

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX, and V-PLEX Plus Human A $\beta$ 42 Kit**Lot Number:** K0082865**Expiration Date:** 28 FEB 2027**Catalog Number:** K151LB-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 $\beta$ 6E10 Antibody	D0082570	2-8°C	30 JUN 2029
AB42 Peptide 96-Well 4-Spot SECTOR Plate	Z0057646	2-8°C	30 JUN 2028
Diluent 35	M0350040	2-8°C	30 NOV 2027
AB1-42 Peptide (6E10)	A0080409	$\leq$ -70°C	28 FEB 2027

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 $\beta$ 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 $\beta$ 42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	A $\beta$ 42	72800	pg/mL

Dilution to top of Curve: 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		20 FEB 2026