



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus A $\beta$  Peptide Panel 1 (4G8) Kit  
**Lot Number:** K0082439  
**Expiration Date:** 31 DEC 2025  
**Catalog Number:** K15199-Series; K150SH-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A $\beta$ 4G8 Antibody	D0082222	2-8°C	31 AUG 2027
A $\beta$ Peptide Panel 1 (4G8) Plate	Z0057456	2-8°C	30 SEP 2026
A $\beta$ 40 Blocker	R0010896	2-8°C	30 SEP 2026
Diluent 100	Y0270334	2-8°C	30 SEP 2026
Diluent 35	M0350037	2-8°C	31 DEC 2025
A $\beta$ 1-38 Peptide (4G8)	A0080391	$\leq$ -70°C	30 SEP 2026
A $\beta$ 1-40 Peptide (4G8)	A0080392	$\leq$ -70°C	30 SEP 2026
A $\beta$ 1-42 Peptide (4G8)	A0080393	$\leq$ -70°C	30 SEP 2026

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

**Precision, Accuracy, and Sensitivity 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	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	A $\beta$ 38	60	75% - 125%	< 20%	7500	80% - 120%	< 20%
4	A $\beta$ 42	2.5	80% - 120%	< 20%	1270	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

Spot	Name	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	656000	pg/mL
2	A $\beta$ 38	592000	pg/mL
4	A $\beta$ 42	80400	pg/mL

Dilution to top of Curve: 40X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		25 JUN 2024