

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Aβ Peptide Panel 1 (4G8) Kit

**Kit Lot Number:** K0081844

**Expiration Date:** 31 Mar 2023

**Kit Catalog Numbers:** K15199-Series; K150SH-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Aβ-40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Aβ-38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	Aβ-42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ-40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	Aβ-38	60.0	75% - 125%	< 20%	7500	80% - 120%	< 20%
4	Aβ-42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

Spot	Name	NSB Spec
1	Aβ-40	< 1.0%
2	Aβ-38	< 1.0%
4	Aβ-42	< 1.0%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	Aβ-40	652000	pg/mL
2	Aβ-38	472000	pg/mL
4	Aβ-42	74800	pg/mL


## CERTIFICATE OF ANALYSIS

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 (4G8) Plate	Z0057047	2-8°C	29 Feb 2024
Anti-Aβ 4G8 Antibody	D0081743	2-8°C	30 Apr 2025
Aβ1-38 Peptide (4G8)	A0080292	≤ -70°C	31 Oct 2023
Aβ1-40 Peptide (4G8)	A0080293	≤ -70°C	31 Oct 2023
Aβ1-42 Peptide (4G8)	A0080294	≤ -70°C	31 Oct 2023
Aβ 40 Blocker	R0010728	2-8°C	31 Aug 2023
Diluent 35	M0350032	2-8°C	31 Mar 2023
Diluent 100	Y0270302	2-8°C	29 Feb 2024

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	Xiomara Talavera	Quality		22 DEC 2021