



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human Tau (pT181) Kit
Catalog Number: K151AGMS-Series; K156AGMS-Series
Lot Number: K00S0056
Expiration Date: 30 APR 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT181) Antibody	B00T0066	2-8°C	30 NOV 2024
Human Tau (pT181) Calibrator	A00S0050	2-8°C	31 MAY 2026
TURBO-BOOST™ Human Tau Antibody	D00T0066	2-8°C	31 DEC 2024
Blocker S1	R0140004	≤ -10°C	30 APR 2023
Diluent 2	M0020115	≤ -10°C	31 MAY 2023
S-PLEX Enhance E1	Y0550006	≤ -10°C	31 OCT 2023
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 OCT 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 AUG 2023
S-PLEX Detect D1 with TURBO-TAG™ reagent	Y0580008	≤ -70°C	30 NOV 2023
S-PLEX Detect D2	Y0590006	≤ -70°C	31 OCT 2023
S-PLEX Enhance E3	Y0570006	≤ -70°C	31 OCT 2023

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 (pT181)	80% - 120%	< 20%

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (pT181)	115 - 1840	460	fg/mL	80% - 120%	< 20%	76.7	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT181)	1X Calibrator	2110000	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		03 JAN 2024