



CERTIFICATE OF ANALYSIS

Product Description: Human Tau (pT181) Calibrator
Catalog Number: C01ARH-2
Lot Number: A00S0115
Expiration Date: 30 APR 2027
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Tau (pT181)	39000000	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	<i>Karen Hamilla</i>	20 JAN 2025