

FLICA® 660 Caspase 3/7 Assay Kit

Detect caspase 3 and 7 activity with the FLICA 660 Caspase 3/7 Assay Kit. This in vitro assay employs the fluorescent inhibitor probe 660-DEVD-FMK to label active caspase 3 and 7 enzyme in living cells. Analyze the fluorescent signal using fluorescence microscopy or flow cytometry.

Ordering Information

[Order Online »](#)

ICT-9125	25 tests
----------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, Fluorescence microscopy

Contents FLICA Caspase-3/7 Inhibitor Reagent (660-DEVD-FMK)
10X Apoptosis Wash Buffer
Fixative

Technical Info / Product Notes For the Manual please [click here](#).



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