



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Human α -Synuclein Control Pack

Catalog Number: C41WK-1

Lot Number: CP000137

Expiration Date: 31 JAN 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
α -Synuclein	16600	1550	182	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human α -Synuclein Control 1	A00C0857	$\leq -70^{\circ}\text{C}$	31 JAN 2025
Human α -Synuclein Control 2	A00C0858	$\leq -70^{\circ}\text{C}$	31 JAN 2025
Human α -Synuclein Control 3	A00C0859	$\leq -70^{\circ}\text{C}$	31 JAN 2025

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		25 OCT 2022