

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Neurology Panel 1 (human) Kit
Lot Number: K00S0244
Expiration Date: 31 JUL 2026
Catalog Number: K15639S-Series; K15640S-Series; K151AKGS-Series; K156AKGS-Series; K151AMPS-Series; K156AMPS-Series; K151APSS-Series; K156APSS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST Hu Neurofilament L Antibody	D00T0207	2-8°C	30 NOV 2027
TURBO-BOOST Human GFAP Antibody	D00T0206	2-8°C	31 AUG 2027
TURBO-BOOST Human Tau Antibody	D00T0208	2-8°C	31 OCT 2027
Blocker S1	R0140008	≤ -10°C	31 JUL 2026
Blocker S2	476747	≤ -10°C	30 APR 2029
S-PLEX Enhance E1	Y0550020	≤ -10°C	30 APR 2027
S-PLEX Enhance E2	Y0560020	≤ -10°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580029	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590021	≤ -70°C	31 MAR 2027
S-PLEX Enhance E3	Y0570022	≤ -70°C	30 APR 2027
S-PLEX Neurology Panel 1 (human) Calibrator Blend	A00S0106	≤ -70°C	30 SEP 2026
S-PLEX Neurology Panel 1 (human) Coating Solution	485774	≤ -70°C	31 AUG 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	464	fg/mL	80% - 120%	< 25%	210	fg/mL
NF-L	420 - 99800	2120	fg/mL	80% - 120%	< 25%	612	fg/mL
Tau(total)	17.5 - 4160	38.8	fg/mL	80% - 120%	< 25%	29.2	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GFAP	20X Calibrator	18540000	fg/mL
NF-L	20X Calibrator	84800000	fg/mL
Tau(total)	20X Calibrator	3100000	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		26 DEC 2024