

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human Tau (total) Kit
Catalog Number: K151AGPS-Series, K156AGPS-Series
Lot Number: K00S0067
Expiration Date: 31 MAY 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (total) Antibody	B00T0076	2-8°C	31 MAR 2025
Human Tau (total) Calibrator	A00S0055	2-8°C	30 SEP 2026
TURBO-BOOST™ Human Tau Antibody	D00T0076	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 (total)	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
Tau (total)	21.7 - 347	86.8	fg/mL	80% - 120%	< 20%	18.6	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (total)	3X Calibrator	657000	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		29 DEC 2023