



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human Tau (pT231) Kit

Catalog Number: K151AGNS-Series; K156AGNS-Series

Lot Number: K00S0110

Expiration Date: 28 Feb 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human Tau Antibody	D00T0114	2-8°C	30 Jun 2026
Biotin Human Tau (pT231) Antibody	B00T0115	2-8°C	31 Aug 2026
Human Tau (pT231) Calibrator	A00S0052	2-8°C	30 Jun 2026
Diluent 2	M0020124	≤ -10°C	28 Feb 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 Aug 2025
Blocker S1	R0010824	≤ -10°C	30 Jun 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 Jul 2025
S-PLEX Enhance E2	Y0560013	≤ -10°C	31 Aug 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580016	≤ -70°C	31 May 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 Jul 2025

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

Precision, Accuracy, and Sensitivity Specifications:

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

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
Tau (pT231)	3680 - 58800	14700	fg/mL	80% - 120%	< 20%	2470	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT231)	1X Calibrator	40000000	fg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 MAY 2023