

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human Aβ42 Kit  
**Kit Lot Number:** K0082100  
**Expiration Date:** 31 Mar 2023  
**Kit Catalog Numbers:** K151LB-Series

**Precision and Accuracy 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	AB42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	AB42	3.00	80% - 120%	< 20%	2000	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	AB42	68400	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
AB42 Peptide Plate	Z0057235	2-8°C	31 Jan 2025
Anti-Aβ 6E10 Antibody	D0081630	2-8°C	31 Jul 2024
AB1-42 Peptide (6E10)	A0080298	≤ -70°C	30 Nov 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 FEB 2024