

## 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

## 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 colorimetric 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 ( $\lambda_{\max}$  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 ( $\lambda_{\max}$  540 nm) precipitate. [BCIP Red](#) produces a red colored ( $\lambda_{\max}$  565 nm) precipitate. Please see our [BCIP Kits](#) that are paired with [NBT \(nitro blue tetrazolium chloride\)](#) for user convenience.

## References

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

This datasheet was generated on January 14, 2026 at 06:20:23 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/bcipnbt-kit/>