

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus A β Peptide Panel 1 (4G8) Kit
Kit Catalog Numbers: K15199-Series, K150SH-Series
Kit Lot Number: K0081147
Expiration Date: 31 May 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A β -40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A β -38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A β -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 -40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	A β -38	60.0	80% - 120%	< 20%	7500	80% - 120%	< 20%
4	A β -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 -40	40X Calibrator	584000	pg/mL
		Neurodegeneration Control 1 (4G8)	8040	pg/mL
		Neurodegeneration Control 2 (4G8)	2700	pg/mL
		Neurodegeneration Control 3 (4G8)	445	pg/mL
2	A β -38	40X Calibrator	480000	pg/mL
		Neurodegeneration Control 1 (4G8)	4020	pg/mL
		Neurodegeneration Control 2 (4G8)	1050	pg/mL
		Neurodegeneration Control 3 (4G8)	256	pg/mL
4	A β -42	40X Calibrator	99200	pg/mL
		Neurodegeneration Control 1 (4G8)	1410	pg/mL
		Neurodegeneration Control 2 (4G8)	515	pg/mL
		Neurodegeneration Control 3 (4G8)	211	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:

A β -40 - Spot 1			A β -38 - Spot 2			A β -42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
14600	795,000	2.9	12000	347,000	2.9	2480	1,120,000	1.8
3650	285,000	3.0	3000	56,200	3.6	620	322,000	2.2
913	62,900	6.7	750	5,800	5.2	155	53,100	4.0
228	8,680	6.2	188	789	3.1	38.8	6,320	4.6
57.0	1,200	4.1	46.9	220	4.9	9.69	948	2.8
14.3	291	3.6	11.7	122	5.5	2.42	226	4.1
3.56	142	4.0	2.93	101	6.2	0.605	114	4.7
0.00	100	4.6	0.00	97	7.7	0.00	81	6.9


Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 Plate	Z0056396	2-8°C	31 May 2021
Anti-A β 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
A β 1-38 Peptide (4G8)	A0080189	≤ -70°C	30 Sep 2021
A β 1-40 Peptide (4G8)	A0080202	≤ -70°C	30 Sep 2021
A β 1-42 Peptide (4G8)	A0080201	≤ -70°C	31 Jul 2021
A β 40 Blocker	R0010657	2-8°C	30 Sep 2021
Diluent 35	M0350030	2-8°C	31 May 2021
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 (4G8)	Not Kit Specific
Neurodegeneration Control 2 (4G8)	Not Kit Specific
Neurodegeneration Control 3 (4G8)	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		18 JUL 2019