



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human GFAP Kit
Catalog Number: K151AJXS-Series; K156AJXS-Series
Lot Number: K00S0111
Expiration Date: 30 Apr 2024

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human GFAP Detection Antibody	D00T0116	2-8°C	31 Mar 2026
Biotin Human GFAP Antibody	B00T0116	2-8°C	31 Mar 2026
Human GFAP Calibrator	A00S0064	2-8°C	31 Mar 2027
Diluent I2	M0120077	≤ -10°C	28 Feb 2025
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 Apr 2024
Blocker S1	R0010811	≤ -10°C	28 Feb 2025
S-PLEX Enhance E1	Y0550009	≤ -10°C	28 Feb 2025
S-PLEX Enhance E2	Y0560011	≤ -10°C	31 Mar 2025
S-PLEX Enhance E3	Y0570011	≤ -70°C	31 Mar 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:

Name	Calibrator	
	Recovery	CV
GFAP	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
GFAP	431 - 13800	3450	fg/mL	80% - 120%	< 20%	549	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
GFAP	1X Calibrator	4350000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 MAR 2023