



RFE MEMO

Page 1 of 1

Product Description(s): Neurodegeneration Control Pack

Catalog Number(s) / Size(s): C41LB-1

Lot Number: CP000070

Based on additional testing, the original date of expiration for this control pack has been extended by 23 months. The expiration date of CP000070 has been extended to 31 DEC 2022.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Neurodegeneration Control 1	A00C0567	31 JAN 2021	31 DEC 2022
Neurodegeneration Control 2	A00C0568	31 JAN 2021	31 DEC 2022
Neurodegeneration Control 3	A00C0569	31 JAN 2021	31 DEC 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		24 JAN 2022

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Neurodegeneration Control Pack

Catalog Number: C41LB-1

Lot Number: CP000070

Expiration Date: 31 JAN 2021

Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration of Materials Tested:

The following concentrations were measured using the A β Peptide Panel 1 (4G8) V-PLEX[®] kit:

Units = pg/mL	4G8 Assays		
	A β 40	A β 38	A β 42
Control 1 (A00C0567)	9060	4330	1550
Control 2 (A00C0568)	3160	1120	558
Control 3 (A00C0569)	521	277	231

The following concentrations were measured using A β Peptide Panel 1 (6E10) V-PLEX[®] kit and Human A β 42 Kit V-PLEX[®] kit:

Units = pg/mL	6E10 Assays		
	A β 40	A β 38	A β 42
Control 1 (A00C0567)	9430	4160	1440
Control 2 (A00C0568)	3190	1050	506
Control 3 (A00C0569)	534	231	203

The following concentrations were measured using the Human Total Tau V-PLEX[®] kit

Units = pg/mL	Assay
	Tau
Control 1 (A00C0567)	1370
Control 2 (A00C0568)	436
Control 3 (A00C0569)	169

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Neurodegeneration Control 1	A00C0567	$\leq -70^{\circ}\text{C}$	31 JAN 2021
Neurodegeneration Control 2	A00C0568	$\leq -70^{\circ}\text{C}$	31 JAN 2021
Neurodegeneration Control 3	A00C0569	$\leq -70^{\circ}\text{C}$	31 JAN 2021

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality		20 MAY 2019