



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human α -Synuclein Calibrator

Catalog Number: C01WK-2

Lot Number: A0080272

Expiration Date: 31 MAY 2023


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

Concentration Assignment:

| Analyte | Stock Concentration | Unit of Measure |
|---------------------|---------------------|-----------------|
| α -Synuclein | 202000 | 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 |  | 22 DEC 2020 |