



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX PLUS A β 40 Peptide (6E10) Kit


Catalog Number(s) / Size(s): K150SK-Series

Lot Number: K0082187

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082187 has been extended to 30 SEP 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
A β 1-40 Peptide (6E10)	A0080277	30 SEP 2023	30 SEP 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		08 NOV 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 Aβ 40 Peptide (6E10) Kit
Kit Lot Number: K0082187
Expiration Date: 30 Sep 2023
Kit Catalog Numbers: K150SK-Series

Precision and Accuracy Specifications:

		Calibrator		Neurodegeneration Control 1 (6E10)		Neurodegeneration Control 2 (6E10)		Neurodegeneration Control 3 (6E10)	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	Aβ 40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Aβ 40	612000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0057297	2-8°C	31 May 2025
Anti-Aβ 6E10 Antibody	D0081958	2-8°C	31 Jul 2026
Aβ1-40 Peptide (6E10)	A0080277	≤ -70°C	30 Sep 2023
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		17 MAR 2023