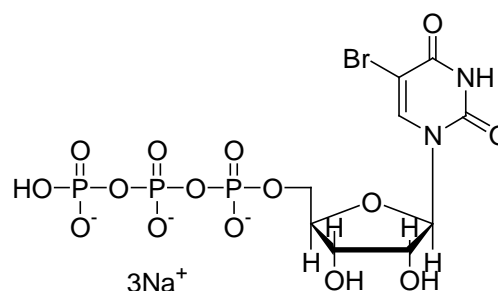


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Bromouridine-5'-triphosphate, 10 mM in TE buffer (BrUTP)**CATALOG #:** 40026**MOLECULAR INFORMATION:** C₉H₁₁BrN₂Na₃O₁₅P₃
MW: 628.98**PROPERTIES:**

Color & Form	Clear colorless liquid
Purity	≥ 97% by HPLC
Solubility	Soluble in water
Absorption/Emission	N/A

STORAGE AND HANDLING:

Store desiccated at -20 °C. Protect from light.

APPLICATION:BrUTP can be enzymatically incorporated into RNA. It is reported to be an excellent substrate for RNA polymerase and has been used to monitor nucleolar transcription *in situ*.Ref.: 1) *J Biochem* (Tokyo) **96**, 1501(1984); 2) *Mol Biol Cell* **10**, 211(1999); 3) *Histochem Cell Biol* **113**, 181 (2000)**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.