

Crystal Violet Cell Cytotoxicity Assay

To study cell proliferation, viability or cytotoxicity.

Crystal violet binding as a measure of cytotoxicity is a well established method for detection of cell viability or drug cytotoxicity. Crystal violet is a triarylmethane dye that binds to ribose containing molecules such as DNA. Typically, dead cells detach from cell culture plates and are removed during the washing steps. Using crystal violet to quantify DNA of the remaining adherent cell population allows a quick measure of cell viability by monitoring its absorbance at 570 nm. AkrivisBio's Crystal Violet Cell Cytotoxicity Assay is a simple, sensitive assay to study cell proliferation, viability or cytotoxicity. This assay provides a quantitative measurement of cell survival and can be used to evaluate the cytotoxic effects of various treatments or compounds.

Ordering Information

[Order Online »](#)

AKR-CPT-0102	1000 wells
--------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

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-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