

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
Kit Lot Number: K0080920
Expiration Date: 30 Sep 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	A β 1-40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A β 1-38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A β 1-42	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 1-40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	A β 1-38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	A β 1-42	3.13	80% - 120%	< 20%	1270	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	A β 1-40	40X Calibrator	608000	pg/mL
		Neurodegeneration Control 1	10200	pg/mL
		Neurodegeneration Control 2	3720	pg/mL
		Neurodegeneration Control 3	555	pg/mL
2	A β 1-38	40X Calibrator	564000	pg/mL
		Neurodegeneration Control 1	4410	pg/mL
		Neurodegeneration Control 2	1130	pg/mL
		Neurodegeneration Control 3	227	pg/mL
4	A β 1-42	40X Calibrator	91200	pg/mL
		Neurodegeneration Control 1	1470	pg/mL
		Neurodegeneration Control 2	494	pg/mL
		Neurodegeneration Control 3	207	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

Standard Curve Data:

A β 1-40 - Spot 1			A β 1-38 - Spot 2			A β 1-42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
15200	781,000	2.4	14100	806,000	3.1	2280	940,000	1.7
3800	241,000	2.9	3530	134,000	4.0	570	246,000	2.8
950	40,100	5.1	881	13,500	4.3	143	39,700	3.4
238	4,310	6.0	220	1,530	4.1	35.6	5,140	4.8
59.4	532	5.1	55.1	398	4.3	8.91	920	3.3
14.8	156	3.3	13.8	228	6.9	2.23	251	3.7
3.71	97	5.0	3.44	196	7.5	0.557	125	3.8
0.00	81	6.6	0.00	189	7.3	0.00	85	6.5


Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 Plate	Z0056109	2-8°C	30 Apr 2020
Anti-A β 6E10 Antibody	D0080985	2-8°C	31 Oct 2019
A β 1-38 Peptide	A0080134	≤ -70°C	30 Nov 2019
A β 1-40 Peptide	A0080129	≤ -70°C	30 Sep 2019
A β 1-42 Peptide	A0080130	≤ -70°C	30 Sep 2019
A β 40 Blocker	R0110015	2-8°C	31 Dec 2019
Diluent 35	M0350028	2-8°C	31 Dec 2019
Diluent 100	Y0270252	2-8°C	29 Feb 2020
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), 1000mL	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		21 DEC 2017