BCIP, Toluidine

The most widely used 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 form is soluble in DMF.



Product attributes

CAS number	6578-06-9
Molecular weight	433.6
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, which forms a dark blue (λ_{max} 615 nm) precipitate on enzymatic hydrolysis. It is often used with the oxidant NBT (nitro blue tetrazolium chloride), which facilitates the precipitation, to detect alkaline phosphatase-activity and -labeled proteins in a variety of applications, such as immunohistochemistry, westerns, and in situ hybridization.

- Blue/Purple colormetric detection of alkaline phosphatase activity and labels
- Compatible with a variety of applications
- Use alone or in combination with NBT
- White solid soluble in DMF

For your convenience, we offer a BCIP/NBT Kit (10003) that contains both reagents. 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.

Molecular Structure:

References

- 1. Histochemistry 58, 203 (1978), DOI: 10.1007/bf00495720

- Histocheritary 36, 203 (1976), <u>DOI: 10.1007/bit0493720</u>
 Biotechniques 12, 656 (1992), <u>PMID: 1381193</u>
 Dev. Dyn., 240, 589 (2011), <u>DOI: 10.1002/dvdy.22544</u>
 Dev Comp Immunol. 65, 41 (2016), <u>DOI: 10.1016/j.dci.2016.06.017</u>
 Molecular Medicine Reports 15, 1455 (2017), <u>DOI: 10.3892/mmr.2017.6162</u>

This datasheet was generated on November 10, 2025 at 03:27:47 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bciptoluidine-5-bromo-4-chloro-3-indoxyl-phosphate-p-toluidine-salt/