

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX<sup>®</sup> Neurology Panel 1 (human) Kit  
**Lot Number:** K00S0211  
**Expiration Date:** 30 JUN 2026  
**Catalog Number:** K151APSS-Series, K151AMPS-Series, K151AKGS-Series, K156APSS-Series, K156AMPS-Series, K15639S-Series, K15640S-Series, K156AKGS-Series

**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	R0140007	≤ -10°C	30 JUN 2026
Blocker S2	438076	≤ -10°C	31 MAY 2028
S-PLEX Enhance E1	Y0550017	≤ -10°C	31 AUG 2026
S-PLEX Enhance E2	Y0560019	≤ -10°C	30 NOV 2026
S-PLEX Detect D1 with TURBO-TAG <sup>™</sup> Reagent	Y0580025	≤ -70°C	30 SEP 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570020	≤ -70°C	31 OCT 2026
S-PLEX Neurology Panel 1 (human) Calibrator Blend	A00S0113	≤ -70°C	31 JAN 2027
S-PLEX Neurology Panel 1 (human) Coating Solution	496767	≤ -70°C	31 DEC 2027

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

**Precision, Accuracy, and Sensitivity Specifications:**

		Calibrator					
Spot Number	Name	Recovery	CV				
1	GFAP	80% - 120%	< 20%				
3	NF-L	80% - 120%	< 20%				
10	Tau (total)	80% - 120%	< 20%				
		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
GFAP	85.0 - 20300	446	fg/mL	80% - 120%	< 25%	172	fg/mL
NF-L	420 - 99800	2270	fg/mL	80% - 120%	< 25%	656	fg/mL
Tau (total)	17.5 - 4160	80.0	fg/mL	80% - 120%	< 25%	30.4	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
GFAP	20X Calibrator	17840000	fg/mL
NF-L	20X Calibrator	90600000	fg/mL
Tau (total)	20X Calibrator	3200000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		07 OCT 2024