



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX<sup>®</sup> Human Tau (pT217) Kit  
**Catalog Number:** K151APFS-series  
**Lot Number:** K00S0173  
**Expiration Date:** 28 FEB 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau Antibody	B00T0180	2-8°C	31 MAY 2027
TURBO-BOOST™ Human Tau Detection Antibody	D00T0180	2-8°C	30 APR 2027
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
Diluent 2	M0020124	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550015	≤ -10°C	31 MAY 2026
S-PLEX Enhance E2	Y0560016	≤ -10°C	31 MAY 2026
Human Tau (pT217) Calibrator	A00S0094	≤ -70°C	30 APR 2026
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580022	≤ -70°C	31 MAY 2026
S-PLEX Detect D2	Y0590016	≤ -70°C	30 APR 2026
S-PLEX Enhance E3	Y0570017	≤ -70°C	31 MAR 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 (pT217)	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
Tau (pT217)	1050 - 16800	2100	fg/mL	80% - 120%	< 20%	695	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
Tau (pT217)	20X Calibrator	72600000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		08 DEC 2023