

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Neurology Panel 1 (human) Kit
Catalog Number: K15639S-Series; K15640S-Series; K151AKGS-Series; K156AKGS-Series;
 K151AMPS-Series; K156AMPS-Series; K151APSS-Series; K156APSS-Series
Lot Number: K00S0157
Expiration Date: 31 OCT 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Hu Neurofilament L Antibody	D00T0167	2-8°C	31 JAN 2027
TURBO-BOOST [™] Human GFAP Antibody	D00T0166	2-8°C	31 MAR 2027
TURBO-BOOST [™] Human Tau Antibody	D00T0168	2-8°C	28 FEB 2027
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
Blocker S2	438077	≤ -10°C	30 SEP 2028
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 NOV 2025
S-PLEX Enhance E2	Y0560014	≤ -10°C	31 OCT 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580019	≤ -70°C	31 DEC 2025
S-PLEX Detect D2	Y0590015	≤ -70°C	31 JAN 2026
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 2025
S-PLEX Neurology Panel 1 (human) Calibrator	A00S0089	≤ -70°C	31 JAN 2026
S-PLEX Neurology Panel 1(human) Coating Solution	452377	≤ -70°C	30 NOV 2026

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
GFAP	85.0 - 20300	880	fg/mL	80% - 120%	< 25%	165	fg/mL
NF-L	420 - 99800	2170	fg/mL	80% - 120%	< 25%	532	fg/mL
Tau (total)	17.5 - 4160	38.3	fg/mL	80% - 120%	< 25%	21.9	fg/mL



CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GFAP	20X Calibrator	17600000	fg/mL
NF-L	20X Calibrator	86600000	fg/mL
Tau (total)	20X Calibrator	3060000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 FEB 2024