

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 of antigen	433.6
Storage Conditions	Store at 2 to 8 °C or below, Protect from light, Desiccate

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.

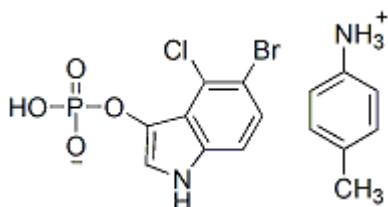
- Blue/Purple colorimetric detection of alkaline phosphatase activity and labels
- Compatible with a variety of applications
- Use alone or in combination with NBT
- White solid soluble in DMF

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](https://doi.org/10.1007/bf00495720)
2. Biotechniques 12, 656 (1992), [PMID: 1381193](https://pubmed.ncbi.nlm.nih.gov/1381193/)
3. Dev. Dyn., 240, 589 (2011), [DOI: 10.1002/dvdy.22544](https://doi.org/10.1002/dvdy.22544)
4. Dev Comp Immunol. 65, 41 (2016), [DOI: 10.1016/j.dci.2016.06.017](https://doi.org/10.1016/j.dci.2016.06.017)
5. Molecular Medicine Reports 15, 1455 (2017), [DOI: 10.3892/mmr.2017.6162](https://doi.org/10.3892/mmr.2017.6162)

This datasheet was generated on June 11, 2026 at 08:33:55 PM. 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/>