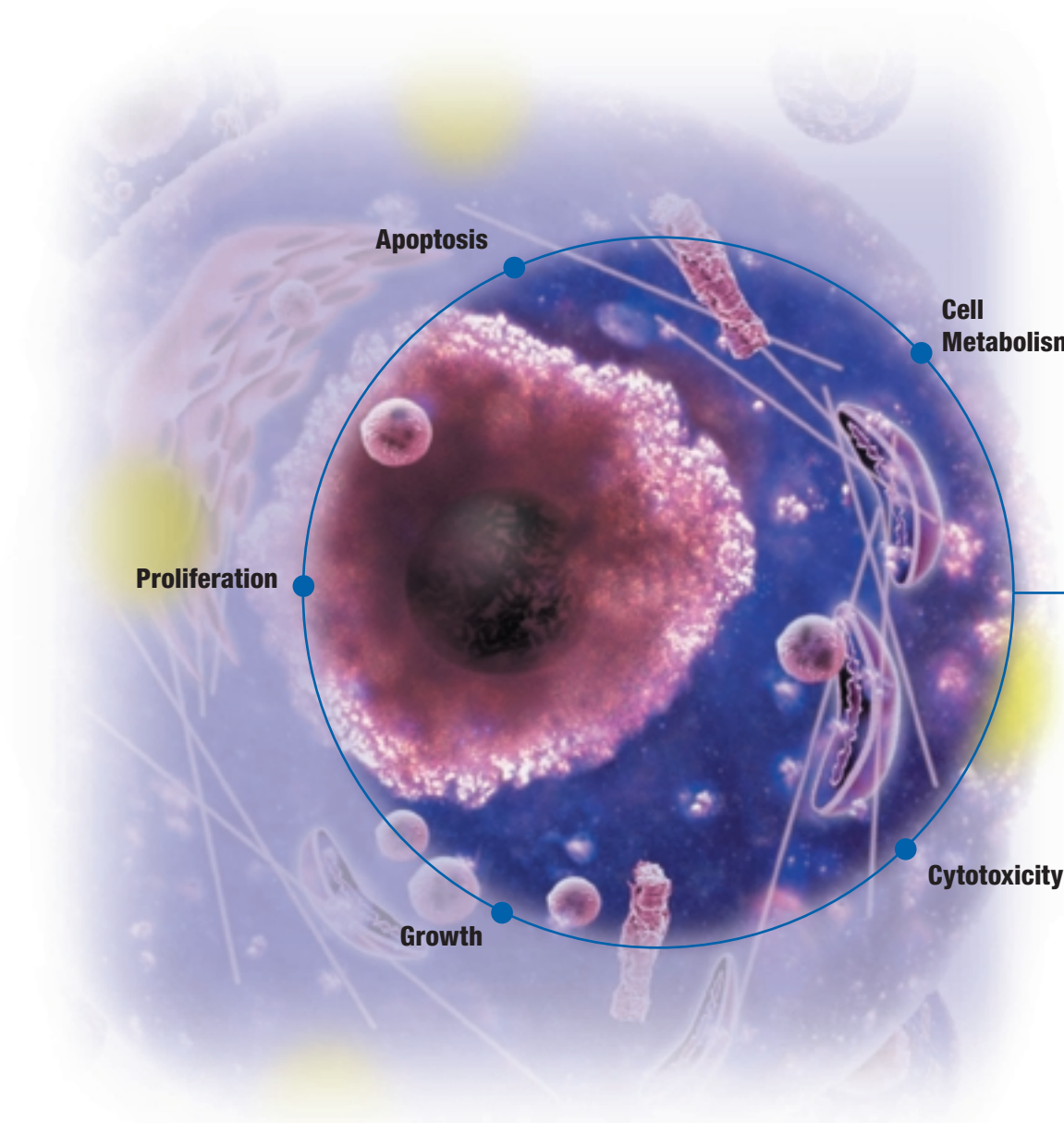


# Tox Suite

Non-radioactive  
assays for cell toxicity testing



# Technologies

**PerkinElmer's Tox Suite portfolio offers** a wide array of sensitive, non-radioactive assays to study different aspects of cell growth and death. ATP<sup>lite</sup>™ assays allow you to assess cell vitality and the level of cell metabolism. DELFIA® cell proliferation assays can be used to enumerate cells by quantitating DNA synthesis of cells. Early events of apoptosis such as DNA fragmentation can be studied using a DELFIA assay or appearance of caspases with TruPoint™ technology.

