

Automation of the Enzo APP Δ C31 ELISA kit

APP Δ C31 ELISA kit (#ADI-900-227, Enzo)

Bettina Plommer¹, Erica Brooks², Miriam Cortes-Camirero³

1. Berthold Detection System GmbH (Titertek-Berthold), Pforzheim, Germany
2. Enzo Life Sciences, Ann Arbor, USA
3. Enzo Life Sciences, Farmingdale, USA

INTRODUCTION

Unlike the β -amyloid and tau fibril formation pathways leading to amyloid plaques and tangles, APP Δ C31 reveals an alternative, unique pro-apoptotic mechanism leading to Alzheimer's disease. APP Δ C31 is the stable amyloid precursor protein fragment created from a caspase cleavage event of the APP695 molecule at Asp664 leaving a smaller 31-residue intracellular fragment. Both the APP Δ C31 and 31-residue fragments are pro-apoptotic and are present in 4-fold greater levels in Alzheimer's disease patients. Given that the smaller 31-residue fragment has a short half-life and is difficult to measure, the use of this APP Δ C31 ELISA provides for the first time a sensitive research tool to measure the levels of the APP caspase cleavage from tissue, biological fluids and cells.

The APP Δ C31 ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of human APP Δ C31 in cell lysate, serum, plasma and cerebral spinal fluid samples with results in just 2 hours.

MATERIALS AND METHODS

Material Name	Company Name
Crocodile miniWorkstation	Titertek-Berthold
MikroWin module, quantitative and qualitative data reduction package	Titertek-Berthold
APP Δ C31 ELISA kit	Enzo Life Sciences
ddH ₂ O	
Pipette and tips	
Sample tubes	
Vortex Mixer	

Table 1: Materials provided with the Δ C31 ELISA kit from Enzo Life Sciences

Step	Component/Condition	Amount
1. Remove	All materials from cold source, allow to come to room temperature	
2. Prepare	Wash Buffer and 8 standards	
3. Transfer	Samples and standards in duplicate to assay plate	50 μ L
4. Follow Steps in Table 3	Dispense antibody	
5. Read	Finished plate at 450 nm absorbance	
6. Fit	Calculate standard curve using MikroWin and fit with the four-parameter algorithm	

Table 2: Procedure Overview for Δ C31 ELISA kit

RESULTS

Standard	Concentration (pM)	OD average (450nm) (Minus Blank OD = 0.063)
Blank	-	0
S0	0	-0.002
S1	1500	2.135
S2	750	1.242
S3	375	0.602
S4	187.5	0.293
S5	93.75	0.139
S6	46.88	0.091
S7	23.44	0.036
S8	11.72	0.020

