

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Ab40 Peptide (6E10) Kit**Lot Number:** K0082743**Expiration Date:** 31 MAR 2027**Catalog Number:** K150SK-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ40 Peptide 96-Well 4-Spot SECTOR Plate	Z0057562	2-8°C	31 AUG 2027
Aβ40 Blocker	R0010925	2-8°C	30 SEP 2027
Diluent 100	Y0270344	2-8°C	30 APR 2027
Diluent 35	M0350039	2-8°C	30 JUN 2027
Human Aβ 6E10 Antibody	D0082394	2-8°C	31 MAY 2028
Aβ1-40 Peptide (6E10)	A0080408	≤ -70°C	31 MAR 2027

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	AB40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	AB40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	AB40	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		07 JUL 2025