



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Ab40 Peptide (6E10) Kit
Lot Number: K0082592
Expiration Date: 30 SEP 2026
Catalog Number: K150SK-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A β 6E10 Antibody	D0082325	2-8°C	31 MAR 2028
A β 40 Peptide 96-Well 4-Spot SECTOR Plate	Z0057531	2-8°C	30 APR 2027
Diluent 100	Y0270342	2-8°C	31 MAR 2027
Diluent 35	M0350038	2-8°C	31 OCT 2026
A β 1-40 Peptide (6E10)	A0080408	$\leq -70^{\circ}\text{C}$	31 MAR 2027
A β 40 Blocker	R0010896	2-8°C	30 SEP 2026

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Neurodegeneration Control 1 (6E10)		Neurodegeneration Control 2 (6E10)		Neurodegeneration Control 3 (6E10)	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	A β 40	50	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	580000	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
Review/Approval	Karen Hamilla	Quality		30 JAN 2025