

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human A β 42 Kit
Kit Catalog Numbers: K151LB-Series
Kit Lot Number: K0080802
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	A β 1-42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	A β 1-42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	A β 1-42	20X Calibrator	91200	pg/mL
		Neurodegeneration Control 1	1470	pg/mL
		Neurodegeneration Control 2	494	pg/mL
		Neurodegeneration Control 3	207	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:

A β 1-42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
4560	781,000	4.9
1140	310,000	8.3
285	74,500	6.9
71.3	12,300	8.8
17.8	2,010	6.0
4.45	463	7.4
1.11	170	11.0
0.278	83	8.1
0.00	81	4.5

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ42 Peptide Plate	Z0055924	2-8°C	31 Oct 2019
Anti-Aβ 6E10 Antibody	D0080970	2-8°C	31 Aug 2019
Aβ1-42 Peptide	A0080130	≤ -70°C	30 Sep 2019
Diluent 35	M0350026	2-8°C	31 Jul 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), 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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	21 JUN 2017