



Extension Memo

Product Description: S-PLEX Human Tau (pT231) Kit
Lot Number: K00S0241
New Expiration Date: 31 AUG 2026
Catalog Number(s): K151AGNS-Series, K156AGNS-Series

Based on additional testing, the original date of expiration for this kit has been extended by 4 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human Tau (pT231) Calibrator	A00S0098	31 MAY 2026	31 AUG 2026
S-PLEX Detect D2	Y0590016	30 APR 2026	31 AUG 2026
S-PLEX Enhance E1	Y0550015	31 MAY 2026	31 AUG 2026
S-PLEX Enhance E2	Y0560017	31 JUL 2026	31 AUG 2026
Blocker S1	R0140007	30 JUN 2026	31 AUG 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		26 Mar 2026

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX[®] Human Tau (pT231) Kit
Lot Number: K00S0241
Expiration Date: 30 APR 2026
Catalog Number: K151AGNS-Series; K156AGNS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT231) Antibody	B00T0241	2-8°C	30 APR 2028
TURBO-BOOST Human Tau Antibody	D00T0241	2-8°C	31 MAY 2028
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 2	M0020133	≤ -10°C	30 JUN 2027
S-PLEX Enhance E1	Y0550015	≤ -10°C	31 MAY 2026
S-PLEX Enhance E2	Y0560017	≤ -10°C	31 JUL 2026
Human Tau (pT231) Calibrator	A00S0098	≤ -70°C	31 MAY 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG [™] Reagent	Y0580030	≤ -70°C	30 APR 2027
S-PLEX Detect D2	Y0590016	≤ -70°C	30 APR 2026
S-PLEX Enhance E3	Y0570022	≤ -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 (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%	1340	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
Review/Approval	Karen Hamilla	Quality		25 MAR 2025