



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus A β 42 Peptide (4G8) Kit


Catalog Number(s) / Size(s): K150SL-Series; K151LB-Series

Lot Number: K0081874

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081874 has been extended to 31 MAR 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
A β 1-42 Peptide (4G8)	A0080294	31 Oct 2023	31 MAR 2024
Diluent 35	M0350032	31 MAR 2023	31 MAR 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		12 JUL 2023

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

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus A β 42 Peptide (4G8) Kit
Kit Lot Number: K0081874
Expiration Date: 31 Mar 2023
Kit Catalog Numbers: K150SL-Series; K151LB-Series

Precision and Accuracy Specifications:

		Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A β 42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	A β 42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	A β 42	74800	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
AB42 Peptide Plate	Z0057054	2-8°C	31 Jul 2024
Anti-A β 4G8 Antibody	D0081743	2-8°C	30 Apr 2025
A β 1-42 Peptide (4G8)	A0080294	\leq -70°C	31 Oct 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 MAY 2022