



## Biotin-7-dATP

N<sup>6</sup>-(6-Aminohexyl)-dATP - Biotin

Bio-7-dATP

N<sup>6</sup>-(6-Aminohexyl)-2'-deoxyadenosine-5'-triphosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-835-BIO-S	200 µl (1 mM)
NU-835-BIO-L	5 x 200 µl (1 mM)

### Applications:

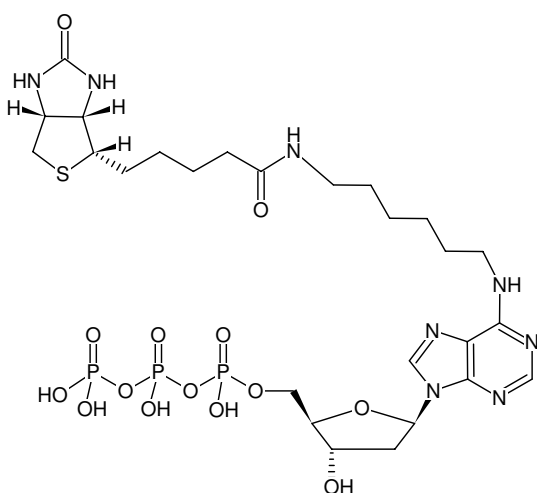
Incorporation by terminal deoxynucleotidyl transferase<sup>[1]</sup>

Incorporation by nick translation<sup>[2]</sup>

### Selected References:

[1] Flickinger *et al.* (1992) Differential incorporation of biotinylated nucleotides by terminal deoxynucleotidyl transferase. *Nucleic Acids Res.* **20**:2382.

[2] Gebeyehu *et al.* (1987) Novel biotinylated nucleotide-analogs for labeling and colorimetric detection of DNA. *Nucleic Acids Res.* **15 (21)**:4513.



Structural formula of Biotin-7-dATP

### For research use only!

**Shipping:** shipped on blue ice

**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>26</sub>H<sub>43</sub>N<sub>8</sub>O<sub>14</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 816.65 g/mol (free acid)

**Exact Mass:** 816.18 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** sterile clear aqueous solution

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 266 nm, ε 16.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)