

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human Tau (pT181) Kit
Catalog Number: K151AGMS-Series, K156AGMS-Series
Lot Number: K00S0109
Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT181) Antibody	B00T0114	2-8°C	31 MAR 2026
Human Tau (pT181) Calibrator	A00S0075	2-8°C	31 OCT 2027
TURBO-BOOST™ Human Tau Antibody	D00T0114	2-8°C	30 JUN 2026
Blocker S1	R0010824	≤ -10°C	30 JUN 2025
Diluent 2	M0020124	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 JUL 2025
S-PLEX Enhance E2	Y0560012	≤ -10°C	31 MAY 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580016	≤ -70°C	31 MAY 2025
S-PLEX Detect D2	Y0590012	≤ -70°C	30 APR 2025
S-PLEX Enhance E3	Y0570012	≤ -70°C	31 MAY 2025

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	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
Tau (pT181)	115 - 1840	460	fg/mL	80% - 120%	< 20%	185	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT181)	1X Calibrator	2210000	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		03 JAN 2024