

# MTT Cell Proliferation Assay

For the quantification of viable cells in proliferation and cytotoxicity assays.

The ability to quantify the number of viable cells in a system under controlled conditions is a requirement in cell biology research. There are several such assays designed to measure metabolic activity of cultured cells, of which MTT is one. The advantages of AkrivisBio's MTT Cell Proliferation Assay are sensitivity (subtle changes can be observed), cost-effectiveness, versatility (adherent or suspension cultures, 96 or 384 well plates) and the ability to monitor metabolic activity over long periods of time. The assay is based on the conversion of MTT to an insoluble formazan. Cells with active metabolism convert MTT while dead cells can't. Absorbance at 590 nm is proportional to metabolic activity.

## Ordering Information

[Order Online »](#)

AKR-CPT-0103	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](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)