



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®] Human Tau (pT181) Kit
Catalog Number: K151AGMS-Series; K156AGMS - Series
Lot Number: K00S0168
Expiration Date: 28 Feb 2025

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau(pT181) Antibody	B00T0077	2-8°C	30 Apr 2025
TURBO-BOOST [™] Human Tau Antibody	D00T0077	2-8°C	30 Apr 2025
Human Tau (pT181) Calibrator	A00S0056	2-8°C	30 Sep 2026
Diluent 2	M0020124	≤ -10°C	28 Feb 2025
S-PLEX Coating Reagent C1	417723	≤ -70°C	31 Aug 2025
Blocker S1	R0010832	≤ -10°C	31 Oct 2025
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 Jul 2025
S-PLEX Enhance E2	Y0560012	≤ -10°C	31 May 2025
S-PLEX Enhance E3	Y0570014	≤ -70°C	31 Oct 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580017	≤ -70°C	31 Aug 2025
S-PLEX Detect D2	Y0590014	≤ -70°C	31 Oct 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(pT181)	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
Tau(pT181)	115 - 1840	230	fg/mL	80% - 120%	< 20%	75.1	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau(pT181)	1X Calibrator	2200000	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		19 SEP 2023