

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8) Kit

Kit Catalog Numbers: K150SJ-Series, K150SK-Series

Kit Lot Number: K0081430

Expiration Date: 30 Sep 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	843,000	8.2
3650	301,000	3.2
913	60,000	3.5
228	7,400	4.3
57.0	1,260	2.8
14.3	404	3.5
3.56	261	2.8
0.00	220	3.1

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056667	2-8°C	31 Jul 2022
Anti-Aβ 4G8 Antibody	D0081432	2-8°C	31 Aug 2023
Aβ1-40 Peptide (4G8)	A0080202	≤ -70°C	30 Sep 2021
Aβ40 Blocker	R0010692	2-8°C	30 Sep 2022
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270270	2-8°C	30 Jun 2022
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		15 JUN 2020