



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human Tau (total) Kit
Catalog Number: K151AGPS-Series; K156AGPS-Series
Lot Number: K00S0058
Expiration Date: 30 APR 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human Tau Antibody	D00T0068	2-8°C	31 Dec 2024
Biotin Human Tau (total) Antibody	B00T0068	2-8°C	30 Nov 2024
Human Tau (total) Calibrator	A00S0051	2-8°C	30 Jun 2026
Diluent 2	M0020115	≤ -10°C	31 May 2023
S-PLEX Coating Reagent C1	329998	≤ -70°C	31 Aug 2023
Blocker S1	R0140004	≤ -10°C	30 Apr 2023
S-PLEX Enhance E1	Y0550006	≤ -10°C	31 Oct 2023
S-PLEX Enhance E2	Y0560007	≤ -10°C	31 Oct 2023
S-PLEX Enhance E3	Y0570006	≤ -70°C	31 Oct 2023
S-PLEX Detect D1	Y0580008	≤ -70°C	30 Nov 2023
S-PLEX Detect D2	Y0590006	≤ -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:

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

Name	LLOQ					LLOD	
	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (total)	21.7 - 347	43.4	fg/mL	80% - 120%	< 20%	11.6	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (total)	3X Calibrator	712000	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	<i>Karen Hamilla</i>	09 FEB 2022