



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human IL-12p70 Kit

Catalog Number: K151G3S - Series

Lot Number: K00S0017

Expiration Date: 31 JAN 2022

Precision, Accuracy, and Sensitivity Specifications:

Calibrator		
Name	Recovery	CV
IL-12p70	80% - 120%	< 20%

LLOQ					
Name	Range Tested	Observed	Units	Recovery	CV
IL-12p70	0.85 - 13.6	3.4	fg/mL	80% - 120%	< 20%

LLOD		
Name	Observed	Units
IL-12p70	2.58	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-12p70	15X Calibrator	496000	fg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0023A	2-8°C	30 SEP 2023
Biotin Human IL-12p70 Antibody	B00T0023A	2-8°C	31 DEC 2023
Human IL-12p70 Calibrator	A00S0024	2-8°C	31 JUL 2025
Diluent 2	M0020107	≤ -10°C	31 JAN 2022
Diluent 3	M0030091	≤ -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	314374	≤ -70°C	28 FEB 2023
Blocker S1	R0140003	≤ -10°C	31 AUG 2022
S-PLEX Enhance E1	Y0550003	≤ -10°C	30 SEP 2022
S-PLEX Enhance E2	Y0560005	≤ -10°C	30 SEP 2022
S-PLEX Enhance E3	Y0570003	≤ -70°C	31 DEC 2022
S-PLEX Detect D1	Y0580003	≤ -70°C	30 NOV 2022
S-PLEX Detect D2	Y0590003	≤ -70°C	31 AUG 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		21 AUG 2020