



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** V-PLEX and V-PLEX Plus A $\beta$ 40 Peptide (4G8) Kit


**Catalog Number(s) / Size(s):** K150SJ-Series

**Lot Number:** K0082125

Based on additional testing, the original date of expiration for this Kit has been extended by 12 months. The expiration date of K0082125 has been extended to 31 OCT 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
A $\beta$ 1-40 Peptide (4G8)	A0080293	31 OCT 2023	31 OCT 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		05 OCT 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 $\beta$ 40 Peptide (4G8) Kit  
**Kit Lot Number:** K0082125  
**Expiration Date:** 31 Oct 2023  
**Kit Catalog Numbers:** K150SJ-Series

**Precision and Accuracy 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 $\beta$ 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	A $\beta$ 40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	A $\beta$ 40	652000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ 40 Peptide Plate	Z0057252	2-8°C	28 Feb 2025
Anti-A $\beta$ 4G8 Antibody	D0081873	2-8°C	31 Dec 2025
A $\beta$ 1-40 Peptide (4G8)	A0080293	$\leq$ -70°C	31 Oct 2023
A $\beta$ 40 Blocker	R0010797	2-8°C	31 Oct 2024
Diluent 100	Y0270313	2-8°C	30 Nov 2024
Diluent 35	M0350033	2-8°C	31 Oct 2024

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	Xiomara Talavera	Quality		01 NOV 2022