



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Human α -Synuclein Control Pack

Catalog Number: C41WK-1

Lot Number: CP000094

Expiration Date: 28 FEB 2022

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
α -Synuclein	17 000	1 610	174	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human α -Synuclein Control 1	A00C0670	$\leq -70^{\circ}\text{C}$	28 FEB 2022
Human α -Synuclein Control 2	A00C0671	$\leq -70^{\circ}\text{C}$	28 FEB 2022
Human α -Synuclein Control 3	A00C0672	$\leq -70^{\circ}\text{C}$	28 FEB 2022

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		11 SEP 2019