

## CERTIFICATE OF ANALYSIS

Page 1 of 2

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

**Kit Catalog Numbers:** K150SL-serise

**Kit Lot Number:** K0081110

**Expiration Date:** 30 Nov 2019

### 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	Aβ1-42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ1-42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ1-42	40X Calibrator	94400	pg/mL
		Neurodegeneration Control 1	1530	pg/mL
		Neurodegeneration Control 2	555	pg/mL
		Neurodegeneration Control 3	237	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:

Aβ1-42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2360	1,140,000	1.0
590	324,000	3.1
148	47,800	2.7
36.9	5,220	3.1
9.22	653	3.1
2.30	168	3.5
0.00	98	4.6

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ 42 Peptide Plate	Z0056387	2-8°C	31 Mar 2021
Anti-A $\beta$ 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
A $\beta$ 1-42 Peptide	A0080140	$\leq$ -70°C	30 Nov 2019
Diluent 35	M0350029	2-8°C	30 Sep 2020
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	Not Kit Specific
Neurodegeneration Control 2	Not Kit Specific
Neurodegeneration Control 3	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>	Dawn Stollar	Quality Assurance		07 NOV 2018