

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Aβ40 Peptide (6E10) Kit
Kit Lot Number: K0081697
Expiration Date: 31 Jan 2022
Kit Catalog Numbers: K150SK-Series

Precision and Accuracy Specifications:

		Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
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	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Aβ40	600000	pg/mL

Lot-Specific 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
Aβ40 Blocker	R0010692	2-8°C	30 Sep 2022
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270266	2-8°C	28 Feb 2022

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	Karen Hamilla	Quality		07 APR 2021