

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: MTT

CATALOG #: 10059

[298-93-1]

Br N—N
N
S
CH
3

PROPERTIES:

Color & Form Yellow solid **Purity** $\geq 98\%$

Soluble in water (upto 10mg/ml)

STORAGE AND HANDLING:

Store desiccated at 4 °C and protect from light.

APPLICATION: MTT 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, see below) are particularly useful for HTS of antiviral and anticancer compounds and for the studies of cytokine effect on cell

proliferation

Ref: 1) Prog Histochem Cytochem. 9(3), 1-56 (1976); 2) Histochemistry, 76(2), 381-405(1982); 3) Antimicro Agents Chemother. 36, 1619(1992); 4) Cancer Res. 48, 4827(1988); 5) Parasitology 107, 175(1993); 6) J. Immunol. Meth. 147,

153(1992)

TOXICITY: Unknown.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on

clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.