



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human Tau (pT231) Kit
Lot Number: K00S0278
Expiration Date: 31 MAY 2026
Catalog Number: K151AGNS-Series, K156AGNS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT231)Antibody	B00T0259	2-8°C	30 NOV 2028
TURBO-BOOST Human Tau Antibody	D00T0259	2-8°C	31 OCT 2028
Blocker S1	R0140015	≤ -10°C	31 OCT 2027
Diluent 2	M0020133	≤ -10°C	30 JUN 2027
S-PLEX Enhance E1	Y0550024	≤ -10°C	31 OCT 2027
S-PLEX Enhance E2	Y0560024	≤ -10°C	31 DEC 2027
Human Tau (pT231) Calibrator	A00S0098	≤ -70°C	31 MAY 2026
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580034	≤ -70°C	31 DEC 2027
S-PLEX Detect D2	Y0590026	≤ -70°C	31 DEC 2027
S-PLEX Enhance E3	Y0570027	≤ -70°C	31 DEC 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Human Tau (pT231) Calibrator	
		Recovery	CV	Recovery	CV
1	Tau (pT231)	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Tau (pT231)	14700	80% - 120%	< 20%	15271844	70% - 130%	< 30%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Tau (pT231)	914000000	fg/mL

Dilution to top of Curve: 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 OCT 2025