



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human Tau (pT231) Kit
Lot Number: K00S0196
Expiration Date: 31 MAR 2026
Catalog Number: K151AGNS-Series; K156AGNS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT231) Antibody	B00T0205	2-8°C	31 AUG 2027
Human Tau (pT231) Calibrator	A00S0099	2-8°C	30 NOV 2028
TURBO-BOOST Human Tau Antibody	D00T0205	2-8°C	31 AUG 2027
Blocker S1	R0140007	≤ -10°C	30 JUN 2026
Diluent 2	M0020128	≤ -10°C	31 MAR 2026
S-PLEX Enhance E1	Y0550016	≤ -10°C	31 JUL 2026
S-PLEX Enhance E2	Y0560018	≤ -10°C	31 AUG 2026
S-PLEX Coating Reagent C1	476750	≤ -70°C	31 MAR 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580025	≤ -70°C	30 SEP 2026
S-PLEX Detect D2	Y0590018	≤ -70°C	31 JUL 2026
S-PLEX Enhance E3	Y0570019	≤ -70°C	31 JUL 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%	1330	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT231)	1X Calibrator	39800000	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		10 JUN 2024