

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Aβ Peptide Panel 1 (6E10) Kit

**Kit Catalog Numbers:** K15200, K15199-Series

**Kit Lot Number:** K0080979

**Expiration Date:** 31 Oct 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	Aβ-40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Aβ-38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	Aβ-42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ-40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	Aβ-38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	Aβ-42	3.13	80% - 120%	< 20%	1270	80% - 120%	< 20%

**Acceptance Data/ Non Specific Binding Specifications:**

Spot	Name	NSB Spec
1	Aβ-40	< 1.0%
2	Aβ-38	< 1.0%
4	Aβ-42	< 1.0%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	Aβ-40	40X Calibrator	608000	pg/mL
		Neurodegeneration Control 1	10100	pg/mL
		Neurodegeneration Control 2	3290	pg/mL
		Neurodegeneration Control 3	571	pg/mL
2	Aβ-38	40X Calibrator	564000	pg/mL
		Neurodegeneration Control 1	4160	pg/mL
		Neurodegeneration Control 2	1060	pg/mL
		Neurodegeneration Control 3	236	pg/mL
4	Aβ-42	40X Calibrator	94400	pg/mL
		Neurodegeneration Control 1	1360	pg/mL
		Neurodegeneration Control 2	476	pg/mL
		Neurodegeneration Control 3	193	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

Page 2 of 2

### Standard Curve Data:

A $\beta$ -40 - Spot 1			A $\beta$ -38 - Spot 2			A $\beta$ -42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
15200	825,000	4.7	14100	814,000	3.9	2360	983,000	2.2
3800	289,000	3.9	3530	158,000	4.1	590	276,000	2.5
950	55,700	4.6	881	16,800	5.4	148	48,800	4.2
238	6,630	4.7	220	1,920	4.9	36.9	6,010	6.4
59.4	777	4.2	55.1	451	5.1	9.22	964	5.9
14.8	203	3.9	13.8	264	5.9	2.30	260	4.8
3.71	112	5.9	3.44	221	6.2	0.576	130	3.7
0.00	92	11.0	0.00	218	8.9	0.00	93	7.8

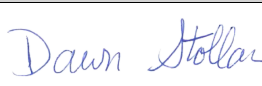
### Components:

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ Peptide Panel 1 Plate	Z0056210	2-8°C	31 Aug 2020
Anti-A $\beta$ 6E10 Antibody	D0081076	2-8°C	31 Oct 2021
A $\beta$ 1-38 Peptide	A0080166	≤ -70°C	30 Sep 2020
A $\beta$ 1-40 Peptide	A0080143	≤ -70°C	31 Dec 2019
A $\beta$ 1-42 Peptide	A0080140	≤ -70°C	30 Nov 2019
A $\beta$ 40 Blocker	R0110016	2-8°C	30 Sep 2020
Diluent 35	M0350028	2-8°C	31 Dec 2019
Diluent 100	Y0270249	2-8°C	31 Oct 2019
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), 100 mL	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
Review/Approval	Dawn Stollar	Quality Assurance		23 MAY 2018