

## Biotin-11-dCTP, 1 mM Solution

Biotin-11-dCTP (Biotin-11-2'-deoxycytidine-5'-triphosphate, tetralithium salt) 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 dCTP.

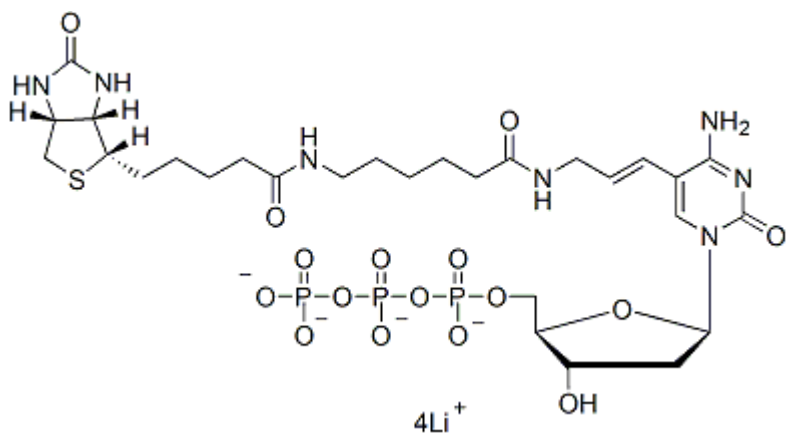


### Product attributes

### Product Description

Biotin-11-dCTP (Biotin-11-2'-deoxycytidine-5'-triphosphate, tetralithium salt) 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 dCTP.

- 1 mM solution in pH 7.5 TE buffer
- Store at -20°C
- $C_{28}H_{42}N_7O_{16}P_3SLi_4$
- MW: 885.42



This datasheet was generated on June 7, 2026 at 01:50:04 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-11-dctp-1-mm-in-ph-7-5-tris-hcl-buffer-biotin-11-2-deoxycytidine-5-triphosphate-tetralithium-salt/>