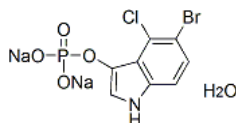
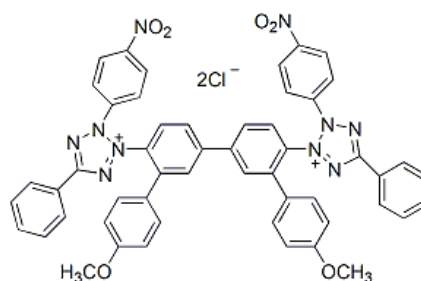


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: BCIP/NBT Kit**CATALOG #:** 10003**MOLECULAR INFORMATION:** **BCIP:**
 $C_8H_4BrClNNa_2O_4P$
Mwt: 370.5
[102185-33-1]**NBT:**
 $C_{40}H_{30}Cl_2N_{10}O_6$
Mwt: 817.7
[298-83-9]**PROPERTIES:**
Color & Form
Purity
Solubility**BCIP:**
White solid
≥ 95% by TLC
Soluble in water**NBT:**
Yellow solid
≥ 95% by TLC
Soluble in water and DMSO**STORAGE AND HANDLING:**

Stored desiccated at ≤ 4 °C upon receipt. Protect components from light, especially when in solution.

APPLICATION:

This kit contains 100 mg each of BCIP disodium salt (#10001) and the precipitation inducer NBT (nitro blue tetrazolium, #10008). BCIP is the most widely used chromogenic phosphatase substrate, which forms a dark blue precipitate on enzymatic hydrolysis. It is often used with the oxidant NBT (nitro blue tetrazolium chloride), which facilitates the precipitation. BCIP/NBT produces a dark blue/violet precipitate product that can be mounted with permanent or aqueous mounting medium.

Ref: *Biotechniques* **12**, 656(1992); *Histochemistry* **58**, 203(1978)**TOXICITY:**

Unknown

FIRST AID:

Potentially harmful; avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.