



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Aβ 42 Peptide (4G8) Kit

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

**Lot Number:** K0040347

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040347 has been extended to 31 Jan 2019.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Aβ42 Peptide Plate	Z0055569	30 Nov 2018	31 JAN 2019
Anti-Aβ 4G8 Antibody	D0040613	30 Sep 2018	31 JAN 2019
Aβ1- 42 Peptide	A0040134	31 Jan 2018	31 JAN 2019
Diluent 35	M0350024	30 Sep 2018	31 JAN 2019
Diluent 100	Y0270232	31 May 2018	31 JAN 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	12 JAN 2018

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

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

COA-07147
-----------

**Product Description:** V-PLEX® and V-PLEX Plus Aβ 42 Peptide (4G8) Kit  
**Kit Catalog Numbers:** K150SLE-1, K150SLE-2, K150SLE-4, K150SLG-1, K150SLG-2, K150SLG-4  
**Kit Lot Number:** K0040347  
**Expiration Date:** 31 Jan 2018

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
		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	55200	pg/mL
		Neurodegeneration Control 1	1680	pg/mL
		Neurodegeneration Control 2	545	pg/mL
		Neurodegeneration Control 3	230	pg/mL

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

**CERTIFICATE OF ANALYSIS**
**Standard Curve Data:**

A $\beta$ 42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1380	1,160,000	1.6
345	287,000	1.7
86.3	54,800	2.2
21.6	8,900	1.3
5.39	1,650	2.5
1.35	500	2.5
0.337	269	3.1
0.00	194	3.3


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ 42 Peptide Plate	Z0055569	2-8°C	30 Nov 2018
Anti-A $\beta$ 4G8 Antibody	D0040613	2-8°C	30 Sep 2018
A $\beta$ 1- 42 Peptide	A0040134	$\leq$ -70°C	31 Jan 2018
Diluent 35	M0350024	2-8°C	30 Sep 2018
Diluent 100	Y0270232	2-8°C	31 May 2018
MSD Read Buffer T (4X)	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	Not Kit Specific
Neurodegeneration Control 2	Not Kit Specific
Neurodegeneration Control 3	Not Kit Specific

**Statement:**

The above kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		02 AUG 2016