

**CERTIFICATE OF ANALYSIS**


**Product Description:** Human BD-Tau (pT181) Calibrator  
**Catalog Number:** C01AXG-2  
**Lot Number:** A00S0139  
**Expiration Date:** 31 OCT 2027  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
BD-Tau (pT181)	36000000	fg/mL

**Dilution to Top of Curve:** 20X

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		30 JUN 2025