

NBT

NBT (Nitro blue tetrazolium, chloride) is an oxidant and is widely used in conjunction with BCIP or BCIP derivatives to detect phosphatase activity. NBT is also a substrate for dehydrogenate and other oxidizes.



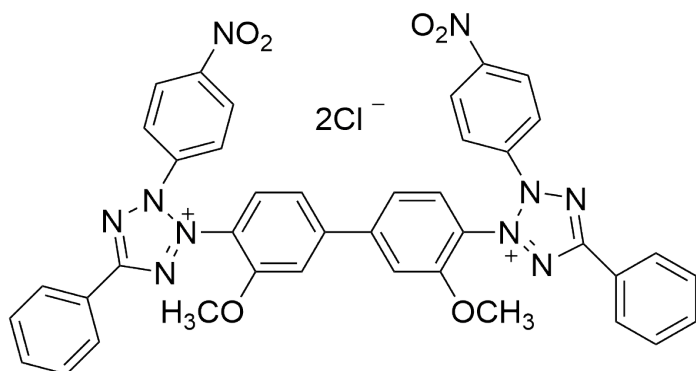
Product attributes

CAS number	298-83-9
------------	----------

Product Description

NBT (Nitro blue tetrazolium, chloride) is an oxidant and is widely used in conjunction with BCIP or BCIP derivatives to detect phosphatase activity. NBT is also a substrate for dehydrogenate and other oxidizes. Thus, it is often used to assess cell viability, proliferation and cytotoxicity, and is also frequently used in histochemistry studies. The colorless NBT is soluble in water, but forms a dark blue-colored, highly insoluble precipitate after being enzymatically reduced.

- White solid soluble in water and DMSO
- Store at 4°C
- C₄₀H₃₀Cl₂N₁₀O₆
- MW: 817.7
- [298-83-9]



References

1. Meth Enzymol 6, 958 (1963).
2. Anal Biochem 56, 353 (1973).
3. J Immunol Meth 179, 95 (1995).
4. J Infect. Dis. 172, 1153 (1995).
5. J Appl Bacteriol 74, 433 (1993).
6. J Immunol Meth 65, 55 (1983).
7. Curr Biol. 27, 2869 (2017), DOI: [10.1016/j.cub.2017.08.019](https://doi.org/10.1016/j.cub.2017.08.019)
8. J Photochem Photobiol B. 106, 24 (2012), DOI: [10.1016/j.jphotobiol.2011.09.006](https://doi.org/10.1016/j.jphotobiol.2011.09.006)
9. J Nanosci Nanotechnol. 5, 3285 (2014) DOI: [10.1166/jnn.2014.8719](https://doi.org/10.1166/jnn.2014.8719)
10. J Photochem Photobiol B. 98, 144 (2010) DOI: [10.1016/j.jphotobiol.2009.11.015](https://doi.org/10.1016/j.jphotobiol.2009.11.015)

This datasheet was generated on June 4, 2026 at 10:42:45 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nbt-nitro-blue-tetrazolium-chloride/>