XTT 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	For live/intact cells
Assay type/options	Endpoint assay, High-throughput assay, Homogeneous assay, Real-time imaging
Detection method/readout	Microplate reader (absorbance)
Storage Conditions	Store at 2 to 8 °C

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

Product Description

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

- Measure cell metabolism by reading absorbance at 475 nm
- Non-radioactive, simple procedure
- XTT is a tetrazolium derivative
- Endpoint assay
- Unlike MTT, does not require solubilization step

Determination of live cell numbers is often used to assess the rate of cell proliferation and cytotoxicity caused by drugs and cytotoxic agents. The non-radioactive MTT assay developed by Mossman is one of the most popular viability assays; however, the product produced is non-soluble, and requires a solubilization step. Unlike MTT, XTT is reduced to a highly water-soluble orange-colored product instead of the insoluble formazan from MTT, thus eliminating the solubilization step required for the MTT assay.

The amount of water-soluble product generated from XTT reduction is proportional to the number of living cells in the sample and can be photometrically quantified at 475 nm. Continuous color development permits measurements at multiple time points for extended detection range.

References

- 1. PLoS ONE (2013) 8(8): e71165 doi:10.1371/journal.pone.0071165 2. ACS Nano (2013) 7, 9, 7524-7533 doi.org/10.1021/nn403722x 3. Oncogene (2014) 6 October doi:10.1038/onc.2014.328
- 5. Chicogene (2014) Sep 26;289(39):26882-94. doi: 10.1074/jbc.M114.567685 5. Environ. Sci Nano (2015) 2, 370-379, doi: 10.1039/C5EN00031A 6. Lab Invest (2015) Mar;95(3):296-307 doi: 10.1038/labinvest.2014.161

- 7. Mol Biol Rep (2016) 43, 229-240 doi:10.1007/s11033-016-3948-4
- 8. Nat Commun (2016) 7, 11551 doi:10.1038/ncomms11551

This datasheet was generated on November 18, 2025 at 01:50:46 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/xtt-cell-viability-assay-kit/