



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Tau Calibrator
Catalog Number: C01LA-2
Lot Number: A0080120
Expiration Date: 31 JUL 2019
Storage Temperature: ≤ -70°C

Concentration of Materials Tested:

	Assay
	Tau
Calibrator (pg/mL)	412000

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	30 MAR 2017