

## XTT

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 ( $\lambda = 475 \text{ nm}$ ).



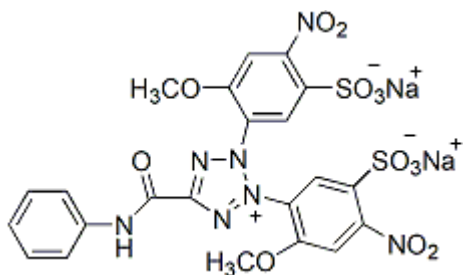
### Product attributes

|            |             |
|------------|-------------|
| CAS number | 111072-31-2 |
|------------|-------------|

## 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 ( $\lambda = 475 \text{ nm}$ ). 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  $4^{\circ}\text{C}$
- $\text{C}_{22}\text{H}_{12}\text{N}_7\text{NaO}_{13}\text{S}_2$
- 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 January 2, 2026 at 01:41:47 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/>