



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX and V-PLEX Plus A β 42 Peptide (4G8) Kit


Catalog Number(s) / Size(s): K150SL-Series

Lot Number: K0082117

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082117 has been extended to 31 MAR 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
A β 1-42 Peptide (4G8)	A0080294	31 OCT 2023	31 MAR 2024
Diluent 35	M0350032	31 MAR 2023	31 MAR 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		15 SEP 2023

This memo supersedes all other statements regarding date of expiration for this product.



Meso Scale Diagnostics, LLC.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX and V-PLEX Plus Aβ42 Peptide (4G8) Kit
Kit Lot Number: K0082117
Expiration Date: 31 Mar 2023
Kit Catalog Numbers: K150SL-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
AB42 Peptide Plate	Z0057246	2-8°C	28 Feb 2025
Anti-Aβ 4G8 Antibody	D0081899	2-8°C	28 Feb 2026
Aβ1-42 Peptide (4G8)	A0080294	≤ -70°C	31 Oct 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023

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 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Aβ42	74800	pg/mL

Dilution to Top of Curve = 40X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 SEP 2023