



Extension Memo

Page 1 of 1

Product Description: S-PLEX® Human Tau (pT217) Kit
Catalog Number(s): K151APFS-series
Lot Number: K00S0173
New Expiration Date: 28 Feb 2026

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 2	M0020124	28 Feb 2025	28 Feb 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		16 Dec 2024



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX[®] 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
Review/Approval	Karen Hamilla	Quality		08 DEC 2023