



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] Human Tau (pT217) Kit
Catalog Number: K151APFS - Series
Lot Number: K00S0152
Expiration Date: 28 FEB 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT217) Antibody	B00T0156	2-8°C	31 JAN 2027
TURBO-BOOST [™] Human Tau Antibody	D00T0156	2-8°C	31 JAN 2027
Blocker S1	R0010832	≤ -10°C	31 OCT 2025
Diluent 2	M0020124	≤ -10°C	28 FEB 2025
S-PLEX Enhance E1	Y0550013	≤ -10°C	30 NOV 2025
S-PLEX Enhance E2	Y0560014	≤ -10°C	31 OCT 2025
Human Tau (pT217) Calibrator	A00S0088	≤ -70°C	31 JAN 2026
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 AUG 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580018	≤ -70°C	31 OCT 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 OCT 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 OCT 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 (pT217)	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (pT217)	1480 - 23600	5900	fg/mL	80% - 120%	< 20%	877	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT217)	20X Calibrator	71800000	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		19 JAN 2024