



CERTIFICATE OF ANALYSIS

Page 1 of 2

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

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

Lot Number: K00S0130

Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human GFAP Antibody	D00T0151	2-8°C	31 Oct 2026
TURBO-BOOST [™] Hu Neurofilament L Antibody	D00T0152	2-8°C	30 Nov 2026
TURBO-BOOST [™] Human Tau Antibody	D00T0153	2-8°C	30 Nov 2026
S-PLEX Neurology Panel 1(human) Coating Solution	417648	≤ -70°C	28 Feb 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	437541	≤ -10°C	31 Dec 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	Y0570012	≤ -70°C	31 May 2025
S-PLEX Detect D1	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:

Name	Calibrator	
	Recovery	CV
GFAP	80% - 120%	< 20%
NF-L	80% - 120%	< 20%
Tau (total)	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
GFAP	53 - 12500	391	fg/mL	80% - 120%	< 25%	211	fg/mL
NF-L	193 - 45900	5980	fg/mL	80% - 120%	< 25%	2630	fg/mL
Tau (total)	15 - 4160	104	fg/mL	80% - 120%	< 25%	31.4	fg/mL

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



CERTIFICATE OF ANALYSIS

Page 2 of 2

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

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		10 JUL 2023