

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Ab Peptide Panel 1 (6E10) Kit**Lot Number:** K0082932**Expiration Date:** 31 DEC 2027**Catalog Number:** K15200-Series, K150SI-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
Anti-A β 6E10 Antibody	D0082484	2-8°C	28 FEB 2029
A β Peptide Panel 1 (6E10) 96-Well 4-Spot SECTOR Plate	Z0057680	2-8°C	30 SEP 2028
A β 40 Blocker	R0010990	2-8°C	31 OCT 2028
Diluent 100	Y0270356	2-8°C	31 MAR 2028
Diluent 35	M0350041	2-8°C	31 DEC 2027
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

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 (6E10)		Neurodegeneration Control 2 (6E10)		Neurodegeneration Control 3 (6E10)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	A β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A β 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A β 42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	A β 40	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	A β 38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	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%
2	A β 38	< 1.0%
4	A β 42	< 1.0%



CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	596000	pg/mL
2	A β 38	552000	pg/mL
4	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		28 JUL 2026