

**CERTIFICATE OF ANALYSIS**

Product Description: S-PLEX Human BD-Tau (total) Kit
Lot Number: K00S0282
Expiration Date: 31 JUL 2027
Catalog Number: K151AXFS-Series, K156AXFS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human BD-Tau Antibody	B00T0262	2-8°C	30 NOV 2028
TURBO-BOOST Hu BD-Tau (total) Antibody	D00T0262	2-8°C	30 NOV 2028
Blocker S1	R0140013	≤ -10°C	31 JUL 2027
S-PLEX Enhance E1	Y0550024	≤ -10°C	31 OCT 2027
S-PLEX Enhance E2	Y0560023	≤ -10°C	31 OCT 2027
Human BD-Tau (total) Calibrator	A00S0137	≤ -70°C	31 OCT 2027
S-PLEX Coating Reagent C1	521329	≤ -70°C	31 DEC 2027
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580033	≤ -70°C	30 NOV 2027
S-PLEX Detect D2	Y0590025	≤ -70°C	30 NOV 2027
S-PLEX Enhance E3	Y0570026	≤ -70°C	31 OCT 2027

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	BD-Tau (total)	80% - 120%	< 20%

Name	Range Tested	Observed	LLOQ			LLOD	
			Units	Recovery	CV	Observed	Units
BD-Tau (total)	118 - 3760	941	fg/mL	80% - 120%	< 20%	47.2	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
BD-Tau (total)	20X Calibrator	18280000	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		17 JUL 2025