

DNA Fragmentation Assays

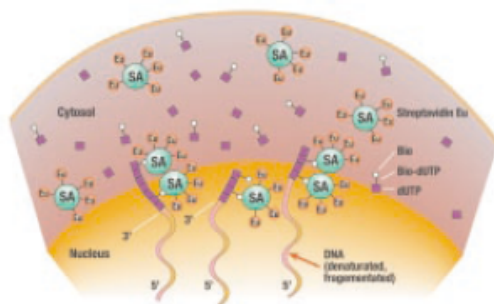
Application Note

A new cell-based DELFIA® DNA Fragmentation assay for testing apoptotic effect of lead compounds

INTRODUCTION

The DELFIA® DNA Fragmentation assay is performed in 96-well microplate format for quantitative detection of apoptosis. It is a simple and upscalable method to test substances for their apoptotic effect on cells.

During the process of apoptosis, DNA fragmentation occurs following the activation of endonucleases. The labeling of the 3' ends of DNA fragments provides an easy measure of cells undergoing apoptosis. Modified nucleotides are incorporated at the 3' ends by the activity of terminal deoxynucleotidyl transferase (TdT).



DNA fragmentation has long been used to distinguish apoptosis from necrosis, and is among the most reliable methods for detection of apoptotic cells. When DNA strands are cleaved or nicked by nucleases, 3'-hydroxyl ends are exposed. These hydroxyl ends can then serve as a starting point for terminal deoxynucleotidyl transferase (TdT), which adds deoxyribonucleotides in a template-independent fashion. Addition of labelled dUTP to the TdT reaction thus provides a means to label the DNA strand breaks. Once incorporated into DNA, the labelled nucleotide can be detected by Europium labeled streptavidin. This method of labeling DNA breaks is referred to as

terminal deoxynucleotide transferase dUTP nick-end labeling, or TUNEL¹.

METHODS

CHO cells, grown in RPMI medium supplemented with 10% FCS and 1% Penicillin & Streptomycin (Life Technologies), were seeded in 96-well plates and grown overnight. The cells were treated with 10 μ M staurosporine (Sigma) for 6 h at 37°C (CO₂ incubator). After treatment, the growth medium was removed and 100 μ L/well Fixation Solution was added. The cells were fixed 30 minutes at room temperature. The Fixation Solution was tapped off and the following reaction was started: 0.01% CHAPS, 5.5 U TdT enzyme (Amersham), 15 μ M dTTP, 5 μ M Bio-dUTP (Roche) and TdT buffer (Amersham, 500mM sodium cacodylate pH 7.2, 10 mM CoCl₂, 1 mM 2-mercapthoethanol), in a final reaction volume of 50 μ L. A control reaction without the enzyme was also started. The reaction was performed at 37°C for 30 minutes. The plate was washed 6 times using the DELFIA Platewash. Eu-labelled streptavidin was added to the wells at a concentration of 100 ng/mL in 50 μ L. The Eu-labelled streptavidin was diluted in DELFIA Assay Buffer. The plate was incubated 1 hour at room temperature and was washed 6 times prior to addition of 200 μ L/well Enhancement Solution. The plate was shaken for 5 minutes and the fluorescence was measured using the VICTOR³™ plate reader.

RESULTS

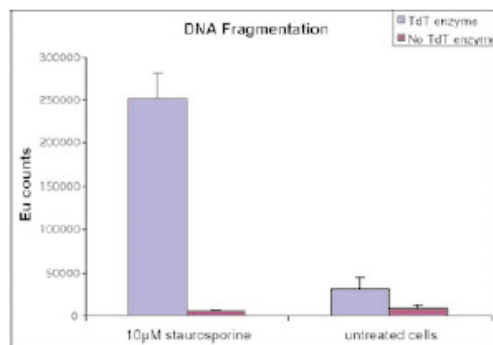


Fig.1. The results show the amount of DNA fragmentation in CHO cells treated 6h with 10µM staurosporin compared to untreated cells. The cells were fixed and the TUNEL reaction was performed at 37°C. The incorporated Bio-dUTP was detected using Eu-labelled streptavidin. The fluorescence was measured using the VICTOR²™ plate reader.

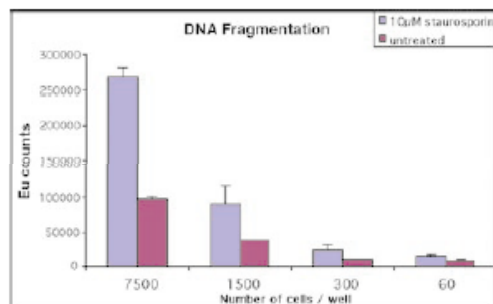


Fig.2. Detection of apoptosis in CHO cells. The amount of fragmented DNA is dependent on the number of cells. The cells were fixed and the TUNEL reaction was performed at 37°C. The incorporated Bio-dUTP was detected using Eu-labelled streptavidin. The fluorescence was measured using the VICTOR²™ plate reader.

The data demonstrate that the DELFIA® DNA Fragmentation Assay is a simple and fast microplate format method, for detection of DNA strand breaks in cells. This is a sensitive and efficient method for screening potential inducers of apoptosis.

REFERENCES:

- Gavrieli, Y., Sherman, Y. and Ben-Sasson S.A. (1992). Identification of programmed cell death in situ via specific labelling of nuclear DNA fragmentation. *J. Cell. Biol.* **119**, 493–501

Catalog Items: (see DELFIA buffers guide)

Cat.No.	Products	Size
1244-360	DELFLA Eu-labelled streptavidin	0.25 mg
1244-114	DELFLA Wash Concentrate	250 ml
1244-111	DELFLA Assay Buffer	250 ml
1244-105	DELFLA Enhancement Solution	250 ml
1450-516	Wallac Isoplates TC	25/100 plets
1450-517		
1296-026	DELFLA Platetwash	
1296-001/002	DELFLA Plateshake	
1296-003/004		
Wallac 1420	VICTOR ² Plate Reader	



Worldwide Headquarters: PerkinElmer Life Sciences, 549 Albany Street, Boston, MA 02118-2512 USA (800) 551-2121

European Headquarters: PerkinElmer Life Sciences, Imperiastraat 8, BE-1930 Zaventem Belgium +32 2 717 7911

Technical Support: in Europe: techsupport.europe@perkinelmer.com in US and Rest of World: techsupport@perkinelmer.com

Belgium: Tel: 0800 94 540 • **France:** Tel: 0800 90 77 62 • **Netherlands:** Tel: 0800 02 23 042 • **Germany:** Tel: 800 1 81 00 32 • **United Kingdom:** Tel: 0800 89 60 46
Switzerland: Tel: 0800 55 50 27 • **Italy:** Tel: 800 79 03 10 • **Sweden:** Tel: 020 79 07 35 • **Norway:** Tel: 800 11 947 • **Denmark:** Tel: 80 88 3477 • **Spain:** Tel: 900 973 255

All trademarks or registered trademarks are the property of PerkinElmer Life Sciences, Inc.

www.perkinelmer.com/lifesciences