

## 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 \(10008\)](#).



### 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 ( $\lambda_{\max}$  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 BCIP Pink alone or in combination with NBT
- White solid soluble in DMF

This kit contains 100 mg each of [BCIP Pink \(10006\)](#) and the precipitation inducer [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.