

MTS Cell Proliferation Assay

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

Among the various cell proliferation assays, there are distinct advantages to using MTS as the detection reagent. One major plus with MTS is that it is a direct measurement of mitochondrial metabolic activity, thus providing a more accurate assessment of cell viability and growth, compared to MTT. AkrivisBio's MTS Cell Proliferation Assay is a one step, optimized reagent for quantification of viable cells in proliferation and cytotoxicity assays. The reduction of MTS by viable cells to a colored formazan is believed to be due to mitochondrial dehydrogenases. The formazan is simply quantified by absorbance at 490 nm.

Ordering Information

[Order Online »](#)

AKR-CPT-0104-2500	2500 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)