



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-Plex Plus Aβ42 Peptide (4G8) Kit


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

**Lot Number:** K0081410

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081410 has been extended to 31 May 2022.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Aβ1-42 Peptide (4G8)	A0080201	31 Jul 2021	31 May 2022
Diluent 35	M0350030	31 May 2021	31 May 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		25 JAN 2021

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

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-Plex Plus Aβ42 Peptide (4G8) Kit

**Kit Catalog Numbers:** K150SL-Series

**Kit Lot Number:** K0081410

**Expiration Date:** 31 May 2021

**Acceptance Data/ Percent Recovery 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 Material Tested :**

Spot	Name	Item	Conc	Units
1	Aβ42	40X Calibrator	99200	0
		Neurodegeneration Control 1 (4G8)	1410	0
		Neurodegeneration Control 2 (4G8)	515	0
		Neurodegeneration Control 3 (4G8)	211	0

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

**Standard Curve Data:**

Aβ42 - Spot 1		
Conc. 0	ECL Signal	%CV
2480	1,250,000	1.1
620	395,000	5.4
155	66,100	3.2
38.8	7,910	5.0
9.69	1,010	4.3
2.42	293	2.1
0.605	196	3.7
0.00	168	3.2

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ42 Peptide Plate	Z0056655	2-8°C	31 Aug 2022
Anti-Aβ 4G8 Antibody	D0081432	2-8°C	31 Aug 2023
Aβ1-42 Peptide (4G8)	A0080201	≤ -70°C	31 Jul 2021
Diluent 35	M0350030	2-8°C	31 May 2021
Diluent 100	Not Kit Specific	Not Kit Specific	N/A
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20x), 100mL	Not Kit Specific
Neurodegeneration Control 1 (4G8)	Not Kit Specific
Neurodegeneration Control 2 (4G8)	Not Kit Specific
Neurodegeneration Control 3 (4G8)	Not Kit Specific

### Statement:

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		14 APR 2020