



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human α -Synuclein Calibrator

Catalog Number: C01WK-2

Lot Number: A0080237

Expiration Date: 30 APR 2022

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
α -Synuclein	210000	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		14 JAN 2020