



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-17A Kit
Lot Number: K00S0312
Expiration Date: 31 MAR 2028
Catalog Number: K151C3S-Series; K156C3S-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-17A Antibody	B00T0283	2-8°C	30 SEP 2029
Human IL-17A Calibrator	A00S0151	2-8°C	28 FEB 2031
TURBO-BOOST Human IL-17A Antibody	D00T0283	2-8°C	31 AUG 2029
Blocker S1	R0140022	≤ -10°C	31 OCT 2028
Diluent 3	M0030149	≤ -10°C	31 AUG 2028
Diluent 43	M0430060	≤ -10°C	30 JUN 2028
S-PLEX Enhance E1	Y0550028	≤ -10°C	31 MAR 2028
S-PLEX Enhance E2	Y0560028	≤ -10°C	31 MAY 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	IL-17A	75% - 125%	< 25%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-17A	7.50 - 120	78.7	fg/mL	75% - 125%	< 25%	20.0	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-17A	10X Calibrator	2820000	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	Michelle Likiam	Quality		14 JUL 2026