Glowing products for science

BCIP Pink/NBT Kit

NBT and BCIP PINK 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 Pink (10006) and the precipitation inducer NBT (10008).

siètum 8

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

Product attributes

CAS number 159954-33-3, 298-83-9

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

Product Description

BCIP Pink (6-Chloro-3-indoxyl phosphate, p-toluidine salt) is a derivative of BCIP. While BCIP yields a blue precipitating product, BCIP Pink produces a pink colored (λ_{max} 540 nm) precipitate. BCIP and the derivatives are the most widely used chromogenic phosphatase substrate. They are often used with the oxidant NBT (nitro blue tetrazolium chloride), which facilitates the precipitation, to detect alkaline phosphatase activity and alkaline phosphatase labeled proteins in a variety of applications, such as immunohistochemistry, westerns, and *in situ* hybridization.

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

This kit contains 100 mg each of <u>BCIP Pink (10006)</u> and the precipitation inducer <u>NBT (10008)</u>. We also offer <u>Alkaline Phosphatase Conjugated Antibodies</u>.

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.

This datasheet was generated on November 15, 2025 at 03:17:37 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bcip-pinknbt-kit/