



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-03723

**Product Description:** A $\beta$ 40 Peptide (4G8) Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K150SJE-1, K150SJE-2, K150SJE-4, K150SJG-1, K150SJG-2, K150SJG-4

**Kit Lot Number:** K0040191

**Expiration Date:** 30SEP2015

### Acceptance Data:

Metric		Specifications	Assay A $\beta$ 40
% Recovery Concentration	Calibrators	80% - 120%	PASS
	Controls	75% - 125%	PASS
	LLOQ	80% - 120%	PASS
	ULOQ	80% - 120%	PASS
% Recovery CV	Calibrators	<20%	PASS
	Controls	<20%	PASS
	LLOQ	<20%	PASS
	ULOQ	<20%	PASS
Non-specific binding		<1.0%	PASS

### Stock Concentration of Materials Tested:

Units = pg/mL	Assay A $\beta$ 40
Calibrator	582 000
Control 1 (A00C0113)	9071
Control 2 (A00C0114)	3036
Control 3 (A00C0115)	570
Control 4 (A00C0145)	8963
Control 5 (A00C0146)	3019
Control 6 (A00C0147)	1870
ULOQ	6000
LLOQ	20.0

### Standard Curve Data from Kit QC:

A $\beta$ 40		
Conc., (pg/ml)	Average Signal	%CV
0	158	4.1
3.55	189	5.1
14.2	290	4.8
56.8	839	6.3
227	4646	3.6
909	34 187	2.8
3638	167 265	3.8
14 550	490 176	2.7




## CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-03723

A $\beta$ 40 Peptide (4G8) Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ 40 Peptide Plate	Z0054560	2-8°C	31OCT2015
Anti-A $\beta$ 4G8 Antibody	D0040351	2-8°C	30NOV2015
A $\beta$ 1-40 Peptide	A0040071	$\leq -70^{\circ}\text{C}$	31OCT2015
A $\beta$ 40 Blocker	R0110003	$\leq -70^{\circ}\text{C}$	31OCT2015
Diluent 35	M0350011	2-8°C	31OCT2015
Diluent 100	Y0270188	2-8°C	30SEP2015
Read Buffer	N/A	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100 mL	Not Kit Specific	Room Temperature	N/A
Plate Seals, pack of 3	Not Kit Specific	Room Temperature	N/A
Neurodegeneration Control 1	Not Kit Specific; See Separate Control Pack CoA		
Neurodegeneration Control 2			
Neurodegeneration Control 3			

**Statement:** The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Paula Eason	Quality Control		23SEP2013