

**Biotin-AMP**
 $\alpha$ -[(6-Aminoheptyl)imido]-AMP - Biotin, Triethylammonium salt

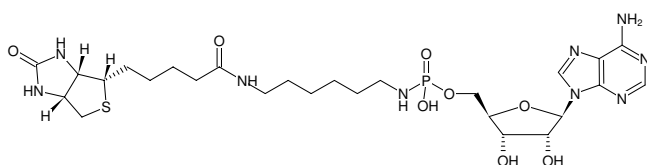
Cat. No.	Amount
NU-894-BIO-S	20 $\mu$ l (100 mM)
NU-894-BIO-L	5 x 20 $\mu$ l (100 mM)

**Applications:**5'-labeling of RNA by in vitro transcription<sup>[1]</sup>

**Please note:** This compound contains a phosphoramidite linkage which is hydrolyzed at pH <7.0.

**Selected References:**

[1] Huang *et al.* (2008) Synthesis of biotin-AMP conjugate for 5' biotin labeling of RNA through one-step in vitro transcription. *Nat. Protoc.* **3** (12):1848.



Structural formula of Biotin-AMP

**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>42</sub>N<sub>9</sub>O<sub>8</sub>PS (free acid)**Molecular Weight:** 671.71 g/mol (free acid)**Exact Mass:** 671.26 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5  $\pm$ 0.5

**Spectroscopic Properties:**  $\lambda_{\text{max}}$  260 nm,  $\epsilon$  15.4 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5)