

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Aβ Peptide Panel 1 (4G8) Kit

**Kit Catalog Numbers:** K15199-Series, K150SH-Series

**Kit Lot Number:** K0081314

**Expiration Date:** 31 Jul 2021

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ 40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	Aβ 38	60.0	80% - 120%	< 20%	7500	80% - 120%	< 20%
4	Aβ 42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

**Acceptance Data/ Non Specific Binding Specifications:**

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

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	Aβ 40	40X Calibrator	584000	pg/mL
		Neurodegeneration Control 1	8040	pg/mL
		Neurodegeneration Control 2	2700	pg/mL
		Neurodegeneration Control 3	445	pg/mL
2	Aβ 38	40X Calibrator	480000	pg/mL
		Neurodegeneration Control 1	4020	pg/mL
		Neurodegeneration Control 2	1050	pg/mL
		Neurodegeneration Control 3	256	pg/mL
4	Aβ 42	40X Calibrator	99200	pg/mL
		Neurodegeneration Control 1	1410	pg/mL
		Neurodegeneration Control 2	515	pg/mL
		Neurodegeneration Control 3	211	pg/mL

Controls tested may not be the same lot as shipped.

## CERTIFICATE OF ANALYSIS

Page 2 of 2

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

### Standard Curve Data:

Aβ 40 - Spot 1			Aβ 38 - Spot 2			Aβ 42 - Spot 4		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
14600	736,000	3.3	12000	288,000	3.2	2480	1,060,000	1.9
3650	262,000	5.3	3000	49,400	5.3	620	323,000	4.2
913	58,200	7.8	750	5,020	7.6	155	53,600	7.0
228	7,640	7.3	188	670	3.7	38.8	6,300	5.9
57.0	890	9.9	46.9	173	6.7	9.69	837	5.6
14.3	190	6.0	11.7	99	6.4	2.42	207	5.5
3.56	97	3.9	2.93	82	7.0	0.605	104	5.7
0.00	79	12.0	0.00	82	15.0	0.00	80	15.0


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 (4G8) Plate	Z0056580	2-8°C	28 Feb 2022
Anti-Aβ 4G8 Antibody	D0081341	2-8°C	30 Jun 2023
Aβ1-38 Peptide (4G8)	A0080189	≤ -70°C	30 Sep 2021
Aβ1-40 Peptide (4G8)	A0080202	≤ -70°C	30 Sep 2021
Aβ1-42 Peptide (4G8)	A0080201	≤ -70°C	31 Jul 2021
Aβ 40 Blocker	R0010657	2-8°C	30 Sep 2021
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270266	2-8°C	28 Feb 2022
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	Karen Hamilla	Quality		06 MAR 2020