



Extension Memo

Page 1 of 1

Product Description: V-PLEX and V-PLEX Plus A β 42 Peptide (4G8) Kit
Catalog Number(s): K150SL-Series
Lot Number: K0082476
New Expiration Date: 28 Feb 2026

Based on additional testing, the original date of expiration for this kit 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
A β 42 Peptide Plate	Z0057246	28 Feb 2025	28 Feb 2026
Diluent 35	M0350036	31 Jul 2025	28 Feb 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		18 Dec 2024

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX and V-PLEX Plus A β 42 Peptide (4G8) Kit
Lot Number: K0082476
Expiration Date: 28 FEB 2025
Catalog Number: K150SL-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A β 4G8 Antibody	D0081899	2-8°C	28 FEB 2026
A β 42 Peptide Plate	Z0057246	2-8°C	28 FEB 2025
Diluent 35	M0350036	2-8°C	31 JUL 2025
A β 1-42 Peptide (4G8)	A0080393	\leq -70°C	30 SEP 2026

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

Precision, Accuracy, and Sensitivity Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	A β 42	2.5	80% - 120%	< 20%	1270	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 42	80400	pg/mL

Dilution to top of Curve: 40X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 JUN 2024