



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human A β 42 Kit**Lot Number:** K0082605**Expiration Date:** 28 FEB 2027**Catalog Number:** K151LB-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A β 6E10 Antibody	D0082394	2-8°C	31 MAY 2028
A β 42 Peptide 96-Well 4-Spot SECTOR Plate	Z0057541	2-8°C	30 JUN 2027
Diluent 35	M0350039	2-8°C	30 JUN 2027
A β 1-42 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:

		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 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 42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 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
Review/Approval	Karen Hamilla	Quality		21 MAR 2025