

CERTIFICATE OF ANALYSIS

Page 1 of 2

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

Kit Catalog Numbers: K15200-Series; K150SI-Series

Kit Lot Number: K0081375

Expiration Date: 30 Sep 2021

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 β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A β 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A β 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 β 40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	A β 38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	A β 42	3.13	80% - 120%	< 20%	1270	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	A β 40	< 1.0%
2	A β 38	< 1.0%
4	A β 42	< 1.0%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	A β 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 β 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 β 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.

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 β 40 - Spot 1			A β 38 - Spot 2			A β 42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
15000	752,000	5.0	12700	905,000	4.2	1840	862,000	2.5
3750	250,000	5.9	3180	174,000	6.0	460	240,000	3.5
938	45,700	8.9	794	18,200	7.6	115	42,000	6.6
234	4,810	10.0	198	1,850	8.0	28.8	5,810	6.2
58.6	543	10.0	49.6	332	6.3	7.19	1,010	5.3
14.6	139	5.6	12.4	141	5.6	1.80	270	3.6
3.66	87	4.7	3.10	106	6.4	0.449	127	4.1
0.00	73	6.7	0.00	92	6.1	0.00	82	7.3


Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 (6E10) Plate	Z0056627	2-8°C	31 May 2022
Anti-A β 6E10 Antibody	D0081402	2-8°C	31 May 2023
A β 1-38 Peptide (6E10)	A0080243	≤ -70°C	31 Jul 2022
A β 1-40 Peptide (6E10)	A0080203	≤ -70°C	31 May 2022
A β 1-42 Peptide (6E10)	A0080245	≤ -70°C	31 Jul 2022
A β 40 Blocker	R0010657	2-8°C	30 Sep 2021
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270270	2-8°C	30 Jun 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		14 APR 2020