



CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: S-PLEX[®] Neurology Panel 1 (human) Kit

Catalog Number: K15639S-series; K15640S-Series; K151AKGS-Series; K156AKGS-Series

Lot Number: K00S0122

Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human GFAP Antibody	D00T0123	2-8°C	31 MAY 2026
TURBO-BOOST [™] Hu Neurofilament L Antibody	D00T0124	2-8°C	31 MAY 2026
TURBO-BOOST [™] Human Tau Antibody	D00T0125	2-8°C	31 MAY 2026
S-PLEX Neurology Panel 1 (human) Calibrator Blend	A00S0073	2-8°C	30 SEP 2027
Diluent 63	M0630002	≤ -10°C	30 APR 2025
S-PLEX Neurology Panel 1 (human) Coating Solution	429071	≤ -70°C	31 MAY 2026
Blocker S1	R0010811	≤ -10°C	28 FEB 2025
Blocker S2	420803	≤ -10°C	30 SEP 2027
S-PLEX Enhance E1	Y0550010	≤ -10°C	31 MAY 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 MAR 2025
S-PLEX Enhance E3	Y0570012	≤ -70°C	31 MAY 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580015	≤ -70°C	30 APR 2025
S-PLEX Detect D2	Y0590011	≤ -70°C	31 MAR 2025

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
GFAP	53.0 - 12500	704	fg/mL	80 - 120%	< 25%	298	fg/mL
NF-L	193 - 45900	19800	fg/mL	80 - 120%	< 25%	2780	fg/mL
Tau (Total)	15.0 - 4160	1900	fg/mL	80 - 120%	< 25%	86.1	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GFAP	1X Calibrator	704000	fg/mL
NF-L	1X Calibrator	2470000	fg/mL
Tau (Total)	1X Calibrator	237000	fg/mL



CERTIFICATE OF ANALYSIS

Page 2 of 2

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	17 JAN 2023