

BCIP Red

A red derivative of BCIP that produces a red (λ_{\max} 565 nm) colored precipitate instead of the blue/violet 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	6769-80-8
Molecular weight	433.6
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 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.

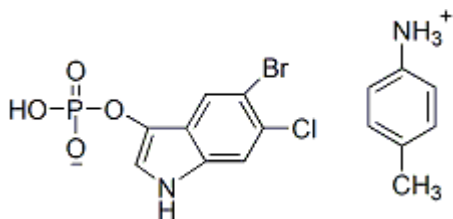
- Red colorimetric detection of alkaline phosphatase activity and labels
- Compatible with a variety of applications
- Use alone or in combination with NBT
- White solid soluble in DMF

For your convenience, we offer a [BCIP Red/NBT kit \(10005\)](#) 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:



References

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

This datasheet was generated on January 8, 2026 at 10:40:53 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/bcip-red-5-bromo-6-chloro-3-indoxyl-phosphate-p-toluidine-salt/>