



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup> Human IL-12p70 Kit

**Catalog Number:** K151G3S-Series

**Lot Number:** K00S0005A

**Expiration Date:** 30 APR 2021

### Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ				
	Range Tested	Observed	Units	Recovery	CV
IL-12p70	0.85 - 13.6	6.8	fg/mL	80% - 120%	< 20%

Name	LLOD	
	Observed	Units
IL-12p70	2.34	fg/mL

### Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-12p70	15X Calibrator	523000	fg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human IL-12p70 Antibody	D00T0009A	2-8°C	31 OCT 2022
Biotin Human IL-12p70 Antibody	B00T0009A	2-8°C	31 OCT 2022
Human IL-12p70 Calibrator	A00S0019	2-8°C	30 SEP 2024
Diluent 2	M0020107	≤ -10°C	31 JAN 2022
Diluent 3	M0030091	≤ -10°C	31 JAN 2022
S-PLEX Coating Reagent C1	265437	≤ -70°C	30 SEP 2021
Blocker S1	R0140001	≤ -10°C	31 MAY 2021
S-PLEX Enhance E1	Y0550001	≤ -10°C	30 APR 2021
S-PLEX Enhance E2	Y0560002	≤ -10°C	31 DEC 2021
S-PLEX Enhance E3	Y0570001	≤ -70°C	31 MAY 2021
S-PLEX Detect D1	Y0580001	≤ -70°C	30 APR 2021
S-PLEX Detect D2	Y0590001	≤ -70°C	30 APR 2021

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		19 MAY 2020