

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Neurology Panel 1 (human) Kit
Lot Number: K00S0270
Expiration Date: 31 OCT 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	D0TB0042	2-8°C	31 JAN 2029
TURBO-BOOST Human GFAP Antibody	D0TB0041	2-8°C	30 NOV 2028
TURBO-BOOST Human Tau Antibody	D0TB0043	2-8°C	31 DEC 2028
Blocker S1	R0140017	≤ -10°C	29 FEB 2028
Blocker S2	562164	≤ -10°C	31 OCT 2030
S-PLEX Enhance E1	Y0550026	≤ -10°C	31 JAN 2028
S-PLEX Enhance E2	Y0560025	≤ -10°C	29 FEB 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580034	≤ -70°C	31 DEC 2027
S-PLEX Detect D2	Y0590027	≤ -70°C	31 JAN 2028
S-PLEX Enhance E3	Y0570028	≤ -70°C	31 JAN 2028
S-PLEX Neurology Panel 1 (human) Calibrator Blend	A00S0133	≤ -70°C	31 OCT 2027
S-PLEX Neurology Panel 1 (human) Coating Solution	528465	≤ -70°C	30 SEP 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	453	fg/mL	80% - 120%	< 25%	159	fg/mL
NF-L	420 - 99800	9140	fg/mL	80% - 120%	< 25%	1850	fg/mL
Tau (total)	17.5 - 4160	40.3	fg/mL	80% - 120%	< 25%	17.2	fg/mL

**CERTIFICATE OF ANALYSIS****Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
GFAP	20X Calibrator	18120000	fg/mL
NF-L	20X Calibrator	91400000	fg/mL
Tau (total)	20X Calibrator	3220000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 DEC 2025