



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Neurology Panel 1 (human) Kit
Lot Number: K00S0268
Expiration Date: 30 JUN 2027
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	D0TB0030	2-8°C	30 NOV 2028
TURBO-BOOST Human GFAP Antibody	D0TB0029	2-8°C	31 OCT 2028
TURBO-BOOST Human Tau Antibody	D0TB0031	2-8°C	30 NOV 2028
Blocker S1	R0140014	≤ -10°C	30 SEP 2027
Blocker S2	476749	≤ -10°C	31 OCT 2029
S-PLEX Enhance E1	Y0550022	≤ -10°C	31 JUL 2027
S-PLEX Enhance E2	Y0560022	≤ -10°C	31 JUL 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580031	≤ -70°C	31 JUL 2027
S-PLEX Detect D2	Y0590024	≤ -70°C	31 AUG 2027
S-PLEX Enhance E3	Y0570024	≤ -70°C	30 JUN 2027
S-PLEX Neurology Panel 1 (human) Calibrator Blend	A00S0126	≤ -70°C	31 AUG 2027
S-PLEX Neurology Panel 1 (human) Coating Solution	527692	≤ -70°C	31 AUG 2028

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	391	fg/mL	80% - 120%	< 25%	142	fg/mL
NF-L	420 - 99800	4350	fg/mL	80% - 120%	< 25%	1440	fg/mL
Tau (total)	17.5 - 4160	38.8	fg/mL	80% - 120%	< 25%	21.1	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GFAP	20X Calibrator	15640000	fg/mL
NF-L	20X Calibrator	87000000	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		02 JUL 2025