XTT

(2,3-Bis-(2-Methoxy-4-nitro-5-sulfophenyl)-2H-tetrazolium-5-carboxanilidisodium salt) is a tetrazolium derivative used widely in cell viability and proliferation testing. The material is readily reduced in viable cells to a highly water-soluble orange colored product (λ = 475 nm).

a sictum

Product attributes

CAS number

111072-31-2

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

Product Description

XTT (2,3-Bis-(2-Methoxy-4-nitro-5-sulfophenyl)-2H-tetrazolium-5-carboxanilide, disodium salt) is a tetrazolium derivative used widely in cell viability and proliferation testing. The material is readily reduced in viable cells to a highly water-soluble orange colored product (λ = 475 nm). Applications of the reagent include studies of antifungal susceptibility, drug sensitivity of cells, parasitic nematode viability and tumor cell cytotoxicity. Please also see our XTT Cell Proliferation Assay Kit.

- Pale yellow solid soluble in water
- Store desiccated at 4°C
- C₂₂H₁₂N₇NaO₁₃S₂
- MW: 675
- [111072-31-2]

References

- 1. European Journal of Pharmacology (2010) doi:10.1016/j.ejphar.2010.06.029
- 2. Toxicology in Vitro (2009) doi:10.1016/j.tiv.2009.07.01
- 3. Parasitology 107, 175 (1993).
- 4. J Immunol Meth 147, 153 (1992).
- 5. Antimicrob Agents Chemother 36, 1619 (1992).
- 6. Cancer Res 48, 4827 (1988).

This datasheet was generated on November 4, 2025 at 07:17:11 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/xtt-23-bis-2-methoxy-4-nitro-5-sulfophenyl-2h-tetrazolium-5-carboxanilide-disodium-salt/