



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8) Kit
Lot Number: K0082585
Expiration Date: 28 FEB 2025
Catalog Number: K150SJ-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-Aβ 4G8 Antibody	D0081873	2-8°C	31 DEC 2025
Aβ40 Peptide Plate	Z0057252	2-8°C	28 FEB 2025
Aβ40 Blocker	R0010896	2-8°C	30 SEP 2026
Diluent 100	Y0270334	2-8°C	30 SEP 2026
Diluent 35	M0350037	2-8°C	31 DEC 2025
Aβ1-40 Peptide (4G8)	A0080359	≤ -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	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	20	80% - 120%	< 20%	6000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Aβ40	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
Review/Approval	Michelle Likiam	Quality		27 NOV 2024