

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus A β Peptide Panel 1 (4G8) Kit; V-PLEX® and V-PLEX Plus A β 38 Peptide (4G8) Kit
Kit Catalog Numbers: K15199-Series, K150SH-Series
Kit Lot Number: K0080978
Expiration Date: 30 Nov 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	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	A β 1-38	60.0	80% - 120%	< 20%	7500	80% - 120%	< 20%
4	A β 1-42	2.50	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	9540	pg/mL
		Neurodegeneration Control 2	3270	pg/mL
		Neurodegeneration Control 3	547	pg/mL
2	A β 1-38	40X Calibrator	592000	pg/mL
		Neurodegeneration Control 1	4120	pg/mL
		Neurodegeneration Control 2	1110	pg/mL
		Neurodegeneration Control 3	281	pg/mL
4	A β 1-42	40X Calibrator	94400	pg/mL
		Neurodegeneration Control 1	1530	pg/mL
		Neurodegeneration Control 2	555	pg/mL
		Neurodegeneration Control 3	237	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:

Aβ1-40 - Spot 1			Aβ1-38 - Spot 2			Aβ1-42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
15200	777,000	6.5	14800	506,000	2.9	2360	1,250,000	1.5
3800	290,000	4.9	3700	87,400	4.3	590	377,000	1.6
950	68,000	6.9	925	8,550	5.5	148	66,400	3.9
238	9,690	4.2	231	975	4.0	36.9	8,280	3.4
59.4	1,300	4.5	57.8	233	4.8	9.22	1,210	5.3
14.8	317	3.3	14.5	127	5.9	2.30	279	4.3
3.71	147	3.1	3.61	102	7.8	0.576	128	5.9
0.00	104	7.4	0.00	98	10.0	0.00	83	8.9


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 Plate	Z0056207	2-8°C	31 Aug 2020
Anti-Aβ 4G8 Antibody	D0081010	2-8°C	31 Dec 2019
Aβ1-38 Peptide	A0080148	≤ -70°C	31 Aug 2020
Aβ1-40 Peptide	A0080143	≤ -70°C	31 Dec 2019
Aβ1-42 Peptide	A0080140	≤ -70°C	30 Nov 2019
Aβ40 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		01 MAY 2018