

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus A $\beta$  Peptide Panel 1 (4G8) Kit

**Kit Lot Number:** K0082066

**Expiration Date:** 31 Oct 2023

**Kit Catalog Numbers:** K15199-Series; K150SH-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	A $\beta$ 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A $\beta$ 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A $\beta$ 42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

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

**Non-Specific Binding Specifications:**

Spot	Name	NSB Spec
1	A $\beta$ 40	< 1.0%
2	A $\beta$ 38	< 1.0%
4	A $\beta$ 42	< 1.0%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	A $\beta$ 40	584000	pg/mL
2	A $\beta$ 38	544000	pg/mL
4	A $\beta$ 42	74800	pg/mL


**CERTIFICATE OF ANALYSIS**

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ Peptide Panel 1 (4G8) Plate	Z0057208	2-8°C	31 Dec 2024
Anti-A $\beta$ 4G8 Antibody	D0081873	2-8°C	31 Dec 2025
A $\beta$ 1-38 Peptide (4G8)	A0080335	$\leq$ -70°C	31 Jan 2025
A $\beta$ 1-40 Peptide (4G8)	A0080336	$\leq$ -70°C	31 Dec 2024
A $\beta$ 1-42 Peptide (4G8)	A0080294	$\leq$ -70°C	31 Oct 2023
A $\beta$ 40 Blocker	R0010797	2-8°C	31 Oct 2024
Diluent 100	Y0270313	2-8°C	30 Nov 2024
Diluent 35	M0350033	2-8°C	31 Oct 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	Karen Hamilla	Quality		29 SEP 2022