



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus A β 40 Peptide (6E10) Kit

Catalog Number(s) / Size(s): K150SK-Series

Lot Number: K0080966

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080966 has been extended to 30 JUN 2020.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
A β 1-40 Peptide	A0080118	30 JUN 2019	30 JUN 2020
A β 40 Blocker	R0110015	31 DEC 2019	30 JUN 2020
Diluent 35	M0350028	31 DEC 2019	30 JUN 2020
Diluent 100	Y0270249	31 OCT 2019	30 JUN 2020
Anti-A β 6E10 Antibody	D0080883	31 JUL 2019	30 JUN 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		12 JUL 2019

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Aβ 40 Peptide (6E10) Kit
Kit Catalog Numbers: K150SK-Series
Kit Lot Number: K0080966
Expiration Date: 30 Jun 2019

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	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	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Aβ40	40X Calibrator	604000	pg/mL
		Neurodegeneration Control 1	9340	pg/mL
		Neurodegeneration Control 2	2890	pg/mL
		Neurodegeneration Control 3	455	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
15100	994,000	2.7
3780	325,000	4.3
944	57,800	5.4
236	6,440	7.9
59.0	807	4.1
14.7	199	2.7
3.69	110	4.9
0.00	82	5.3

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056181	2-8°C	31 Aug 2020
Anti-Aβ 6E10 Antibody	D0080883	2-8°C	31 Jul 2019
Aβ1-40 Peptide	A0080118	≤ -70°C	30 Jun 2019
Aβ40 Blocker	R0110015	2-8°C	31 Dec 2019
Diluent 35	M0350028	2-8°C	31 Dec 2019
Diluent 100	Y0270249	2-8°C	31 Oct 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), 100 mL	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	<i>Dawn Stollar</i>	23 MAY 2018