

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: K0081219
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

Page 2 of 2

Standard Curve Data:

A β 1-40 - Spot 1			A β 1-38 - Spot 2			A β 1-42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
15200	716,000	3.1	14800	381,000	4.1	2360	1,070,000	2.2
3800	238,000	3.9	3700	66,800	6.0	590	304,000	2.7
950	50,900	5.5	925	6,700	4.1	148	50,400	4.1
238	6,730	4.5	231	793	2.4	36.9	5,900	4.3
59.4	1,050	3.4	57.8	208	3.0	9.22	876	3.2
14.8	264	4.5	14.5	115	3.3	2.30	221	3.6
3.71	131	3.2	3.61	94	5.2	0.576	117	4.1
0.00	95	4.2	0.00	90	7.7	0.00	83	7.1

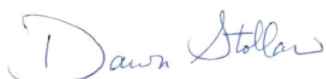
Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 Plate	Z0056284	2-8°C	30 Nov 2020
Anti-A β 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
A β 1-38 Peptide	A0080148	≤ -70°C	31 Aug 2020
A β 1-40 Peptide	A0080143	≤ -70°C	31 Dec 2019
A β 1-42 Peptide	A0080140	≤ -70°C	30 Nov 2019
A β 40 Blocker	R0110016	2-8°C	30 Sep 2020
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), 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		28 MAR 2019