



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human BD-Tau (pT181) Kit
Lot Number: K00S0284
Expiration Date: 31 OCT 2027
Catalog Number: K151AXHS-Series, K156AXHS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human BD-Tau Antibody	B00T0267	2-8°C	30 NOV 2028
TURBO-BOOST Hu BD-Tau (pT181) Antibody	D00T0264	2-8°C	30 NOV 2028
S-PLEX Enhance E1	Y0550024	≤ -10°C	31 OCT 2027
S-PLEX Enhance E2	Y0560023	≤ -10°C	31 OCT 2027
Blocker S1	R0140015	≤ -70°C	31 OCT 2027
Human BD-Tau (pT181) Calibrator	A00S0139	≤ -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:

		Calibrator	
Spot Number	Name	Recovery	CV
1	BD-Tau (pT181)	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
BD-Tau (pT181)	363 - 11600	2900	fg/mL	80% - 120%	< 20%	209	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
BD-Tau (pT181)	20X Calibrator	36000000	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		23 JUL 2025