



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX Ab40 Peptide (4G8) Kit  
**Lot Number:** K0082584  
**Expiration Date:** 28 FEB 2025  
**Catalog Number:** K150SJ-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A $\beta$ 4G8 Antibody	D0081873	2-8°C	31 DEC 2025
A $\beta$ 40 Peptide 96-Well 4-Spot SECTOR Plate	Z0057252	2-8°C	28 FEB 2025
A $\beta$ 40 Blocker	R0010829	2-8°C	30 JUN 2025
Diluent 100	Y0270334	2-8°C	30 SEP 2026
Diluent 35	M0350037	2-8°C	31 DEC 2025
A $\beta$ 1-40 Peptide (4G8)	A0080359	$\leq$ -70°C	31 JUL 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 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		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	20	80% - 120%	< 20%	6000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	A $\beta$ 40	616000	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		12 NOV 2024