



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human Tau (pT217) Kit
Lot Number: K00S0318
Expiration Date: 30 APR 2027
Catalog Number: K151APFS-Series; K156APFS-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT217) Antibody	B00T0289	2-8°C	31 AUG 2029
TURBO-BOOST™ Human Tau Antibody	D00T0289	2-8°C	30 SEP 2029
Blocker S1	R0140019	≤ -10°C	31 MAY 2028
Diluent 2	M0020139	≤ -10°C	30 NOV 2028
S-PLEX Enhance E1	Y0550027	≤ -10°C	29 FEB 2028
S-PLEX Enhance E2	Y0560027	≤ -10°C	31 MAR 2028
Human Tau (pT217) Calibrator	A00S0121	≤ -70°C	30 APR 2027
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580036	≤ -70°C	30 APR 2028
S-PLEX Detect D2	Y0590029	≤ -70°C	31 MAY 2028
S-PLEX Enhance E3	Y0570030	≤ -70°C	31 MAR 2028

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	Tau (pT217)	80% - 120%	< 20%

LLOQ						LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (pT217)	1050 - 16800	3670	fg/mL	80% - 120%	< 20%	735	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT217)	20X Calibrator	74600000	fg/mL

This lot meets MSD's non-specific binding specifications.

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		04 AUG 2026