

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: K0080934
Expiration Date: 31 Jul 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	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	608000	pg/mL
		Neurodegeneration Control 1	10200	pg/mL
		Neurodegeneration Control 2	3720	pg/mL
		Neurodegeneration Control 3	555	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
15200	705,000	3.1
3800	213,000	1.8
950	36,000	3.6
238	3,900	3.6
59.4	526	2.5
14.8	164	4.5
3.71	104	5.2
0.00	92	9.7

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056130	2-8°C	30 Jun 2020
Anti-Aβ 6E10 Antibody	D0080883	2-8°C	31 Jul 2019
Aβ1-40 Peptide	A0080143	≤ -70°C	31 Dec 2019
Aβ40 Blocker	R0110015	2-8°C	31 Dec 2019
Diluent 35	M0350028	2-8°C	31 Dec 2019
Diluent 100	Y0270249	2-8°C	31 Oct 2019
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), 1000mL	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		31 JAN 2018