



## RFE MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus A $\beta$ 40 Peptide (6E10) Kit

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

**Lot Number:** K0081733

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081733 been extended to 31 JAN 2023.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
A $\beta$ 40 Peptide Plate	Z0056593	31 MAR 2022	31 JAN 2023
Anti-A $\beta$ 6E10 Antibody	D0081119	31 JAN 2022	31 JAN 2023
A $\beta$ 1-40 Peptide (6E10)	A0080203	31 MAY 2022	31 JAN 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		22 FEB 2022

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 Aβ40 Peptide (6E10) Kit  
**Kit Lot Number:** K0081733  
**Expiration Date:** 31 Jan 2022  
**Kit Catalog Numbers:** K150SK-Series

**Precision and Accuracy Specifications:**

		Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
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	600000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056593	2-8°C	31 Mar 2022
Anti-Aβ 6E10 Antibody	D0081119	2-8°C	31 Jan 2022
Aβ1-40 Peptide (6E10)	A0080203	≤ -70°C	31 May 2022
Aβ40 Blocker	R0010728	2-8°C	31 Aug 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023
Diluent 100	Y0270290	2-8°C	31 Jul 2023

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		10 MAY 2021