



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human α -Synuclein Calibrator
Catalog Number: C01WK-2
Lot Number: A00U0270
Expiration Date: 30 NOV 2024
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

| Analyte | Stock Concentration | Unit of Measure |
|---------------------|---------------------|-----------------|
| α -Synuclein | 200000 | 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 | | 16 JUN 2022 |