



CERTIFICATE OF ANALYSIS

COA-5885

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus Aβ Peptide Panel 1 (4G8) Kit
Kit Catalog Numbers: K15199E-1, K15199E-2, K15199E-4, K15199G-1, K15199G-2, K15199G-4
Kit Lot Number: K0040309
Expiration Date: 31 Oct 2016

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	ND Matrix Control 1		ND Matrix Control 2		ND Matrix Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ40	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Aβ38	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	Aβ42	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	Aβ38	60.0	80% - 120%	< 20%	7500	80% - 120%	< 20%
4	Aβ42	2.50	80% - 120%	< 20%	1271	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSBSpec
1	Aβ40	NA
1	Aβ40	< 1.0%
2	Aβ38	NA
2	Aβ38	< 1.0%
4	Aβ42	NA
4	Aβ42	< 1.0%



CERTIFICATE OF ANALYSIS

COA-5885

Page 2 of 4

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	A β 40	40X Calibrator	620000	pg/mL
		Neurodegeneration Control 1	9690	pg/mL
		Neurodegeneration Control 2	3480	pg/mL
		Neurodegeneration Control 3	534	pg/mL
		ND Matrix Control 1	7747	pg/mL
		ND Matrix Control 2	4054	pg/mL
		ND Matrix Control 3	2073	pg/mL
2	A β 38	40X Calibrator	431000	pg/mL
		Neurodegeneration Control 1	3550	pg/mL
		Neurodegeneration Control 2	820	pg/mL
		Neurodegeneration Control 3	160	pg/mL
		ND Matrix Control 1	4160	pg/mL
		ND Matrix Control 2	4964	pg/mL
		ND Matrix Control 3	455	pg/mL
4	A β 42	40X Calibrator	53300	pg/mL
		Neurodegeneration Control 1	1680	pg/mL
		Neurodegeneration Control 2	545	pg/mL
		Neurodegeneration Control 3	230	pg/mL
		ND Matrix Control 1	1854	pg/mL
		ND Matrix Control 2	615	pg/mL
		ND Matrix Control 3	296	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

COA-5885

Standard Curve Data:

Aβ40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
15500	512,233	4.1
3875	170,773	7.4
969	37,359	6.6
242	5,103	7.7
60.5	815	4.7
15.1	261	4.8
3.78	165	3.0
0	141	3.4

Aβ38 - Spot 2		
Conc. pg/mL	ECL Signal	%CV
10775	633,316	2.4
2694	100,791	3.8
673	9,032	2.5
168	890	2.5
42.1	240	3.1
10.5	150	4.6
2.63	133	4.4
0	136	5.5

Aβ42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1333	819,801	4.4
333	201,758	3.5
83.3	30,222	3.9
20.8	3,516	5.3
5.21	631	3.9
1.30	200	4.2
0.325	130	5.0
0	109	3.9

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 Plate	Z0055385	2-8°C	31 Dec 2017
Anti-Aβ 4G8 Antibody	D0040541	2-8°C	31 Dec 2017
Aβ1-38 Peptide	A0040099	≤ -70°C	31 Oct 2016
Aβ1-40 Peptide	A0040118	≤ -70°C	31 Oct 2017
Aβ1-42 Peptide	A0040101	≤ -70°C	31 Oct 2016
Diluent 35	M0350021	2-8°C	30 Sep 2017
Diluent 100	Y0270221	2-8°C	30 Nov 2017
Aβ40 Blocker	R0110009	2-8°C	30 Nov 2017
MSD Read Buffer T (4X)	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



CERTIFICATE OF ANALYSIS

COA-5885

Page 4 of 4

Statement:

The above kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Monika Jain	Quality Assurance	<i>Monika Jain</i>	19 AUG 2015