

Pyroptosis/Caspase 1 Assay, Green (25 – 50 tests)

ICT's Pyroptosis/Caspase-1 Assay Kit utilizes our popular FLICA technology to detect caspase-1 activation. Using this kit, researchers can easily assess pyroptotic cells and utilize nigericin as a positive control. Analyze the green fluorescent signal using fluorescence microscopy, a fluorescent plate reader, or by flow cytometry.

Ordering Information

[Order Online »](#)

ICT-9145	25 tests
----------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, Fluorescence microscopy, Fluorescent detection

Contents FLICA Caspase-1 Inhibitor Reagent (FAM-YVAD-FMK), 10X Cellular Wash Buffer, Fixative, Hoechst 33342, Nigericin



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)