



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Tau Calibrator  
**Catalog Number:** C01LA-2  
**Lot Number:** A0080179  
**Expiration Date:** 28 FEB 2021  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration of Materials Tested:

	Assay
	Tau
Calibrator (pg/mL)	250000

Concentration: 20X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	08 OCT 2018