

**CERTIFICATE OF ANALYSIS**


Product Description: Human BD-Tau (total) Calibrator
Catalog Number: C01AXE-2
Lot Number: A00S0137
Expiration Date: 31 OCT 2027
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

| Analyte | Stock Concentration | Unit of Measure |
|----------------|---------------------|-----------------|
| BD-Tau (total) | 18280000 | 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 |  | 06 JUN 2025 |