



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus A β 40 Peptide (4G8) Kit, QuickPlex
Lot Number: K0082986
Expiration Date: 31 DEC 2027
Catalog Number: K150AYS-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
A β 4G8 Antibody	D0082580	2-8°C	31 JUL 2029
A β 40 Peptide (4G8) 96-Well 10-Spot QuickPlexUltra Plate	Z0057728	2-8°C	31 DEC 2028
Diluent 100	Y0270356	2-8°C	31 MAR 2028
Diluent 35	M0350041	2-8°C	31 DEC 2027
A β 40 Blocker	R0010958	2-8°C	30 APR 2028
A β 1-40 Peptide (4G8)	A0080449	$\leq -70^{\circ}\text{C}$	31 AUG 2028

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

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	576000	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		28 JUL 2026