

MTT

MTT (3-(4,5-Dimethylthiazol-2-yl)-2,5-diphenyltetrazolium bromide) is a tetrazolium salt that is commonly used to detect reductive metabolism in cells for viability, proliferation and cytotoxicity assays. On reduction tetrazolium salts turn into deeply colored products that can be easily quantified by colorimetric measurement.



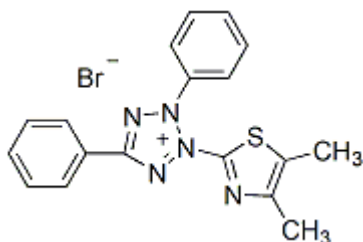
Product attributes

CAS number	298-93-1
------------	----------

Product Description

MTT (3-(4,5-Dimethylthiazol-2-yl)-2,5-diphenyltetrazolium bromide) is a tetrazolium salt that is commonly used to detect reductive metabolism in cells for viability, proliferation and cytotoxicity assays. On reduction tetrazolium salts turn into deeply colored products that can be easily quantified by colorimetric measurement. MTT forms a purple formazan product that needs to be solubilized prior to quantification. MTT and XTT ([10060](#)) are particularly useful for HTS of antiviral and anticancer compounds and for the studies of cytokine effect on cell proliferation. Please also see the [MTT Cell Proliferation Assay Kit](#).

- Yellow solid soluble in water or DMSO
- C₁₈H₁₆BrN₅S
- MW: 414.32
- [298-93-1]



References

1. Biomaterials (2010) doi:10.1016/j.biomaterials.2010.08.035
2. Parasitology 107, 175 (1993).
3. Antimicro Agents Chemother. 36, 1619 (1992).
4. J Immunol Meth 147, 153 (1992).
5. Cancer Res 48, 4827 (1988).

This datasheet was generated on June 12, 2026 at 12:10:29 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mtt-3-45-dimethylthiazol-2-yl-25-diphenyltetrazolium-bromide/>