



## EXTENSION MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX PLUS A $\beta$  Peptide Panel 1 (6E10) Kit


**Catalog Number(s) / Size(s):** K15200-Series, K150SI-Series

**Lot Number:** K0082181

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082181 has been extended to 30 Sep 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
A $\beta$ 1-38 Peptide (6E10)	A0080287	30 Sep 2023	30 Sep 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		15 JUN 2023

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

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX PLUS A $\beta$  Peptide Panel 1 (6E10) Kit  
**Kit Lot Number:** K0082181  
**Expiration Date:** 30 Sep 2023  
**Kit Catalog Numbers:** K15200-Series, K150SI-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Neurodegeneration Control 1 (6E10)		Neurodegeneration Control 2 (6E10)		Neurodegeneration Control 3 (6E10)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A $\beta$ 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A $\beta$ 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A $\beta$ 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 $\beta$ 40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	A $\beta$ 38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	A $\beta$ 42	3.13	75% - 125%	< 20%	1270	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

Spot	Name	NSB Spec
1	A $\beta$ 40	< 1.0%
2	A $\beta$ 38	< 1.0%
4	A $\beta$ 42	< 1.0%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	A $\beta$ 40	624000	pg/mL
2	A $\beta$ 38	488000	pg/mL
4	A $\beta$ 42	72400	pg/mL


**CERTIFICATE OF ANALYSIS**

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 (6E10) Plate	Z0057294	2-8°C	30 Apr 2025
Anti-Aβ 6E10 Antibody	D0081958	2-8°C	31 Jul 2026
AB1-42 Peptide (6E10)	A0080350	≤ -70°C	31 May 2025
Aβ1-38 Peptide (6E10)	A0080287	≤ -70°C	30 Sep 2023
Aβ1-40 Peptide (6E10)	A0080349	≤ -70°C	31 May 2025
Aβ40 Blocker	R0010817	2-8°C	31 Mar 2025
Diluent 100	Y0270316	2-8°C	31 Jan 2025
Diluent 35	M0350034	2-8°C	31 Jan 2025

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		24 FEB 2023