Glowing products for science

## MTT Cell Viability Assay Kit

A simple, colorimetric assay for determining live cell numbers by absorbance on a microplate reader.



## Product attributes

Apoptosis/viability marker	Metabolic activity
For live or fixed cells	Cell lysis required
Assay type/options	Endpoint assay, Homogeneous assay, No-wash staining, Real-time imaging
Detection method/readout	Microplate reader (absorbance)
Storage Conditions	Store at -10 to -35 °C

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

The MTT Cell Viability Assay Kit provides a simple method for determining live cell numbers by absorbance on a microplate reader. MTT is a tetrazolium salt that is turned into a purple formazan product after reduction by mitochondrial enzymes that are only present in metabolically active live cells. The amount of formazan product generated is proportional to the number of living cells in the sample.

- Measure cell metabolism by reading absorbance at 570 nm
- Non-radioactive, simple procedure
- Endpoint assay, requires cell lysis

Determination of live cell numbers is often used to assess the rate of cell proliferation and cytotoxicity caused by drugs and cytotoxic agents. Among all non-radioactive viability assays, MTT assay developed by Mossman is one of the most versatile and popular assays. At the end of the assay, the cells containing the formazan product are solubilized and then photometrically quantified at 570 nm.

Also see the XTT Cell Viability Assay Kit, which does not require solubilization before absorbance measurement.

## References

- 1. PLoS ONE (2011) 6(8): e22635. doi:10.1371/journal.pone.0022635 2. Cell (2011) 145, 941-955. doi:10.1016/j.cell.2011.05.004 3. PLoS ONE (2012) 7(4): e35866. doi:10.1371/journal.pone.0035866 4. J. Virol (2013) Jul;87(14):8179-94. doi:10.1128/JVI.00443-13 5. J Am Chem Soc. (2014) Apr 9;136(14):5295-300. doi: 10.1021/ja408505n.
- 6. Endocrinology. (2015) Dec;156(12):4695-706. doi: 10.1210/en.2015-1500
- 7. J Mol Neurosci. (2015) May;56(1):122-30. doi: 10.1007/s12031-014-0473-6
- 8. Nat Commun (2016) Feb 16;7:10686. doi: 10.1038/ncomms10686

This datasheet was generated on November 6, 2025 at 07:27:41 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mtt-cell-viability-assay-kit/">https://biotium.com/product/mtt-cell-viability-assay-kit/</a>