2-8°C	31 MAY 2029
Blocker S1	R0140017	≤ -10°C	29 FEB 2028
Diluent 2	M0020133	≤ -10°C	30 JUN 2027
Diluent 3	M0030145	≤ -10°C	29 FEB 2028
S-PLEX Enhance E1	Y0550026	≤ -10°C	31 JAN 2028
S-PLEX Enhance E2	Y0560026	≤ -10°C	29 FEB 2028
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580036	≤ -70°C	30 APR 2028
S-PLEX Detect D2	Y0590028	≤ -70°C	29 FEB 2028
S-PLEX Enhance E3	Y0570028	≤ -70°C	31 JAN 2028

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	IL-6	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
IL-6	0.325 - 5.20	5.08	fg/mL	80% - 120%	< 20%	0.339	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-6	120X Calibrator	744000	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	Karen Hamilla	Quality		19 MAR 2026