



CERTIFICATE OF ANALYSIS


Product Description: Human α -Synuclein Calibrator
Catalog Number: C01WK-2
Lot Number: A0080441
Expiration Date: 29 FEB 2028
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: 20 X

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		03 OCT 2025