

## 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.



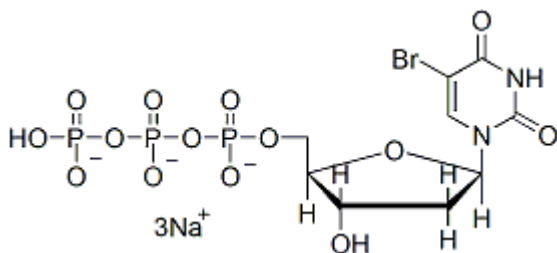
### Product attributes

|            |             |
|------------|-------------|
| CAS number | 102212-99-7 |
|------------|-------------|

## 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
- $C_9H_{11}BrN_2Na_3O_{14}P_3$
- 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 September 10, 2026 at 05:56:46 PM. 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/>