

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human A β 42 Kit
Kit Lot Number: K0081758
Expiration Date: 31 Mar 2023
Kit Catalog Numbers: K151LB-Series; K150SL-Series

Precision and Accuracy Specifications:

		Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
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	68400	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β 42 Peptide Plate	Z0056970	2-8°C	30 Nov 2023
Anti-A β 6E10 Antibody	D0081509	2-8°C	30 Sep 2023
A β 1-42 Peptide (6E10)	A0080298	\leq -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		08 SEP 2021