

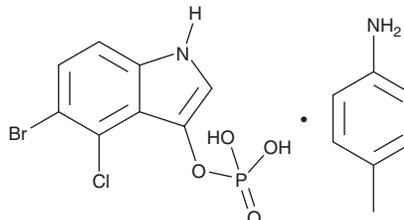
PRODUCT INFORMATION



BCIP (*p*-toluidine salt)

Item No. 34423

CAS Registry No.: 6578-06-9
Formal Name: 5-bromo-4-chloro-1H-indol-3-ol
3-(dihydrogen phosphate) compd. with
4-methylbenzenamine
Synonym: 5-Bromo-4-chloro-3-indoyl phosphate
MF: C₈H₆BrClNO₄P • C₇H₉N
FW: 433.6
Purity: ≥95%
UV/Vis.: λ_{max}: 232 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BCIP (*p*-toluidine salt) is supplied as a solid. A stock solution may be made by dissolving the BCIP (*p*-toluidine salt) in the solvent of choice, which should be purged with an inert gas. BCIP (*p*-toluidine salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BCIP (*p*-toluidine salt) in these solvents is approximately 30 mg/ml.

BCIP (*p*-toluidine salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BCIP (*p*-toluidine salt) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BCIP (*p*-toluidine salt) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BCIP is a colorimetric substrate for alkaline phosphatase.^{1,2} Upon enzymatic cleavage by alkaline phosphatase, an indoxyl intermediate is released and subsequently oxidized to produce water-insoluble indigo. It can be used to detect alkaline phosphatase activity by immunohistochemistry (IHC) or Western blot (WB).

References

1. Epstein, E., Wolf, P.L., Horwitz, J.P., *et al.* An indigogenic reaction for alkaline phosphatase in disk electrophoresis. *Am. J. Clin. Pathol.* **48(5)**, 530-534 (1967).
2. Wolf, P.L., Horwitz, J.P., Vasquez, J., *et al.* The indigogenic reaction for histochemical demonstration of alkaline and acid phosphatase. *Experientia* **23(3)**, 183-185 (1967).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/16/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM