



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus A β 40 Peptide (4G8) Kit


Catalog Number(s) / Size(s): K150SJ-Series

Lot Number: K0082050

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082050 has been extended to 31 OCT 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 (4G8)	A0080293	31 OCT 2023	31 OCT 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		09 OCT 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 Aβ40 Peptide (4G8) Kit
Kit Lot Number: K0082050
Expiration Date: 31 Oct 2023
Kit Catalog Numbers: K150SJ-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ1-40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ1-40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Aβ1-40	652000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0057187	2-8°C	30 Nov 2024
Anti-Aβ 4G8 Antibody	D0081795	2-8°C	30 Sep 2025
Aβ1-40 Peptide (4G8)	A0080293	≤ -70°C	31 Oct 2023
Diluent 35	M0350033	2-8°C	31 Oct 2024
Diluent 100	Y0270311	2-8°C	31 Oct 2024
Aβ40 Blocker	R0010797	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		09 AUG 2022