



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Human  $\alpha$ -Synuclein Control pack

**Catalog Number:** C41WK-1

**Lot Number:** CP000106

**Expiration Date:** 31 DEC 2022

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
$\alpha$ -Synuclein	18400	1740	200	pg/mL

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human $\alpha$ -Synuclein Control 1	A00C0723	$\leq -70^{\circ}\text{C}$	31 DEC 2022
Human $\alpha$ -Synuclein Control 2	A00C0724	$\leq -70^{\circ}\text{C}$	31 DEC 2022
Human $\alpha$ -Synuclein Control 3	A00C0725	$\leq -70^{\circ}\text{C}$	31 DEC 2022

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		19 AUG 2020