

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8) Kit  
**Kit Catalog Numbers:** K150SJ-Series  
**Kit Lot Number:** K0081361  
**Expiration Date:** 30 Sep 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	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ40	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

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

### Standard Curve Data:

Aβ40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
14600	790,000	4.5
3650	302,000	3.9
913	67,200	2.8
228	8,620	6.3
57.0	1,110	5.4
14.3	216	5.4
3.56	102	6.2
0.00	75	3.9

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ 40 Peptide Plate	Z0056621	2-8°C	30 Apr 2022
Anti-A $\beta$ 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
A $\beta$ 1-40 Peptide (4G8)	A0080202	≤ -70°C	30 Sep 2021
A $\beta$ 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 (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
<b>Review/Approval</b>	Xiomara Talavera	Quality		12 DEC 2019