

Annexin V-FITC

apoptosis detection

kit

Easy, quick assay to distinguish apoptotic cells vs necrotic cells

FITC-labelled recombinant human annexin V shows bright green fluorescence (Ex(max): 488nm, Em(max): 530nm).

Citations: 22

[View Online »](#)

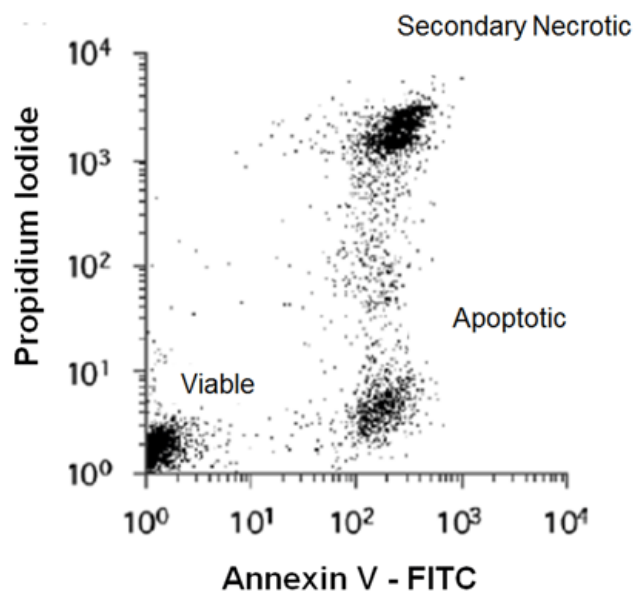
Ordering Information

[Order Online »](#)

ALX-850-0200	300 tests
--------------	-----------

Manuals, SDS & CofA

[View Online »](#)



Typical flow cytometry results. The Annexin V-FITC apoptosis detection kit allows the detection and discrimination between apoptotic and necrotic cells.

Handling & Storage

Handling	After opening annexin V-FITC, aliquot contents and freeze at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Flow Cytometry, Fluorescence microscopy, Fluorescent detection

Application Notes For the detection of apoptotic cells by flow cytometry or microscopy

Contents

Annexin V-FITC: 100 tests/500µl; 300 tests/1.5ml; Liquid. In 50mM TRIS, pH 7.4, 100mM sodium chloride, 1% BSA and 0.02% sodium azide.

Binding Buffer: 50ml (4x concentrated) Concentration after dilution: 10mM HEPES/NaOH, pH 7.4, 140mM sodium chloride, 2.5mM CaCl₂.

Propidium Iodide: 20µg/ml (100 tests: 1.6ml, 300 tests: 2 x 1.6ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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