



### Extension Memo

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8)Kit  
**Catalog Number(s):** K150SJ-Series  
**Lot Number:** K0082585  
**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
Aβ40 Peptide 96-Well 4-Spot SECTOR Plate	Z0057252	28 FEB 2025	28 Feb 2026
Aβ1-40 Peptide (4G8)	A0080359	31 JUL 2025	28 Feb 2026
Anti-Aβ 4G8 Antibody	D0081873	31 DEC 2025	28 Feb 2026
Diluent 35	M0350037	31 DEC 2025	28 Feb 2026

Function	Name	Signature	Date
Quality	Michelle Likiam		02 Dec 2024



### CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8) Kit  
**Lot Number:** K0082585  
**Expiration Date:** 28 FEB 2025  
**Catalog Number:** K150SJ-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-Aβ 4G8 Antibody	D0081873	2-8°C	31 DEC 2025
Aβ40 Peptide Plate	Z0057252	2-8°C	28 FEB 2025
Aβ40 Blocker	R0010896	2-8°C	30 SEP 2026
Diluent 100	Y0270334	2-8°C	30 SEP 2026
Diluent 35	M0350037	2-8°C	31 DEC 2025
Aβ1-40 Peptide (4G8)	A0080359	≤ -70°C	31 JUL 2025

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	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	20	80% - 120%	< 20%	6000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	Aβ40	616000	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
<b>Review/Approval</b>	Michelle Likiam	Quality		27 NOV 2024