

Live vs. Dead Cell Differentiation Staining Kit

To stain live cells. Stained cells are visualized by fluorescence microscopy.

Differentiating between live and dead cells is common but a very important task when investigating growth regulation controls, cell death mechanisms and pathways. AkrivisBio's Live vs. Dead Cell Differentiation Staining Kit provides ready-to-use reagents. The kit combines a proprietary cell-permeable green fluorescent dye to stain live cells and propidium iodide, a red non-permeable dye, to stain all cells. Stained cells are visualized by fluorescence microscopy.

Ordering Information

[Order Online »](#)

AKR-COS-0103	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescence microscopy

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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

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

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

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

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

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