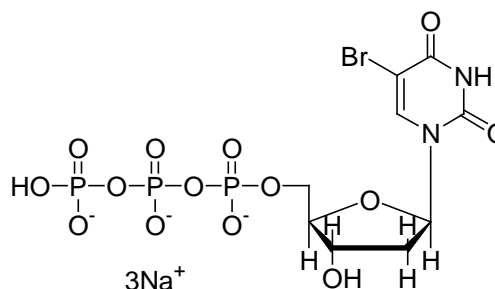


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Bromo-2'-deoxyuridine-5'-triphosphate, 10 mM in TE buffer (5-Bromo-dUTP)

CATALOG # 40025

MOLECULAR INFORMATION: C₉H₁₁BrN₂Na₃O₁₄P₃
MWt: 612.99
[102212-99-7]

**PROPERTIES:**

Color & Form Colorless solution
Solubility Soluble in water
Purity ≥ 95% by HPLC
Absorption/Emission
Extinction
Coefficient

STORAGE AND HANDLING: Store at -20 °C and protect from light

APPLICATION: 5-Bromo-dUTP is widely used in TUNEL assay to detect apoptosis cells,^{1,2} and it is also a good substrate for reverse transcriptase.^{3,4} Note: for PCR applications, Taq polymerase should be used with dUTP conjugates, because dUTP inhibits archaeal polymerases such as Pfu and Vent.^{5,6}

Ref.: 1) *Exp Cell Res* **222**, 28(1996); 2) *Cell Prolif* **28**, 571 (1995); 3) *Biotechnol Appl Biochem* **29**, 241(1999); 4) *Biotechnol Appl Biochem* **23**, 95(1996); 5) *Anal Biochem.* **211**,164(1993); 6) *PNAS* **99**,596(2002).

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

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.