

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Aβ Peptide Panel 1 (4G8) Kit
Kit Lot Number: K0082376
Expiration Date: 30 JUN 2025
Kit Catalog Numbers: K15199-Series, K150SH-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Aβ Peptide Panel 1 (4G8) Plate	Z0057410	2-8°C	28 FEB 2026
Anti-Aβ 4G8 Antibody	D0081965	2-8°C	30 JUN 2026
Aβ1-38 Peptide (4G8)	A0080358	≤ -70°C	31 JUL 2025
Aβ1-40 Peptide (4G8)	A0080359	≤ -70°C	31 JUL 2025
Aβ1-42 Peptide (4G8)	A0080360	≤ -70°C	30 SEP 2025
Aβ40 Blocker	R0010829	2-8°C	30 JUN 2025
Diluent 100	Y0270330	2-8°C	30 SEP 2025
Diluent 35	M0350036	2-8°C	31 JUL 2025

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

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	Aβ40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Aβ38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	Aβ42	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Aβ40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	Aβ38	60.0	75% - 125%	< 20%	7500	80% - 120%	< 20%
4	Aβ42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

Non-Specific Binding Specifications:


Spot	Name	NSB Spec
1	Aβ40	< 1.0%
2	Aβ38	< 1.0%
4	Aβ42	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Aβ40	616000	pg/mL
2	Aβ38	528000	pg/mL
4	Aβ42	79600	pg/mL

Dilution to Top of Curve = 40X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 NOV 2023