

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human Tau (pT217) Kit
Lot Number: K00S0232
Expiration Date: 31 MAR 2026
Catalog Number: K151APFS-Series; K156APFS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT217) Antibody	B00T0235	2-8°C	30 APR 2028
TURBO-BOOST Human Tau Antibody	D00T0235	2-8°C	31 MAR 2028
Blocker S1	R0140012	≤ -10°C	30 APR 2027
Diluent 2	M0020128	≤ -10°C	31 MAR 2026
S-PLEX Enhance E1	Y0550020	≤ -10°C	30 APR 2027
S-PLEX Enhance E2	Y0560020	≤ -10°C	31 MAR 2027
Human Tau (pT217) Calibrator	A00S0114	≤ -70°C	28 FEB 2027
S-PLEX Coating Reagent C1	446866	≤ -70°C	31 MAR 2026
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580029	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590022	≤ -70°C	30 APR 2027
S-PLEX Enhance E3	Y0570023	≤ -70°C	30 APR 2027

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)	1050 - 16800	4200	fg/mL	80% - 120%	< 20%	559	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT217)	20X Calibrator	70000000	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		02 JAN 2025