



## Extension Memo

Page 1 of 1

**Product Description:** Neurodegeneration Control Pack 1 (4G8)  
**Catalog Number(s):** C40RQ-1  
**Lot Number:** CP000089  
**New Expiration Date:** 31 Mar 2026

Based on additional testing, the original date of expiration for this control pack has been extended by 48 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Neurodegeneration Control 1 (4G8)	A00C0643	31 Mar 2022	31 Mar 2026
Neurodegeneration Control 2 (4G8)	A00C0644	31 Mar 2022	31 Mar 2026
Neurodegeneration Control 3 (4G8)	A00C0645	31 Mar 2022	31 Mar 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		12 Feb 2025



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Control Pack Description:** Neurodegeneration Control Pack 1 (4G8)

**Catalog Number:** C40RQ-1

**Lot Number:** CP000089

**Expiration Date:** 31 MAR 2022

**Concentration Assignment:** The following concentrations were measured using the A $\beta$  Peptide Panel 1 (4G8) V-PLEX® kit:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
A $\beta$ 40	7790	2650	417	pg/mL
A $\beta$ 38	3840	961	223	pg/mL
A $\beta$ 42	1490	516	205	pg/mL


**The following concentrations were measured using the Human Total Tau V-PLEX® kit:**

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Tau	1330	415	161	pg/mL

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Neurodegeneration Control 1 (4G8)	A00C0643	$\leq -70^{\circ}\text{C}$	31 MAR 2022
Neurodegeneration Control 2 (4G8)	A00C0644	$\leq -70^{\circ}\text{C}$	31 MAR 2022
Neurodegeneration Control 3 (4G8)	A00C0645	$\leq -70^{\circ}\text{C}$	31 MAR 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		05 MAR 2024