

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus A $\beta$  Peptide Panel 1 (4G8) Kit  
**Kit Lot Number:** K0081632  
**Expiration Date:** 31 Mar 2022  
**Kit Catalog Numbers:** K15199-Series; K150SH-Series

**Precision and Accuracy Specifications:**

		Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
Spot	Name	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%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	A $\beta$ 40	20.0	80% - 120%	< 20%	6000	80% - 120%	< 20%
2	A $\beta$ 38	60.0	75% - 125%	< 20%	7500	80% - 120%	< 20%
4	A $\beta$ 42	2.50	80% - 120%	< 20%	1270	80% - 120%	< 20%

**Non-Specific Binding Specifications:**

Spot	Name	NSB 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	584000	pg/mL
2	A $\beta$ 38	480000	pg/mL
4	A $\beta$ 42	99200	pg/mL


## CERTIFICATE OF ANALYSIS

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
A $\beta$ Peptide Panel 1 (4G8) Plate	Z0056856	2-8°C	30 Jun 2023
Anti-A $\beta$ 4G8 Antibody	D0081432	2-8°C	31 Aug 2023
A $\beta$ 1-38 Peptide (4G8)	A0080189	$\leq$ -70°C	30 Sep 2022
A $\beta$ 1-40 Peptide (4G8)	A0080202	$\leq$ -70°C	30 Sep 2022
A $\beta$ 1-42 Peptide (4G8)	A0080201	$\leq$ -70°C	31 May 2022
A $\beta$ 40 Blocker	R0010692	2-8°C	30 Sep 2022
Diluent 35	M0350031	2-8°C	31 Mar 2022
Diluent 100	Y0270270	2-8°C	30 Jun 2022

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		10 JUN 2021



**RFE MEMO**


Page 1 of 1

**Product Description(s):** A $\beta$ 1-38 Peptide (4G8)

**Catalog Number(s) / Size(s):** C002Y-2

**Lot Number:** A0080189

Based on additional testing, the original date of expiration for this calibrator has been extended by 12 months. The expiration date of A0080189 has been extended to 30 SEP 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		10 JUN 2021

This memo supersedes all other statements regarding date of expiration for this product.



**RFE MEMO**


Page 1 of 1

**Product Description(s):** A $\beta$ 1-42 Peptide (4G8)

**Catalog Number(s) / Size(s):** C003A-2

**Lot Number:** A0080201

Based on additional testing, the original date of expiration for this calibrator has been extended by 10 months. The expiration date of A0080201 has been extended to 31 MAY 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		08 APR 2021

This memo supersedes all other statements regarding date of expiration for this product.



**RFE MEMO**


Page 1 of 1

**Product Description(s):** A $\beta$ 1-40 Peptide (4G8)

**Catalog Number(s) / Size(s):** C002Z-2

**Lot Number:** A0080202

Based on additional testing, the original date of expiration for this calibrator has been extended by 12 months. The expiration date of A0080202 has been extended to 30 SEP 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		10 JUN 2021

This memo supersedes all other statements regarding date of expiration for this product.