



CERTIFICATE OF ANALYSIS

Product Description: Human α -Synuclein Calibrator
Catalog Number: C01WK-2
Lot Number: A0080362
Expiration Date: 31 AUG 2025
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
α -Synuclein	214000	pg/mL

Dilution to Top of Curve: 20X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 APR 2023