



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus A β 40 Peptide (4G8) Kit**Lot Number:** K0082706**Expiration Date:** 30 APR 2027**Catalog Number:** K150SJ-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
A β 40 Peptide 96-Well 4-Spot SECTOR Plate	Z0057575	2-8°C	31 OCT 2027
Diluent 100	Y0270344	2-8°C	30 APR 2027
Diluent 35	M0350039	2-8°C	30 JUN 2027
Anti-A β 4G8 Antibody	D0082222	2-8°C	31 AUG 2027
A β 1-40 Peptide (4G8)	A0080425	$\leq -70^{\circ}\text{C}$	31 AUG 2027
A β 40 Blocker	R0010910	2-8°C	30 APR 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	A β 40	20	80% - 120%	< 20%	6000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 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
Review/Approval	Karen Hamilla	Quality		18 JUN 2025