

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Aβ40 Peptide (4G8) Kit
Kit Catalog Numbers: K150SJ-Series
Kit Lot Number: K0081312
Expiration Date: 30 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	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ40	40X Calibrator	584000	pg/mL
		Neurodegeneration Control 1 (4G8)	8040	pg/mL
		Neurodegeneration Control 2 (4G8)	2700	pg/mL
		Neurodegeneration Control 3 (4G8)	445	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:

Aβ40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
14600	871,000	2.7
3650	299,000	3.5
913	61,200	5.6
228	7,680	5.9
57.0	1,090	4.4
14.3	260	3.5
3.56	124	4.9
0.00	84	7.5

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056576	2-8°C	31 Jan 2022
Anti-Aβ 4G8 Antibody	D0081102	2-8°C	30 Nov 2021
Aβ1-40 Peptide (4G8)	A0080202	≤ -70°C	30 Sep 2021
Aβ40 Blocker	R0010657	2-8°C	30 Sep 2021
Diluent 35	M0350030	2-8°C	30 May 2021
Diluent 100	Y0270260	2-8°C	30 Jun 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
Was Buffer (20x), 100 ml	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 OCT 2019