



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human α -Synuclein Calibrator

Catalog Number: C01WK-2

Lot Number: A0080128

Expiration Date: 30 SEP 2019

Storage Temperature: -70°C

Concentration of Materials Tested:

	Assay
	hu α-Synuclein
Calibrator (pg/mL)	198 200

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>	21 APR 2017