



## CERTIFICATE OF ANALYSIS

**Product Description:** S-PLEX Human Tau (pT181) Kit  
**Catalog Number:** K151AGMS-Series; K156AGMS-Series  
**Lot Number:** K00S0068  
**Expiration Date:** 31 MAY 2023

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT181) Antibody	B00T0077	2-8°C	30 APR 2025
Human Tau (pT181 ) Calibrator	A00S0056	2-8°C	30 SEP 2026
TURBO-BOOST™ Human Tau Antibody	D00T0077	2-8°C	30 APR 2025
Blocker S1	R0140005	≤ -10°C	31 AUG 2023
Diluent 2	M0020115	≤ -10°C	31 MAY 2023
S-PLEX Enhance E1	R0140006	≤ -10°C	31 JAN 2024
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	Y0590007	≤ -70°C	29 FEB 2024
S-PLEX Enhance E3	Y0570007	≤ -70°C	29 FEB 2024

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	920	fg/mL	80% - 120%	< 20%	114	fg/mL

**Stock Concentration of Calibrator:**

Name	Item	Concentration	Units
Tau (pT181)	1X Calibrator	2200000	fg/mL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		04 JAN 2024