



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IL-29/IFN- λ 1 Kit
Lot Number: K00S0280
Expiration Date: 31 OCT 2027
Catalog Number: K151AXCS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human IL-29/IFN- λ 1 Antibody	B00T0260	2-8°C	30 NOV 2028
TURBO-BOOST Human IL-29/IFN- λ 1 Antibody	D00T0260	2-8°C	31 DEC 2028
Blocker S1	R0140015	$\leq -10^{\circ}\text{C}$	31 OCT 2027
S-PLEX Enhance E1	Y0550024	$\leq -10^{\circ}\text{C}$	31 OCT 2027
S-PLEX Enhance E2	Y0560024	$\leq -10^{\circ}\text{C}$	31 DEC 2027
Human IL-29/IFN- λ 1 Calibrator	A00S0135	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Coating Reagent C1	521329	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580034	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Detect D2	Y0590026	$\leq -70^{\circ}\text{C}$	31 DEC 2027
S-PLEX Enhance E3	Y0570027	$\leq -70^{\circ}\text{C}$	31 DEC 2027

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-29/IFN- λ 1	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IL-29/IFN- λ 1	22.7 - 726	182	fg/mL	80% - 120%	< 20%	10.4	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-29/IFN- λ 1	20X Calibrator	3700000	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		02 SEP 2025