

## 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
------------	------------

## 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
- C<sub>28</sub>H<sub>41</sub>N<sub>6</sub>O<sub>17</sub>P<sub>3</sub>SLi<sub>4</sub>
- MW: 886.5
- [86303-25-5]

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

