

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Aβ42 Peptide (4G8) Kit
Kit Catalog Numbers: K150SL-Series; K151LB-Series
Kit Lot Number: K0081342
Expiration Date: 31 May 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ1-42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ1-42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ1-42	40X Calibrator	99200	pg/mL
		Neurodegeneration Control 1 (4G8)	1410	pg/mL
		Neurodegeneration Control 2 (4G8)	515	pg/mL
		Neurodegeneration Control 3 (4G8)	211	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:

Aβ1-42 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2480	1,170,000	1.4
620	334,000	3.1
155	52,800	5.8
38.8	6,270	4.5
9.69	896	3.2
2.42	197	7.2
0.605	100	5.0
0.00	73	3.5

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ42 Peptide Plate	Z0056598	2-8°C	31 Mar 2022
Anti-Aβ 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
Aβ1-42 Peptide (4G8)	A0080201	≤ -70°C	31 Jul 2021
Diluent 35	M0350030	2-8°C	31 May 2021
Diluent 100	Not Kit Specific	Not Kit Specific	N/A
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 (4G8)	Not Kit Specific
Neurodegeneration Control 2 (4G8)	Not Kit Specific
Neurodegeneration Control 3 (4G8)	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		11 DEC 2019