



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** S-PLEX<sup>®</sup>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:** K00S0159

**Expiration Date:** 31 Oct 2025

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST <sup>™</sup> Human GFAP Antibody	D00T0151	2-8°C	31 Oct 2026
TURBO-BOOST <sup>™</sup> Hu Neurofilament L Antibody	D00T0152	2-8°C	30 Nov 2026
TURBO-BOOST <sup>™</sup> Human Tau Antibody	D00T0153	2-8°C	30 Nov 2026
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
Blocker S1	R0010832	≤ -10°C	31 Oct 2025
Blocker S2	437541	≤ -10°C	31 Dec 2027
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 Nov 2025
S-PLEX Enhance E2	Y0560014	≤ -10°C	31 Oct 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 Oct 2025
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> reagent	Y0580019	≤ -70°C	31 Dec 2025
S-PLEX Detect D2	Y0590015	≤ -70°C	31 Jan 2026

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	85.0 - 20300	220	fg/mL	80% - 120%	< 25%	179	fg/mL
NF-L	420 - 99800	4330	fg/mL	80% - 120%	< 25%	715	fg/mL
Tau (total)	17.5 - 4160	38.3	fg/mL	80% - 120%	< 25%	25.2	fg/mL

#### 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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 JAN 2024