



CERTIFICATE OF ANALYSIS

**Control Pack Description:** Human  $\alpha$ -Synuclein Control Pack  
**Catalog Number:** C41WK-1  
**Lot Number:** CP000167  
**Expiration Date:** 31 JAN 2027

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human $\alpha$ -Synuclein Control 1	A00C0946	$\leq -70^{\circ}\text{C}$	31 JAN 2027
Human $\alpha$ -Synuclein Control 2	A00C0947	$\leq -70^{\circ}\text{C}$	31 JAN 2027
Human $\alpha$ -Synuclein Control 3	A00C0948	$\leq -70^{\circ}\text{C}$	31 JAN 2027

**Concentration Assignment:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
$\alpha$ -Synuclein	18400	1920	218	pg/mL

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 SEP 2024