

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human Tau (pT181) Kit
Catalog Number: K151AGMS-Series; K156AGMS-Series
Lot Number: K00S0167
Expiration Date: 30 NOV 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT181) Antibody	B00T0176	2-8°C	31 JAN 2027
TURBO-BOOST™ Human Tau Antibody	D00T0176	2-8°C	31 MAR 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 2	M0020127	≤ -10°C	30 NOV 2025
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
Human Tau (pT181) Calibrator	A00S0097	≤ -70°C	31 MAY 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580023	≤ -70°C	31 JUL 2026
S-PLEX Detect D2	Y0590017	≤ -70°C	30 JUN 2026
S-PLEX Enhance E3	Y0570018	≤ -70°C	30 JUN 2026

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	Tau (pT181)	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (pT181)	115 - 1840	920	fg/mL	80% - 120%	< 20%	118	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT181)	20X Calibrator	53800000	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		16 APR 2024