These assays combined with the fast EnVision™ multilabel plate reader allow hits from HTS screens to be conveniently and efficiently evaluated. Screening compounds for their toxicological properties early in the drug discovery process is vital to lead optimization.

## Tox Suite: The complete portfolio of platforms for all your cell toxicity assays

### Cell Metabolism



Cellular damage will result in loss of ability to maintain and provide energy for metabolic cell function and growth. Luminescent ATP<sup>lite</sup> is a highly sensitive way to measure cellular ATP levels. ATP<sup>lite</sup> allows you to put together the perfect assay in less than 10 minutes! Since ATP<sup>lite</sup> is a homogenous assay with one plate preparation (from seeding cells to detection), it offers significant time and labor savings.

#### Products

ATP<sup>lite</sup>, 300 points  
6016943

ATP<sup>lite</sup>, 1000 points  
6016941

### Cell Proliferation

#### DELFIA Cell Proliferation Kit

Cell proliferation assays encompass a characterization of compounds for their capability to either promote or retard cell growth. The DELFIA Cell Proliferation kit is based on quantitating changes in the level of DNA synthesis in cell populations to reflect the number of proliferated cells. During cell growth, 5-bromo-2'-deoxyuridine (BrdU) is incorporated into newly synthesized DNA. It is then immunochemically detected with the extraordinarily sensitive DELFIA assay. There is no need for radiochemicals, dyes or microscopes.

#### Products

DELFIA Proliferation Kit,  
960 wells  
AD0200

# you trust



## Cell Mediated Cytotoxicity

### DELFIA Cytotoxicity Kit

Immunotoxicological effects of hits can be studied with the DELFIA Cytotoxicity Kit. The non-radioactive DELFIA method mimics the principle of the well-established  $^{51}\text{Cr}$ -release test. Compared to radiometric assays, release of the fluorescence enhancing ligand is fast, typically shortening assay times.

#### Products

EuTDA Cytotoxicity Kit,  
contains 10 plates  
AD0116

DELFIA BATDA  
labeling reagent, 50  $\mu\text{L}$   
C136-100

Europium solution, 200 mL  
C135-100

Lysis buffer, 30 mL  
4005-0010

## Apoptosis

### TruPoint Caspase Reagents

#### DELFIA Eu-labeled Streptavidin

TruPoint Caspase-3 reagents can be used for measurement of caspase activity from apoptotic cell lysates (fresh or frozen). The assay is homogenous, sensitive and robust and applicable to HTS screens. TruPoint Assay Kits for caspase-6 and caspase-8 are also available. For other TruPoint substrates, please email at [Labellingservices@perkinelmer.com](mailto:Labellingservices@perkinelmer.com).

Various sensitive, simple DELFIA assays using Europium (Eu)-labeled streptavidin have been used to study different apoptotic pathways. Complicated DNA fragment analysis can be simplified indirectly with Eu-labeled streptavidin (PerkinElmer application note #1234-9864).

#### Products

TruPoint Caspase-3 Assay Kit,  
2 plates  
AD0125

TruPoint Caspase-3 substrate,  
5 vials  
CR93-100

TruPoint Caspase reaction  
buffer, 1000 mL  
CR94-100

Lysis buffer, 30 mL  
4005-0010

DELFIA Eu-labelled  
streptavidin, 0.25 mg  
1244-360

**PerkinElmer Life and  
Analytical Sciences**  
710 Bridgeport Avenue  
Shelton, CT 06484-4794 USA  
Phone: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



---

For a complete listing of our global offices, visit [www.perkinelmer.com/lasoffices](http://www.perkinelmer.com/lasoffices)

©2003 PerkinElmer, Inc. All rights reserved. PerkinElmer is a registered trademark of PerkinElmer, Inc. All trademarks or registered trademarks are the property of PerkinElmer, Inc. PerkinElmer reserves the right to change this document at any time and disclaims liability for editorial, pictorial or typographical errors.