

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: K0080679
Expiration Date: 30 Nov 2018

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	hu 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	hu A β 42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu A β 42	20X Calibrator	69200	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

CERTIFICATE OF ANALYSIS

Page 2 of 2

Standard Curve Data:

hu A β 42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3460	714,000	10.0
865	268,000	7.2
216	63,400	11.0
54.1	10,700	11.0
13.5	1,890	16.0
3.38	467	7.9
0.845	167	5.5
0.211	104	4.5
0.00	83	6.8


Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β 42 Peptide Plate	Z0055864	2-8°C	31 Aug 2019
Anti-A β 6E10 Antibody	D0080970	2-8°C	31 Aug 2019
A β 1-42 Peptide	A0040169	\leq -70°C	30 Nov 2018
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)	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		28 JUN 2017