

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX and V-PLEX Plus A β 42 Peptide (4G8) Kit
Lot Number: K0082476
Expiration Date: 28 FEB 2025
Catalog Number: K150SL-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A β 4G8 Antibody	D0081899	2-8°C	28 FEB 2026
A β 42 Peptide Plate	Z0057246	2-8°C	28 FEB 2025
Diluent 35	M0350036	2-8°C	31 JUL 2025
A β 1-42 Peptide (4G8)	A0080393	\leq -70°C	30 SEP 2026

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 β 42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%


Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	A β 42	2.5	80% - 120%	< 20%	1270	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 42	80400	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		12 JUN 2024