

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus A β 40 Peptide (6E10) Kit
Kit Catalog Numbers: K150SK-Series
Kit Lot Number: K0081335
Expiration Date: 30 Sep 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 40	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 40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	A β 40	40X Calibrator	600000	pg/mL
		Neurodegeneration Control 1	9430	pg/mL
		Neurodegeneration Control 2	3190	pg/mL
		Neurodegeneration Control 3	534	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 40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
15000	836,000	3.5
3750	268,000	7.9
938	51,800	12.0
234	6,000	15.0
58.6	642	10.0
14.6	145	5.0
3.66	84	6.4
0.00	67	3.6

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056593	2-8°C	31 Mar 2022
Anti-Aβ 6E10 Antibody	D0081119	2-8°C	31 Jan 2022
Aβ1-40 Peptide (6E10)	A0080203	≤ -70°C	31 May 2022
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270266	2-8°C	28 Feb 2022
Aβ40 Blocker	R0010657	2-8°C	30 Sep 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		06 MAR 2020