



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus A β Peptide Panel 2 (6E10) QuickPlex Kit**Lot Number:** K0082903**Expiration Date:** 30 NOV 2027**Catalog Number:** K15765-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 Peptide Panel 2 (6E10) 96-Well 10-Spot QuickPlexUltra Plate	Z0057669	2-8°C	31 AUG 2028
Diluent 100	Y0270355	2-8°C	29 FEB 2028
Diluent 35	M0350040	2-8°C	30 NOV 2027
Human A β 6E10 Antibody	D0082579	2-8°C	31 JUL 2029
A β 1-42 Peptide (6E10)	A0080453	$\leq -70^{\circ}\text{C}$	30 JUN 2028
A β 1-38 Peptide (6E10)	A0080451	$\leq -70^{\circ}\text{C}$	30 APR 2028
A β 1-40 Peptide (6E10)	A0080452	$\leq -70^{\circ}\text{C}$	30 APR 2028
A β 40 Blocker	R0010958	2-8°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%
3	A β 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
10	A β 42	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	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
3	A β 38	60.0	80% - 120%	< 20%	8480	75% - 125%	< 20%
10	A β 42	3.13	75% - 125%	< 20%	1270	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	Spec
1	A β 40	< 1.0%
3	A β 38	< 1.0%
10	A β 42	< 1.0%



CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	596000	pg/mL
3	A β 38	552000	pg/mL
10	A β 42	73200	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		26 JUL 2026