



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human GFAP Kit
Catalog Number: K151AJXS-Series; K156AJXS-Series
Lot Number: K00S0080
Expiration Date: 31 AUG 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST™ Human GFAP Detection Antibody	D00T0083	2-8°C	30 SEP 2025
Biotin Human GFAP Antibody	B00T0083	2-8°C	30 SEP 2025
Human GFAP Calibrator	A00S0064	2-8°C	31 MAR 2027
Diluent I2	M0120073	≤ -10°C	31.AUG 2024
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 APR 2024
Blocker S1	R0140005	≤ -10°C	31 AUG 2023
S-PLEX Enhance E1	Y0550007	≤ -10°C	31 AUG 2024
S-PLEX Enhance E2	Y0560008	≤ -10°C	30 APR 2024
S-PLEX Enhance E3	Y0570008	≤ -70°C	30 Jun 2024
S-PLEX Detect D1	Y0580011	≤ -70°C	31 May 2024
S-PLEX Detect D2	Y0590008	≤ -70°C	31 Jul 2024

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	LLOQ				LLOD		
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
GFAP	431 - 13800	987	fg/mL	80% - 120%	< 20%	443	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		18 APR 2022