

BCIP Pink

A pink derivative of BCIP that produces a pink colored (λ_{\max} 540 nm) precipitate instead of the dark blue precipitate of BCIP. BCIP and derivatives are the most widely used chromogenic phosphatase substrate for the detection of alkaline phosphatase labeled proteins and are compatible with a variety of applications.



Product attributes

CAS number	159954-33-3
Molecular weight of antigen	354.7
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.

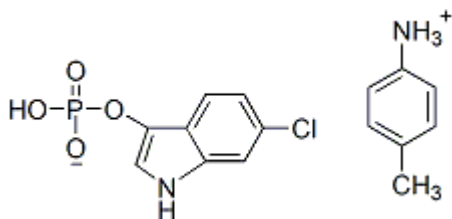
- Pink colorimetric detection of alkaline phosphatase activity and labels
- Compatible with a variety of applications
- Use alone or in combination with NBT
- White solid soluble in DMF

For your convenience, we offer a [BCIP Pink/NBT kit \(10007\)](#) 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.

Chemical Structure:



This datasheet was generated on June 13, 2026 at 08:51:41 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/bcip-pink-6-chloro-3-indoxyl-phosphate-p-toluidine-salt/>