

Biotin-16-dUTP, Lyophilized Powder

Biotin-16-dUTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3' end terminal labeling. The terminal deoxynucleotidyl transferase (TdT)-mediated biotin-dUTP nick end-labeling (TUNEL) method has been commonly used for apoptosis studies.



Product attributes

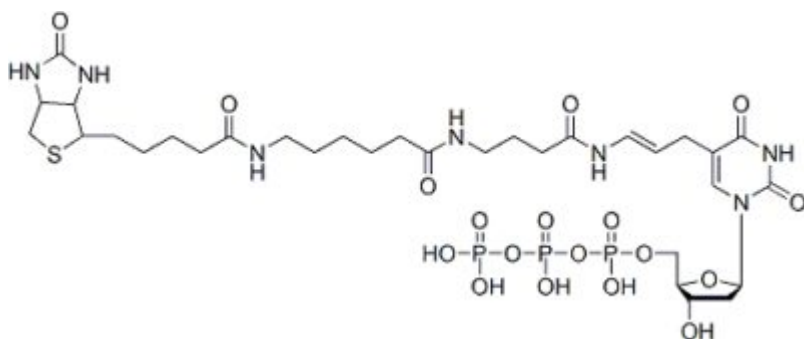
CAS number	136632-31-0
------------	-------------

Product Description

Biotin-16-dUTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3' end terminal labeling. The terminal deoxynucleotidyl transferase (TdT)-mediated biotin-dUTP nick end-labeling (TUNEL) method has been commonly used for apoptosis studies. The number '16' is the number of atoms in the linker between biotin and dUTP. Biotium also offers biotin-11-dUTP (catalog no. [40029-1](#)) and biotin-20-dUTP (catalog no. [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 suitable for long term storage; contains lyophilized TE buffer
- Store at -20 °C
- $C_{32}H_{48}N_7O_{18}P_3S Li_4$
- MW: 971.5
- [136632-31-0]

Biotin-16-dUTP is also available as a 1 mM solution in pH 7.5 Tris-HCl buffer (catalog no. [40022](#)).



References

1. Journal of Virological Methods (2013), <http://dx.doi.org/doi:10.1016/j.jviromet.2013.12.019>
2. The Open Genomics Journal 5, 18-29 (2012)

This datasheet was generated on December 19, 2025 at 05:04:29 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-16-dutp-lyophilized-powder/>