

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: K0080899
Expiration Date: 31 Aug 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	608000	pg/mL
		Neurodegeneration Control 1	9220	pg/mL
		Neurodegeneration Control 2	3550	pg/mL
		Neurodegeneration Control 3	548	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
15200	565,000	5.5
3800	192,000	8.9
950	38,400	5.0
238	4,660	3.7
59.4	770	4.4
14.8	217	4.0
3.71	115	7.1
0.00	85	4.6

Components:

CERTIFICATE OF ANALYSIS


Page 2 of 2

Description	Lot Number	Storage Temperature	Expiration Date
A β 40 Peptide Plate	Z0056077	2-8°C	31 Mar 2020
Anti-A β 4G8 Antibody	D0080947	2-8°C	30 Sep 2019
A β 1-40 Peptide	A0080129	\leq -70°C	30 Sep 2019
A β 40 Blocker	R0110015	2-8°C	31 Dec 2019
Diluent 35	M0350028	2-8°C	31 Dec 2019
Diluent 100	Y0270248	2-8°C	31 Aug 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X)	Not Kit Specific
Plate Seal	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	Dawn Stollar	Quality Assurance		01 NOV 2017