

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, K150SI-Series

Kit Lot Number: K0081053

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	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	608000	pg/mL
		Neurodegeneration Control 1	10100	pg/mL
		Neurodegeneration Control 2	3290	pg/mL
		Neurodegeneration Control 3	571	pg/mL
2	Aβ38	40X Calibrator	564000	pg/mL
		Neurodegeneration Control 1	4160	pg/mL
		Neurodegeneration Control 2	1060	pg/mL
		Neurodegeneration Control 3	236	pg/mL
4	Aβ42	40X Calibrator	94400	pg/mL
		Neurodegeneration Control 1	1360	pg/mL
		Neurodegeneration Control 2	476	pg/mL
		Neurodegeneration Control 3	193	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:

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
15200	796777	3.1	14100	777188	4.3	2360	959369	2.4
3800	281348	3.4	3530	147560	4.4	590	270935	3.4
950	50867	4.8	881	15179	5.3	148	46057	4.4
238	5467	7.4	220	1576	6.2	36.9	5597	6.2
59.4	697	5.1	55.1	344	6.8	9.22	1001	4.7
14.8	183	7.6	13.8	158	8.5	2.30	279	7.7
3.71	106	13.6	3.44	118	9.1	0.576	132	11.7
0.00	81	9.4	0.00	108	8.1	0.00	87	5.6


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 Plate	Z0056313	2-8°C	31 Dec 2020
Anti-Aβ 6E10 Antibody	D0081119	2-8°C	31 Jan 2022
Aβ1-38 Peptide	A0080166	≤ -70°C	30 Sep 2020
Aβ1-40 Peptide	A0080143	≤ -70°C	31 Dec 2019
Aβ1-42 Peptide	A0080140	≤ -70°C	30 Nov 2019
Aβ40 Blocker	R0010642	2-8°C	31 Jan 2021
Diluent 35	M0350029	2-8°C	30 Sep 2020
Diluent 100	Y0270255	2-8°C	30 Sep 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), 100 mL	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		19 DEC 2018