

## BrUTP, 10 mM Solution

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



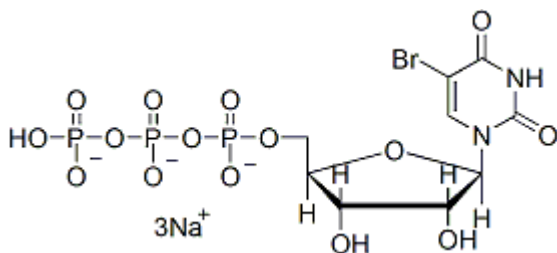
### Product attributes

CAS number	161848-60-8
------------	-------------

## 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
- $C_9H_{11}BrN_2Na_3O_{15}P_3$
- MW: 628.98
- [161848-60-8]



## References

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

This datasheet was generated on January 9, 2026 at 07:09:38 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/>