

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Aβ40 Peptide (4G8) Kit**Lot Number:** K0082631**Expiration Date:** 30 SEP 2026**Catalog Number:** K150SJ-Series, K150SK-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-Aβ 4G8 Antibody	D0082222	2-8°C	31 AUG 2027
Aβ40 Peptide 96-Well 4-Spot SECTOR Plate	Z0057548	2-8°C	31 JUL 2027
Diluent 100	Y0270342	2-8°C	31 MAR 2027
Diluent 35	M0350039	2-8°C	30 JUN 2027
Aβ1-40 Peptide (4G8)	A0080392	≤ -70°C	30 SEP 2026
Aβ40 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:

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	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Aβ40	656000	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		31 MAR 2025



CERTIFICATE OF ANALYSIS