## Biotin-11-dUTP, Lyophilized Powder

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, 40022-1) and biotin-20-dUTP (catalog nos. 40030, 40030-1). 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, a longer linker should facilitate interaction between biotin and avidin or streptavidin.

- Lyophilized solid more suitable for long term storage. The product contains lyophilized TE buffer.
- Store at -20°C
- C28H41N6O17P3SLi4
- MW: 886.5
- [86303-25-5]

Biotin-11-dUTP is also available as a 1 mM solution in Tris buffer (40029).

This datasheet was generated on November 3, 2025 at 09:32:40 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-11-dutp-lyophilized-powder/">https://biotium.com/product/biotin-11-dutp-lyophilized-powder/</a>