5-Bromo-dUTP, 10 mM solution

5-Bromo-dUTP (5-Bromo-2'-deoxyuridine-5'-triphosphate) is widely used in TUNEL assay to detect apoptosis cells, and also is a good substrate for reverse transcriptase.



Product attributes

CAS number

102212-99-7

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

Product Description

5-Bromo-dUTP (5-Bromo-2'-deoxyuridine-5'-triphosphate) is widely used in TUNEL assay to detect apoptotic cells (1,2), and it is also a good substrate for reverse transcriptase (3,4).

- 10 mM solution in TE buffer
- Store at -20°C
- C₉H₁₁BrN₂Na₃O₁₄P₃
- MW: 612.99
- [102212-99-7]

References

- 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).

This datasheet was generated on November 5, 2025 at 12:03:33 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-bromo-2-deoxyuridine-5-triphosphate-10-mm-in-te-buffer-5-bromo-dutp/