



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX® Human IFN-γ Kit

Catalog Number: K151X9S-Series

Lot Number: K00S0032

Expiration Date: 31 AUG 2022

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IFN-γ	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IFN-γ	3.96 - 63.4	15.85	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IFN-γ	5.26	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IFN-γ	1X Calibrator	25400	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IFN-γ Antibody	D00T0037A	2-8°C	30APR 2024
Biotin Human IFN-γ Antibody	B00T0037A	2-8°C	29 FEB 2024
Human IFN-γ Calibrator	A00S0026	2-8°C	30 SEP 2025
Diluent 58	M0580001	≤ -10°C	28 FEB 2023
S-PLEX Coating Reagent C1	314374	≤ -70°C	28 FEB 2023
Blocker S1	R0140003	≤ -10°C	31AUG 2022
S-PLEX Enhance E1	Y0550004	≤ -10°C	28 FEB 2023
S-PLEX Enhance E2	Y0560004	≤ -10°C	31 JAN 2023
S-PLEX Enhance E3	Y0570004	≤ -70°C	31 JAN 2023
S-PLEX Detect D1	Y0580006	≤ -70°C	31MAR 2023
S-PLEX Detect D2	Y0590003	≤ -70°C	31AUG 2022

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 FEB 2021