



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human α -synuclein Calibrator

Catalog Number: C01WK-2

Lot Number: A0080115

Expiration Date: 30 JUN 2019

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

Concentration of Materials Tested:

	Assay
	hu α -synuclein
Calibrator (pg/mL)	210 000

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>	27 JAN 2017