



**CERTIFICATE OF ANALYSIS**

**Product Description:** Human Tau (pT217) Calibrator  
**Catalog Number:** C01APF-2  
**Lot Number:** A00S0121  
**Expiration Date:** 30 APR 2027  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
Tau (pT217)	74600000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		26 DEC 2024