

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Ab40 Peptide (6E10) Kit  
**Lot Number:** K0082554  
**Expiration Date:** 31 MAR 2025  
**Catalog Number:** K150SK-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A $\beta$ 6E10 Antibody	D0081958	2-8°C	31 JUL 2026
A $\beta$ 40 Blocker	R0010817	2-8°C	31 MAR 2025
A $\beta$ 40 Peptide Plate	Z0057297	2-8°C	31 MAY 2025
Diluent 100	Y0270334	2-8°C	30 SEP 2026
Diluent 35	M0350037	2-8°C	31 DEC 2025
A $\beta$ 1-40 Peptide (6E10)	A0080349	$\leq$ -70°C	31 MAY 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot	Name	Calibrator		Neurodegeneration Control 1 (6E10)		Neurodegeneration Control 2 (6E10)		Neurodegeneration Control 3 (6E10)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A $\beta$ 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 $\beta$ 40	50	80% - 120%	< 20%	7000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	A $\beta$ 40	624000	pg/mL

Dilution to top of Curve: 40X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		16 SEP 2024