

**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:** K0080970

**Expiration Date:** 30 Jun 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	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	604000	pg/mL
		Neurodegeneration Control 1	9540	pg/mL
		Neurodegeneration Control 2	3270	pg/mL
		Neurodegeneration Control 3	547	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
15100	899,000	2.4
3780	314,000	4.1
944	66,100	3.3
236	8,380	5.0
59.0	1,190	4.9
14.7	331	2.3
3.69	203	2.3
0.00	163	3.4

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056180	2-8°C	31 Aug 2020
Anti-Aβ 4G8 Antibody	D0080871	2-8°C	30 Jun 2019
Aβ1-40 Peptide	A0080118	≤ -70°C	30 Jun 2019
Aβ40 Blocker	R0110015	2-8°C	31 Dec 2019
Diluent 35	M0350028	2-8°C	31 Dec 2019
Diluent 100	Y0270252	2-8°C	29 Feb 2020
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
<b>Review/Approval</b>	Jessica Drolet	Quality Assurance		18 May 2018