

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Aβ42 Peptide (4G8) Kit
Kit Lot Number: K0081759
Expiration Date: 31 Mar 2023
Kit Catalog Numbers: K150SL-Series; K151LB-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

Concentration of Calibrator:


Spot	Name	Conc.	Units
1	Aβ42	74800	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ42 Peptide Plate	Z0056971	2-8°C	30 Nov 2023
Anti-Aβ 4G8 Antibody	D0081432	2-8°C	31 Aug 2023
Aβ1-42 Peptide (4G8)	A0080294	≤ -70°C	31 Oct 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