

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus A β Peptide Panel 1 (6E10) Kit
Kit Lot Number: K0082067
Expiration Date: 31 May 2023
Kit Catalog Numbers: K15200-Series; K150SI-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3	
		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	50.0	80% - 120%	< 20%	7000	80% - 120%	< 20%
2	A β 38	60.0	80% - 120%	< 20%	8480	80% - 120%	< 20%
4	A β 42	3.13	75% - 125%	< 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	612000	pg/mL
2	A β 38	488000	pg/mL
4	A β 42	68400	pg/mL


CERTIFICATE OF ANALYSIS

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 (6E10) Plate	Z0056874	2-8°C	31 Jul 2023
Anti-A β 6E10 Antibody	D0081402	2-8°C	31 May 2023
A β 1-42 Peptide (6E10)	A0080298	\leq -70°C	30 Nov 2023
A β 1-38 Peptide (6E10)	A0080287	\leq -70°C	30 Sep 2023
A β 1-40 Peptide (6E10)	A0080277	\leq -70°C	30 Sep 2023
A β 40 Blocker	R0010728	2-8°C	31 Aug 2023
Diluent 100	Y0270312	2-8°C	31 Oct 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		14 JUL 2022