BrUTP, 10 mM Solution

BrUTP (5-Bromouridine-5'-triphosphate) can be enzymatically incorporated into RNA and detected with anti-BrdU antibodies.



Product attributes

CAS number

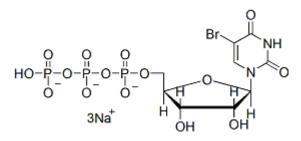
161848-60-8

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

Product Description

BrUTP (5-Bromouridine-5'-triphosphate) can be enzymatically incorporated into RNA and detected with anti-BrdU antibodies.

- 10 mM solution in TE buffer
- Store at -20°C
- C₉H₁₁BrN₂Na₃O₁₅P₃
- MW: 628.98
- [161848-60-8]



References

1) J Biochem (Tokyo) 96, 1501 (1984).

This datasheet was generated on November 8, 2025 at 11:38:24 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-bromouridine-5-triphosphate-10-mm-in-te-buffer-brutp/