

**CERTIFICATE OF ANALYSIS**

Product Description: Human BD-Tau (pT217) Calibrator
Catalog Number: C01AXF-2
Lot Number: A00S0153
Expiration Date: 30 NOV 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
BD-Tau (pT217)	72200000	fg/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 JUN 2026