

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus A $\beta$  Peptide Panel 1 (6E10) Kit

**Kit Catalog Numbers:** K15200-Series, K150SI-Series

**Kit Lot Number:** K0081483

**Expiration Date:** 31 Mar 2022

**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	A $\beta$ 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A $\beta$ 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A $\beta$ 42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

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

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

Spot	Name	NSB Spec
1	A $\beta$ 40	< 1.0%
2	A $\beta$ 38	< 1.0%
4	A $\beta$ 42	< 1.0%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	A $\beta$ 40	40X Calibrator	600000	pg/mL
		Neurodegeneration Control 1	9430	pg/mL
		Neurodegeneration Control 2	3190	pg/mL
		Neurodegeneration Control 3	534	pg/mL
2	A $\beta$ 38	40X Calibrator	508000	pg/mL
		Neurodegeneration Control 1	4160	pg/mL
		Neurodegeneration Control 2	1050	pg/mL
		Neurodegeneration Control 3	231	pg/mL
4	A $\beta$ 42	40X Calibrator	73600	pg/mL
		Neurodegeneration Control 1	1440	pg/mL
		Neurodegeneration Control 2	506	pg/mL
		Neurodegeneration Control 3	203	pg/mL

Controls tested may not be the same lot as shipped.

**CERTIFICATE OF ANALYSIS**

Page 2 of 2

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:**

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
15000	715,000	3.8	12700	948,000	3.3	1840	829,000	2.5
3750	241,000	7.1	3180	195,000	5.8	460	228,000	4.1
938	40,500	8.7	794	22,200	5.6	115	37,900	4.4
234	4,470	7.1	198	2,540	5.9	28.8	4,970	5.0
58.6	633	5.1	49.6	513	4.6	7.19	934	2.7
14.6	188	2.5	12.4	216	5.0	1.80	266	4.3
3.66	115	6.3	3.10	155	8.6	0.449	128	4.8
0.00	94	7.2	0.00	144	9.2	0.00	83	8.8

**Components:**


Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ Peptide Panel 1 (6E10) Plate	Z0056699	2-8°C	31 Oct 2022
Anti- A $\beta$ 6E10 Antibody	D0081509	2-8°C	30 Sep 2023
A $\beta$ 1-38 Peptide (6E10)	A0080243	≤ -70°C	31 Jul 2022
A $\beta$ 1-40 Peptide (6E10)	A0080203	≤ -70°C	31 May 2022
A $\beta$ 1-42 Peptide (6E10)	A0080245	≤ -70°C	31 Jul 2022
A $\beta$ 40 Blocker	R0010692	2-8°C	30 Sep 2022
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270273	2-8°C	30 Sep 2022
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
Review/Approval	Karen Hamilla	Quality		18 JUN 2020