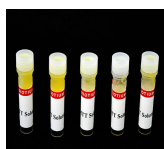


## MTT Cell Viability Assay Kit

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



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

This datasheet was generated on June 5, 2026 at 04:45:25 AM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/mtt-cell-viability-assay-kit/>

### Product attributes

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