

## 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:** K0081259  
**Expiration Date:** 31 Mar 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	pg/mL
		Neurodegeneration Control 1 (4G8)	1410	pg/mL
		Neurodegeneration Control 2 (4G8)	515	pg/mL
		Neurodegeneration Control 3 (4G8)	211	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

### Standard Curve Data:

Aβ42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2480	1,120,000	1.9
620	333,000	3.4
155	53,300	2.5
38.8	6,550	4.6
9.69	930	1.6
2.42	234	3.7
0.605	115	6.2
0.00	78	8.2

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ42 Peptide Plate	Z0056387	2-8°C	31 Mar 2021
Anti-Aβ 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
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		19 JUL 2019