



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human IFN- ω Kit
Lot Number: K00S0320
Expiration Date: 29 FEB 2028
Catalog Number: K151AXDS-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 IFN- ω Antibody	B00T0291	2-8°C	31 JUL 2029
TURBO-BOOST Human IFN- ω Antibody	D00T0291	2-8°C	31 JUL 2029
Blocker S1	R0140018	$\leq -10^{\circ}\text{C}$	30 APR 2028
S-PLEX Enhance E1	Y0550027	$\leq -10^{\circ}\text{C}$	29 FEB 2028
S-PLEX Enhance E2	Y0560026	$\leq -10^{\circ}\text{C}$	29 FEB 2028
Human IFN- ω Calibrator	A00S0149	$\leq -70^{\circ}\text{C}$	31 AUG 2028
S-PLEX Coating Reagent C1	521330	$\leq -70^{\circ}\text{C}$	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG TM Reagent	Y0580036	$\leq -70^{\circ}\text{C}$	30 APR 2028
S-PLEX Detect D2	Y0590028	$\leq -70^{\circ}\text{C}$	29 FEB 2028
S-PLEX Enhance E3	Y0570029	$\leq -70^{\circ}\text{C}$	29 FEB 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	IFN- ω	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
IFN- ω	6.88 - 220	27.4	fg/mL	80% - 120%	< 20%	10.6	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN- ω	20X Calibrator	1158000	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		06 AUG 2026