



## Extension Memo

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human A $\beta$ 42 Kit  
**Catalog Number(s):** K151LB - Series  
**Lot Number:** K0082421  
**New Expiration Date:** 31 Jan 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 $\beta$ 42 Peptide Plate	Z0057235	31 Jan 2025	31 Jan 2026
A $\beta$ 1-42 Peptide (6E10)	A0080350	31 May 2025	31 Jan 2026
Diluent 35	M0350036	31 Jul 2025	31 Jan 2026

Function	Name	Signature	Date
Quality	Karen Hamilla		20 Nov 2024

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human Aβ42 Kit  
**Kit Lot Number:** K0082421  
**Expiration Date:** 31 Jan 2025  
**Kit Catalog Numbers:** K151LB - Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
AB42 Peptide Plate	Z0057235	2-8°C	31 Jan 2025
Anti-Aβ 6E10 Antibody	D0081958	2-8°C	31 Jul 2026
AB1-42 Peptide (6E10)	A0080350	≤ -70°C	31 May 2025
Diluent 35	M0350036	2-8°C	31 Jul 2025

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

**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	AB42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%


Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	AB42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	AB42	72400	pg/mL

Dilution to Top of Curve = 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 Mar 2024