

ApoSENSOR™ Cell viability assay kit

A simple no-wash, chemiluminescent cell viability assay

The assay utilizes bioluminescent detection of the ATP level for a rapid screening of apoptosis and cell proliferation simultaneously in mammalian cells. The assay utilizes the enzyme luciferase to catalyze the formation of light from ATP and luciferin. The light can be measured using a luminometer or Beta Counter. The assay can be fully automatic for high throughput (10 seconds/sample) and is highly sensitive (detects 10-100 mammalian cells/well).

Ordering Information

[Order Online »](#)

| | |
|------------------|-----------|
| ALX-850-247-KI01 | 200 tests |
|------------------|-----------|

Manuals, SDS & CofA

[View Online »](#)

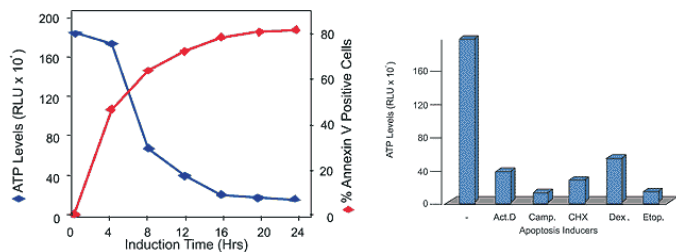


Figure: ApoSENSOR™ Cell Viability Assay effectively detects apoptosis. Apoptosis was induced in Jurkat cells and ATP levels were analyzed according to the kit instructions. **Left:** Time course showing decrease in ATP levels following induction of apoptosis with camptothecin. The decrease in ATP level correlates with the increase in annexin V positive staining cells detected by Annexin V-FITC Apoptosis Detection Kit according to the kit instructions. **Right:** Jurkat cells were treated with various apoptosis inducers as indicated for 12 hours and ATP levels decrease dramatically for all treatments.

Handling & Storage

Handling Protect from light.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Chemiluminescence

Contents Nucleotide Releasing Buffer, ATP Monitoring Enzyme, Enzyme Reconstitution Buffer, ATP.

Technical Info / Product Notes **Application Note:**
[Automated Monitoring of Protein Expression and Metastatic Cell Migration Using Magnetic 3D Bioprinted Colorectal Cancer Cells](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)