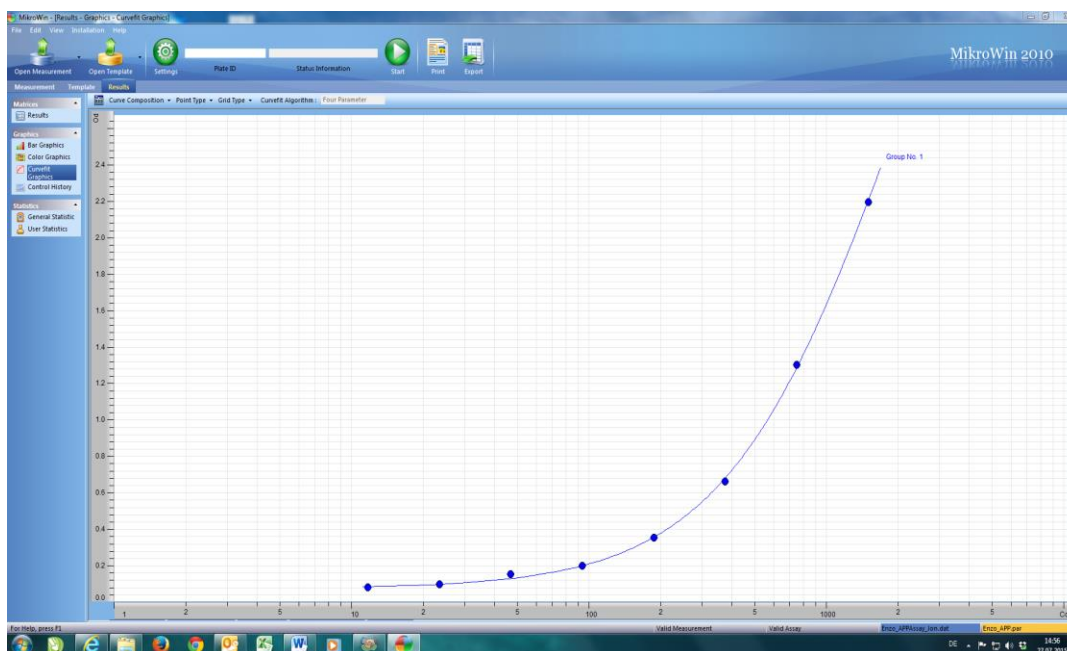


Fig. 1 Standard curve fitted with four parameter algorithm. Y-axis linear, X-axis logarithmic.

CONCLUSION

The standard curve showed excellent fitting so Crocodile provides a convenient and easy-to-use method for the automation of the Enzo APP Δ C31 ELISA kit. The assay procedure is extremely simple and involves only the addition of standards and controls while the instrument is processing all necessary dispense, wash, incubation and reading steps.

Changing the first steps in the assay procedure (pipetting antibody and standards/samples) doesn't affect the assay results at all.

#	Step Name	Description
1	Prime Antibody	Dispensing Volume 1000 μ l Inlet 1 Label "Antibody" Method: Priming Well Count: 1
2	Dispense Antibody	Dispensing Volume 50ul Inlet 1 Label "Antibody" Method: Standard Well Count: 20
3	Incubate 1	Shaking for 01:00:00 at Incubator with 2mm Amplitude at 10 Hz
4	Wash 1	Washing Method: Standard Wash Solution Inlet: 1 Wash buffer Cycles: 3 Volume: 300 μ l Delay: 2s Wait: 500ms Dispenser Depth: 1300 (Plate Offset 50) Aspiration Depth 2725 (Plate Offset 20) Sweep 5mm @1mm/s Well Count 96

5	Prime Conjugate	Dispensing Volume 1000 µl Inlet 2 Label "Conjugate" Method: Priming Well Count: 1
6	Dispense Conjugate	Dispensing Volume 100ul Inlet 2 Label "Conjugate" Method: Standard Well Count: 20
7	Incubate 2	Shaking for 00:30:00 at Incubator with 2mm Amplitude at 10 Hz
8	Wash 2	Washing Method: Standard Wash Solution Inlet: 1 Wash buffer Cycles: 3 Volume: 300 µl Delay: 2s Wait: 500ms Dispenser Depth: 1300 (Plate Offset 50) Aspiration Depth 2725 (Plate Offset 20) Sweep 5mm @1mm/s Well Count 96
9	Prime Substrate	Dispensing Volume 1000 µl Inlet 3 Label "Substrate" Method: Priming Well Count: 1
10	Dispense Substrate	Dispensing Volume 100ul Inlet 3 Label "Substrate" Method: Standard Well Count: 20
11	Incubate 3	Shaking for 00:30:00 at Incubator with 2mm Amplitude at 10 Hz
12	Prime Stop	Dispensing Volume 1000 µl Inlet 4 Label "Stop" Method: Priming Well Count: 1
13	Dispense Stop	Dispensing Volume 100ul Inlet 4 Label "Stop" Method: Standard Well Count: 20
14	Measure	Reading Single Wavelength Filter 1: 450nm (Pos:2) Well Count: 96

Table 3: Assay program with Crocodile Control Software