



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-1 β Kit
Lot Number: K00S0324
Expiration Date: 31 MAR 2028
Catalog Number: K151ADSS-Series; K156ADSS-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-1 β Antibody	B00T0294	2-8°C	30 NOV 2029
Human IL-1 β Calibrator	A00S0155	2-8°C	30 APR 2031
TURBO-BOOST Human IL-1 β Antibody	D00T0294	2-8°C	31 OCT 2029
Blocker S1	R0140023	$\leq -10^{\circ}\text{C}$	30 NOV 2028
Diluent 58	M0580019	$\leq -10^{\circ}\text{C}$	30 NOV 2028
S-PLEX Enhance E1	Y0550028	$\leq -10^{\circ}\text{C}$	31 MAR 2028
S-PLEX Enhance E2	Y0560028	$\leq -10^{\circ}\text{C}$	31 MAY 2028
S-PLEX Coating Reagent C1	521331	$\leq -70^{\circ}\text{C}$	31 OCT 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580038	$\leq -70^{\circ}\text{C}$	31 AUG 2028
S-PLEX Detect D2	Y0590030	$\leq -70^{\circ}\text{C}$	30 JUN 2028
S-PLEX Enhance E3	Y0570031	$\leq -70^{\circ}\text{C}$	30 JUN 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	IL-1 β	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-1 β	48.7 - 780	84.6	fg/mL	80% - 120%	< 20%	15.0	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-1 β	1X Calibrator	111000	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		17 JUL 2026