

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus A β Peptide Panel 1 (4G8) Kit
Kit Lot Number: K0082150
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 β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	A β 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	A β 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 β 40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	A β 38	60.0	75% - 125%	< 20%	7500	80% - 120%	< 20%
4	A β 42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB Spec
1	A β 40	< 1.0%
2	A β 38	< 1.0%
4	A β 42	< 1.0%

Stock Concentration of Calibrator:


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

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 (4G8) Plate	Z0057270	2-8°C	31 Mar 2025
Anti-A β 4G8 Antibody	D0081899	2-8°C	28 Feb 2026
A β 1-38 Peptide (4G8)	A0080335	\leq -70°C	31 Jan 2025
A β 1-40 Peptide (4G8)	A0080336	\leq -70°C	31 Dec 2024
A β 1-42 Peptide (4G8)	A0080294	\leq -70°C	31 Oct 2023
A β 40 Blocker	R0010817	2-8°C	31 Mar 2025
Diluent 100	Y0270315	2-8°C	31 Dec 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		07 DEC 2022