

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: K0081325
Expiration Date: 31 May 2021

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 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 42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	A β 42	20X Calibrator	70000	pg/mL
		Neurodegeneration Control 1	1440	pg/mL
		Neurodegeneration Control 2	506	pg/mL
		Neurodegeneration Control 3	203	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 42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3500	1,090,000	3.8
875	412,000	4.5
219	100,000	6.3
54.7	17,700	8.4
13.7	3,090	6.5
3.42	697	7.1
0.854	227	4.9
0.214	117	9.6
0.00	101	8.3

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ42 Peptide Plate	Z0056583	2-8°C	28 Feb 2022
Anti-Aβ 6E10 Antibody	D0081076	2-8°C	31 Oct 2021
Aβ1-42 Peptide	A0080200	≤ -70°C	30 Jun 2021
Diluent 35	M0350030	2-8°C	31 May 2021
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	Karen Hamilla	Quality		19 DEC 2019