

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Aβ40 Peptide (6E10) Kit

**Kit Catalog Numbers:** K150SK-series

**Kit Lot Number:** K0081137

**Expiration Date:** 30 Sep 2020

**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%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ-40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	Aβ-40	40X Calibrator	604000	pg/mL
		Neurodegeneration Control 1	10100	pg/mL
		Neurodegeneration Control 2	3290	pg/mL
		Neurodegeneration Control 3	571	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:**

Aβ-40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
15100	704,000	5.2
3780	208,000	5.4
944	35,200	6.1
236	3,930	9.5
59.0	585	9.6
14.7	179	13.0
3.69	110	13.0
0.00	108	35.0

**CERTIFICATE OF ANALYSIS**

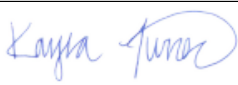
**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056400	2-8°C	30 Apr 2021
Anti-Aβ 6E10 Antibody	D0081076	2-8°C	31 Oct 2021
Aβ1-40 Peptide	A0080190	≤ -70°C	30 Apr 2021
Aβ40 Blocker	R0110016	2-8°C	30 Sep 2020
Diluent 35	M0350029	2-8°C	30 Sep 2020
Diluent 100	Y0270255	2-8°C	30 Sep 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), 100mL	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>	Kayla Turner	Quality Assurance		23 Jan 2019