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

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

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 colormetric 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 <u>BCIP Red/NBT kit (10005)</u> that contains both reagents. 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.

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 November 10, 2025 at 03:36:13 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/