

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Aβ40 Peptide (6E10) Kit  
**Kit Lot Number:** K0081997  
**Expiration Date:** 31 Mar 2023  
**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	7000	80% - 120%	< 20%	50.0	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	Aβ40	612000	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide Plate	Z0057151	2-8°C	30 Sep 2024
Anti-Aβ 6E10 Antibody	D0081630	2-8°C	31 Jul 2024
Aβ1-40 Peptide (6E10)	A0080277	≤ -70°C	30 Sep 2023
Aβ40 Blocker	R0010728	2-8°C	31 Aug 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023
Diluent 100	Y0270309	2-8°C	31 Aug 2024

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		03 AUG 2022