

BCIP Red/NBT Kit

A red derivative of BCIP that produces a red (λ_{\max} 565 nm) colored 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 in a variety of applications.



Product attributes

CAS number	298-83-9, 6769-80-8
Storage Conditions	Store at 2 to 8 °C or below, Protect from light, Desiccate

Product Description

BCIP Red (5-Bromo-6-chloro-3-indoxyl phosphate, *p*-toluidine salt) is a derivative of BCIP. While BCIP yields a dark blue precipitating product, BCIP Red produces a red (λ_{\max} 565 nm) colored 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.

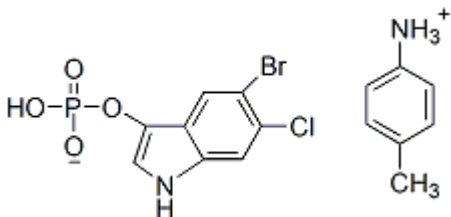
- Red colorimetric detection of alkaline phosphatase activity and labels
- Compatible with a variety of applications
- Use BCIP Red alone or in combination with NBT
- White solid soluble in DMF

This kit contains 100 mg each of [BCIP Red \(10004\)](#) 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 (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 BCIP Red:



References

1. Dev Dyn. 236, 1475 (2007) [DOI: 10.1002/dvdy.21152](https://doi.org/10.1002/dvdy.21152)
2. Dev Dyn. 238, 2179 (2009) [DOI: 10.1002/dvdy.21883](https://doi.org/10.1002/dvdy.21883)