



Extension Memo

Page 1 of 1

Product Description: Neurodegeneration Control Pack 1 (6E10)
Catalog Number(s): C41LB-1
Lot Number: CP000134
New Expiration Date: 30 Jun 2026

Based on additional testing, the original date of expiration for this control pack has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Neurodegeneration Control 1 (6E10)	A00C0828	30 Jun 2025	30 Jun 2026
Neurodegeneration Control 2 (6E10)	A00C0829	30 Jun 2025	30 Jun 2026
Neurodegeneration Control 3 (6E10)	A00C0830	30 Jun 2025	30 Jun 2026

Function	Name	Signature	Date
Quality	Michelle Likiam		25 Apr 2025



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