

**CERTIFICATE OF ANALYSIS**

**Product Description:** S-PLEX Human Tau (pT231) Kit  
**Catalog Number:** K151AGNS-Series  
**Lot Number:** K00S0183  
**Expiration Date:** 28 FEB 2025

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT231) Antibody	B00T0115	2-8°C	31 AUG 2026
TURBO-BOOST™ Human Tau Antibody	D00T0114	2-8°C	30 JUN 2026
Blocker S1	R0010871	≤ -10°C	31 MAR 2026
Diluent 2	M0020124	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550014	≤ -10°C	28 FEB 2026
S-PLEX Enhance E2	Y0560016	≤ -10°C	31 MAY 2026
Human Tau (pT231) Calibrator	A00S0098	≤ -70°C	31 MAY 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 (pT231)	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (pT231)	3680 - 58800	7350	fg/mL	80% - 120%	< 20%	2880	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
Tau (pT231)	20X Calibrator	914000000	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		27 FEB 2024