



RFE MEMO

Page 1 of 1

Product Description(s): Human α -Synuclein Control Pack

Catalog Number(s) / Size(s): C41WK-1

Lot Number: CP000057

Based on additional testing, the original date of expiration for this Control Pack has been extended by 41 months. The expiration date of CP000057 has been extended to 30 APR 2023

The expiration dates for the following controls have been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human α -Synuclein Control 1	A00C0501	31 OCT 2019	30 APR 2023
Human α -Synuclein Control 2	A00C0502	31 OCT 2019	30 APR 2023
Human α -Synuclein Control 3	A00C0503	31 OCT 2019	30 APR 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		12 MAY 2022

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human α -Synuclein Controls

Catalog Number: C41WK-1

Lot Number: CP000057

Expiration Date: 31 OCT 2019

Concentration of Materials Tested:

Units = pg/mL	Assay
	Human α -Synuclein
Control 1 (A00C0501)	17923.23
Control 2 (A00C0502)	1690.28
Control 3 (A00C0503)	178.47

Control concentrations were assigned across a minimum of three kit lots.

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human α -Synuclein Control 1	A00C0501	$\leq -70^{\circ}\text{C}$	31 OCT 2019
Human α -Synuclein Control 2	A00C0502	$\leq -70^{\circ}\text{C}$	31 OCT 2019
Human α -Synuclein Control 3	A00C0503	$\leq -70^{\circ}\text{C}$	31 OCT 2019

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	20JUL2017