Glowing products for science

BCIP/NBT Kit

BCIP/NBT are widely used together as a chromogenic phosphatase substrate for the detection of alkaline phosphatase labeled proteins in a variety of applications, such as immunohistochemistry, westerns, and in situ hybridization. This kit contains 100 mg each of BCIP sodium salt (10001) and the oxidant NBT (Nitro blue tetrazolium, 10008).

Product attributes

CAS number 102185-33-1, 298-83-9

Storage Conditions Store at 2 to 8 °C or below. Protect from light, Desiccate

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

Product Description

BCIP (5-Bromo-4-chloro-3-indoxyl phosphate) is the most widely used chromogenic phosphatase substrate that is often used with the oxidant NBT (nitro blue tetrazolium chloride, 10008), which facilitates the precipitation. NBT/BCIP produces a dark blue/violet precipitate product in the presence of alkaline phosphatase activity and alkaline phosphatase labeled proteins in a variety of applications, such as immunohistochemistry, westerns, and in situ hybridization. The stain can be mounted with permanent or aqueous mounting medium.

- Dark blue colormetric detection of alkaline phosphatase activity and labels
- Compatible with a variety of applications
- Use BCIP alone or in combination with NBT
- Can be mounted with permanent or aqueous mounting medium

This kit contains 100 mg each of BCIP sodium salt (10001) and the oxidant NBT (10008). We also offer Alkaline Phosphatase Conjugated Antibodies.

Find the Right Stain for your Application

The original BCIP forms a dark blue (λ_{max} 615 nm) precipitate and is available in two different salt formulations; BCIP, toluidine salt is soluble in DMF while BCIP, sodium salt is soluble in water. We also offer a Pink BCIP derivative, which produces a pink colored (λ_{max} 540 nm) precipitate. BCIP Red produces a red colored (λ_{max} 565 nm) precipitate. Please see our BCIP Kits that are paired with NBT (nitro blue tetrazolium chloride) for user convenience.

References

- Histochemistry 58, 203 (1978), <u>DOI: 10.1007/bf00495720</u>
 Biotechniques 12, 656 (1992), <u>PMID: 1381193</u>
 J Photochem Photobiol B. 98, 144 (2010) <u>DOI: 10.1016/j.jphotobiol.2009.11.015</u>
- Dev. Dyn., 240, 589 (2011), DOI: 10.1002/dvdy.22544
 J Photochem Photobiol B. 106, 24 (2012), DOI: 10.1016/j.jphotobiol.2011.09.006
- J Nanosci Nanotechnol. 5, 3285 (2014) DOI: 10.1166/jnn.2014.8719
- 7. Dev Comp Immunol. 65, 41 (2016), <u>DOI: 10.1016/j.dci.2016.06.017</u> 8. Molecular Medicine Reports 15, 1455 (2017), <u>DOI: 10.3892/mmr.2017.6162</u>
- 9. Curr Biol. 27, 2869 (2017), DOI: 10.1016/j.cub.2017.08.019

This datasheet was generated on November 15, 2025 at 04:20:28 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bcipnbt-kit/