



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus A β 40 Peptide (6E10) Kit, QuickPlex
Lot Number: K0082987
Expiration Date: 30 APR 2028
Catalog Number: K150AYT-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 40 Blocker	R0010958	2-8°C	30 APR 2028
A β 40 Peptide (6E10) 96-Well 10-Spot QuickPlexUltra Plate	Z0057729	2-8°C	31 DEC 2028
Diluent 100	Y0270361	2-8°C	30 SEP 2028
Diluent 35	M0350042	2-8°C	31 AUG 2028
Human A β 6E10Antibody	D0082579	2-8°C	31 JUL 2029
A β 1-40 Peptide (6E10)	A0080452	$\leq -70^{\circ}\text{C}$	30 APR 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 (6E10)		Neurodegeneration Control 2 (6E10)		Neurodegeneration Control 3 (6E10)	
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	50	80% - 120%	< 20%	7000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	596000	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		07 AUG 2026