



CERTIFICATE OF ANALYSIS

Page 1 of 1

Control Pack Description: Neurodegeneration Control Pack 1 (6E10)

Catalog Number: C41LB-1

Lot Number: CP000134

Expiration Date: 30 JUN 2025

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
A β 40	9960	3320	563	pg/mL
A β 38	4190	1030	209	pg/mL
A β 42	1430	497	201	pg/mL

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Neurodegeneration Control 1 (6E10)	A00C0828	$\leq -70^{\circ}\text{C}$	30 JUN 2025
Neurodegeneration Control 2 (6E10)	A00C0829	$\leq -70^{\circ}\text{C}$	30 JUN 2025
Neurodegeneration Control 3 (6E10)	A00C0830	$\leq -70^{\circ}\text{C}$	30 JUN 2025

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 AUG 2022