

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, K15199-Series

Kit Lot Number: K0081344

Expiration Date: 30 Sep 2020

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	Aβ1-40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Aβ1-38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	Aβ1-42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ1-40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	Aβ1-38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	Aβ1-42	3.13	80% - 120%	< 20%	1270	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	Aβ1-40	< 1.0%
2	Aβ1-38	< 1.0%
4	Aβ1-42	< 1.0%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ1-40	40X Calibrator	604000	pg/mL
		Neurodegeneration Control 1	9430	pg/mL
		Neurodegeneration Control 2	3190	pg/mL
		Neurodegeneration Control 3	534	pg/mL
2	Aβ1-38	40X Calibrator	516000	pg/mL
		Neurodegeneration Control 1	4160	pg/mL
		Neurodegeneration Control 2	1050	pg/mL
		Neurodegeneration Control 3	231	pg/mL
4	Aβ1-42	40X Calibrator	70000	pg/mL
		Neurodegeneration Control 1	1440	pg/mL
		Neurodegeneration Control 2	506	pg/mL
		Neurodegeneration Control 3	203	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:

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
15100	569,000	4.4	12900	597,000	5.5	1750	583,000	4.3
3780	185,000	4.8	3230	101,000	4.6	438	141,000	5.4
944	32,200	10.0	806	10,500	7.3	109	21,600	8.2
236	3,840	7.5	202	1,260	7.4	27.3	2,980	8.6
59.0	578	7.2	50.4	311	6.2	6.84	598	7.4
14.7	168	4.7	12.6	148	5.2	1.71	181	4.9
3.69	99	6.9	3.15	124	9.0	0.427	100	5.5
0.00	77	6.9	0.00	111	9.5	0.00	71	7.0


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	A0080181	≤ -70°C	28 Feb 2021
Aβ1-40 Peptide	A0080190	≤ -70°C	30 Apr 2021
Aβ1-42 Peptide (6E10)	A0080200	≤ -70°C	30 Jun 2021
Aβ40 Blocker	R0010642	2-8°C	31 Jan 2021
Diluent 35	M0350029	2-8°C	30 Sep 2020
Diluent 100	Y0270260	2-8°C	30 Jun 2021
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	Karen Hamilla	Quality		26 NOV 2019