



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus A β Peptide Panel 1 (4G8) Kit
Lot Number: K0082528
Expiration Date: 30 SEP 2026
Catalog Number: K15199-Series; K150SH-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-A β 4G8 Antibody	D0082302	2-8°C	30 APR 2028
A β Peptide Panel 1 (4G8) 96-Well 4-Spot SECTOR Plate	Z0057488	2-8°C	31 JAN 2027
A β 40 Blocker	R0010896	2-8°C	30 SEP 2026
Diluent 100	Y0270335	2-8°C	30 NOV 2026
Diluent 35	M0350038	2-8°C	31 OCT 2026
A β 1-38 Peptide (4G8)	A0080391	$\leq -70^{\circ}\text{C}$	30 SEP 2026
A β 1-40 Peptide (4G8)	A0080392	$\leq -70^{\circ}\text{C}$	30 SEP 2026
A β 1-42 Peptide (4G8)	A0080393	$\leq -70^{\circ}\text{C}$	30 SEP 2026

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
Spot	Name	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%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	A β 40	20	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	A β 38	60	75% - 125%	< 20%	7500	80% - 120%	< 20%
4	A β 42	2.5	80% - 120%	< 20%	1270	80% - 120%	< 20%

Non-Specific Binding Specifications:

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



CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	656000	pg/mL
2	A β 38	592000	pg/mL
4	A β 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		09 DEC 2024