



RFE MEMO

Page 1 of 1

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


Catalog Number(s) / Size(s): K150SJ-Series

Lot Number: K0081734

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081734 has been extended to 30 SEP 2022.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Aβ40 Peptide Plate	Z0056621	30 APR 2022	30 SEP 2022
Aβ1-40 Peptide (4G8)	A0080202	30 SEP 2021	30 SEP 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		28 MAR 2022

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

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8) Kit
Kit Lot Number: K0081734
Expiration Date: 30 Sep 2021
Kit Catalog Numbers: K150SJ-Series

Precision and Accuracy Specifications:

		Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

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

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Aβ40	584000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0056621	2-8°C	30 Apr 2022
Anti-Aβ 4G8 Antibody	D0081432	2-8°C	31 Aug 2023
Aβ1-40 Peptide (4G8)	A0080202	≤ -70°C	30 Sep 2021
Aβ40 Blocker	R0010728	2-8°C	31 Aug 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023
Diluent 100	Y0270290	2-8°C	31 Jul 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		21 MAY 2021