



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: S-PLEX[®]Neurology Panel 1 (human) Kit

Catalog Number: K15640S-Series; K15639S-Series; K151AKGS-Series; K156AKGS-Series

Lot Number: K00S0112

Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human GFAP Antibody	D00T0117	2-8°C	30 APR 2026
TURBO-BOOST [™] Hu Neurofilament L Antibody	D00T0118	2-8°C	31 MAY 2026
TURBO-BOOST [™] Human Tau Antibody	D00T0119	2-8°C	30 APR 2026
S-PLEX Neurology Panel 1 (human) Coating Solution	429070	≤ -70°C	30 APR 2026
S-PLEX Neurology Panel 1 (human) Calibrator Blend	A00S0074	2-8°C	31 OCT 2027
Diluent 63	M0630002	≤ -10°C	30 APR 2025
Blocker S1	R0010811	≤ -10°C	28 FEB 2025
Blocker S2	420803	≤ -10°C	30 SEP 2027
S-PLEX Enhance E1	Y0550010	≤ -10°C	31 MAY 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 MAR 2025
S-PLEX Enhance E3	Y0570011	≤ -70°C	31 MAR 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580015	≤ -70°C	30 APR 2025
S-PLEX Detect D2	Y0590011	≤ -70°C	31 MAR 2025

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator	
		Recovery	CV
1	GFAP	80% - 120%	< 20%
3	NF-L	80% - 120%	< 20%
10	Tau (Total)	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
GFAP	53.0 - 12500	391	fg/mL	80% - 120%	< 25%	252	fg/mL
NF-L	193 - 45900	23900	fg/mL	80% - 120%	< 25%	2480	fg/mL
Tau (total)	15.0 - 4160	104	fg/mL	80% - 120%	< 25%	49.8	fg/mL




CERTIFICATE OF ANALYSIS

Page 2 of 2

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GFAP	1X Calibrator	781000	fg/mL
NF-L	1X Calibrator	2990000	fg/mL
Tau (total)	1X Calibrator	208000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		27 DEC 2022