



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-Series

Kit Lot Number: K0080759

Expiration Date: 30 Jun 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	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Aβ38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	Aβ42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	Aβ38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	Aβ42	3.13	80% - 120%	< 20%	1270	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	Aβ40	< 1.0%
2	Aβ38	< 1.0%
4	Aβ42	< 1.0%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ40	40X Calibrator	604000	pg/mL
		Neurodegeneration Control 1	10200	pg/mL
		Neurodegeneration Control 2	3720	pg/mL
		Neurodegeneration Control 3	555	pg/mL
2	Aβ38	40X Calibrator	584000	pg/mL
		Neurodegeneration Control 1	4410	pg/mL
		Neurodegeneration Control 2	1130	pg/mL
		Neurodegeneration Control 3	227	pg/mL
4	Aβ42	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

Page 2 of 2

Standard Curve Data:

Aβ40 - Spot 1			Aβ38 - Spot 2			Aβ42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
15100	290,000	5.3	14600	971,000	2.9	2280	861,000	2.4
3780	84,200	5.3	3650	194,000	3.4	570	240,000	1.8
944	14,200	7.2	913	20,800	3.1	143	40,000	2.8
236	1,670	8.9	228	2,150	5.3	35.6	5,250	3.4
59.0	306	6.5	57.0	412	4.8	8.91	873	4.4
14.7	131	5.5	14.3	192	5.1	2.23	244	4.6
3.69	99	5.0	3.56	147	5.5	0.557	118	5.0
0.00	92	5.9	0.00	135	4.8	0.00	81	4.5


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 Plate	Z0055876	2-8°C	31 Aug 2019
Anti-Aβ 6E10 Antibody	D0080970	2-8°C	31 Aug 2019
Aβ1-38 Peptide	A0080117	≤ -70°C	30 Jun 2019
Aβ1-40 Peptide	A0080118	≤ -70°C	30 Jun 2019
Aβ1-42 Peptide	A0080130	≤ -70°C	30 Sep 2019
Aβ40 Blocker	R0110013	2-8°C	30 Jun 2019
Diluent 35	M0350026	2-8°C	31 Jul 2019
Diluent 100	Y0270248	2-8°C	31 Aug 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), 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	Dawn Stollar	Quality Assurance		04 MAY 2017