



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Tau Calibrator

Catalog Number: C01LA-2

Lot Number: A0080278

Expiration Date: 31 JUL 2023


Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Tau	224000	pg/mL

Dilution to Top of Curve: 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		29 APR 2021