## Biotin-11-dUTP, 1 mM Solution

Biotin-11-dUTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3'-end terminal labeling. The number '11' is the number of atoms in the linker between biotin and dUTP.



## **Product attributes**

CAS number

86303-25-5

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

## **Product Description**

Biotin-11-dUTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3'-end terminal labeling. The number '11' is the number of atoms in the linker between biotin and dUTP. Biotium also offers biotin-16-dUTP (40022) and biotin-20-dUTP (40030). The length of the linker affects the incorporation efficiency of the biotin-dUTP probe into DNA using DNA polymerases, and it also affects biotin/avidin or biotin/streptavidin. In general, the shorter the linker, the more efficiently the biotin-dUTP is incorporated into DNA by DNA polymerases. On the other hand, the longer the linker, the better biotin can interact with avidin or streptavidin. Biotin-11-dUTP is also available as a lyophilized solid (40029-1).

- 1 mM solution in pH 7.5 Tris-HCl buffer
- Store at -20°C
- C28H41N6O17P3S Li4
- MW: 886.5
- [86303-25-5]

This datasheet was generated on November 4, 2025 at 04:45:53 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-11-dutp-1-mm-in-ph-7-5-tris-hcl-buffer/