



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: K0082115

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082115 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
Anti-A β 6E10 Antibody	D0081630	31 JUL 2024	30 SEP 2024
A β 1-40 Peptide (6E10)	A0080277	30 SEP 2023	30 SEP 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		10 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 A β 40 Peptide (6E10) Kit
Kit Lot Number: K0082115
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	A β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	A β 40	612000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β 40 PeptideMULTI-SPOT® 96-Well	Z0057242	2-8°C	28 Feb 2025
Anti-A β 6E10 Antibody	D0081630	2-8°C	31 Jul 2024
A β 1-40 Peptide (6E10)	A0080277	\leq -70°C	30 Sep 2023
A β 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	Karen Hamilla	Quality		18 OCT 2022