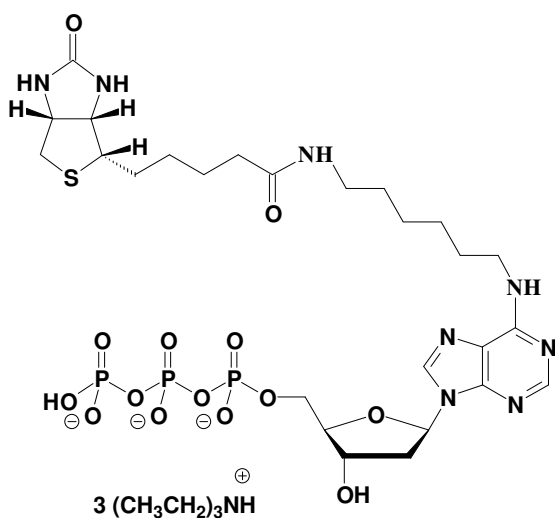


## N<sup>6</sup>-(6-Amino)hexyl-dATP - Biotin

N<sup>6</sup>-(6-Amino)hexyl-2'-deoxy-adenosine-5'-triphosphate - Biotin, Triethylammonium salt

| Cat. No.   | Amount            |
|------------|-------------------|
| NU-835-BIO | 50 $\mu$ l / 5 mM |



**Cat. No.:** NU-835-BIO

**Molecular Formula:** C<sub>26</sub>H<sub>40</sub>N<sub>8</sub>O<sub>14</sub>P<sub>3</sub>S (Anion)

**Molecular Weight:** 813.63 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 6 months after date of delivery.

**For research use only!**