

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus A β Peptide Panel 2 (4G8) SECTOR Kit**Lot Number:** K0082905**Expiration Date:** 30 NOV 2027**Catalog Number:** K15764-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
A β 4G8 Antibody	D0082580	2-8°C	31 JUL 2029
A β Peptide Panel 2 (4G8) 96-Well 10-Spot SECTOR Plate	Z0057671	2-8°C	31 AUG 2028
Diluent 100	Y0270355	2-8°C	29 FEB 2028
Diluent 35	M0350040	2-8°C	30 NOV 2027
A β 1-38 Peptide (4G8)	A0080448	$\leq -70^{\circ}\text{C}$	30 JUN 2028
A β 1-40 Peptide (4G8)	A0080449	$\leq -70^{\circ}\text{C}$	31 AUG 2028
A β 1-42 Peptide (4G8)	A0080450	$\leq -70^{\circ}\text{C}$	30 JUN 2028
A β 40 Blocker	R0010958	2-8°C	30 APR 2028

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 β 40	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	A β 38	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
10	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%
3	A β 38	60.0	75% - 125%	< 20%	7500	75% - 125%	< 20%
10	A β 42	2.5.0	80% - 120%	< 20%	1270	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	Spec
1	A β 40	< 1.0%
3	A β 38	< 1.0%
10	A β 42	< 1.0%



CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	A β 40	576000	pg/mL
3	A β 38	504000	pg/mL
10	A β 42	76400	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		28 JUL 2026