

Annexin V-EGFP

apoptosis detection

kit

EGFP (enhanced green fluorescent protein) has emerged as an ideal fluorochrome for FACS analysis and fluorescent microscopy as it offers a more intense signal than wild-type GFP. It has Ex(max): 488nm and Em(max): 530nm. The green signal produced makes EGFP ideal for dual staining with propidium iodide (PI).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-253-KI02	100 tests
------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

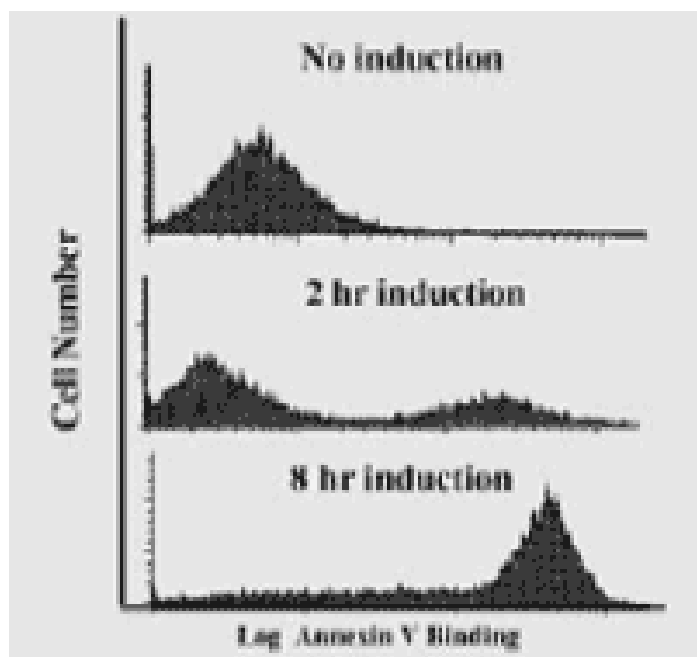


Figure: Detection of apoptosis using Annexin V-EGFP Apoptosis Detection Kit (Prod. No. ALX-850-253).

Method: Apoptosis was induced in Jurkat cells by camptothecin ($2\mu\text{M}$) for various times as indicated. Cells were collected and incubated with Annexin V-EGFP for 5 min. acc. to kit instructions. Results were analyzed by flow cytometry.

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, Fluorescence microscopy

Application Notes For the detection of apoptotic cells by flow cytometry.

Contents Annexin V-EGFP, propidium iodide solution and binding buffer.
An assay protocol is included.



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)