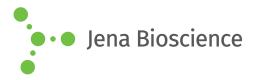
# **DATA SHEET**

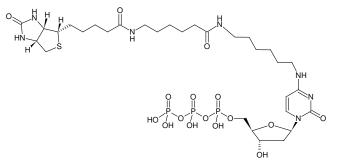




## Biotin-14-dCTP

Bio-14-dCTP, Triethylammonium salt

Cat. No.	Amount
NU-956-BIO14-S	200 µl (1 mM)
NU-956-BIO14-L	5 x 200 μl (1 mM)



Structural formula of Biotin-14-dCTP

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C<sub>31</sub>H<sub>54</sub>N<sub>7</sub>O<sub>16</sub>P<sub>3</sub>S (free acid)

Molecular Weight: 905.78 g/mol (free acid)

Exact Mass: 905.26 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: colorless to slightly yellow

Concentration: 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{max}$  272 nm,  $\epsilon$  13.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

#### Applications:

Incorporation into DNA/cDNA by

- Nick Translation with DNAse I/ DNA Polymerase I <sup>[1] & in-house data</sup> - Primer Extension with Klenow fragment <sup>[2]</sup>

### **Description**:

Biotin-14-dCTP is enzymatically incorporated into DNA/cDNA as substitute for its natural counterpart dCTP. The resulting Biotin-labeled DNA/cDNA probes are subsequently detected using streptavidin conjugated with horseradish peroxidase (HRP), alkaline phosphatase (AP), a fluorescent dye or agarose/magnetic beads. Optimal substrate properties for Nick Translation are ensured by a 14-atom linker attached to the N4 position of cytosine. For PCR incorporation experiments e.g. with *Taq* polymerase Biotin-11-dCTP (#NU-809-BIOX) or Biotin-16-dCTP (#NU-809-BIO16) are recommended whose Biotin moiety is attached to C5 position of cytidine via a 11-atom or 16-atom linker, respectively.

Recommended Biotin-14-dCTP/dCTP ratio for Nick Translation: 50% Biotin-14-dCTP/ 50% dCTP

Please note: The optimal final concentration of Biotin-14-dCTP may very depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Biotin-14-dCTP/dCTP ratio is recommended.

#### Selected References:

 Gebeyehu *et al.* (1987) Novel biotinylated nucleotide-analogs for labeling and colorimetric detection of DNA. *Nucleic Acids Res.* **15 (21)**:4513.
Lieberman-Aiden *et al.* (2009) Comprehensive mapping of long range interactions reveals folding principles of the human genome. *Science* **326** (5959):289.

