

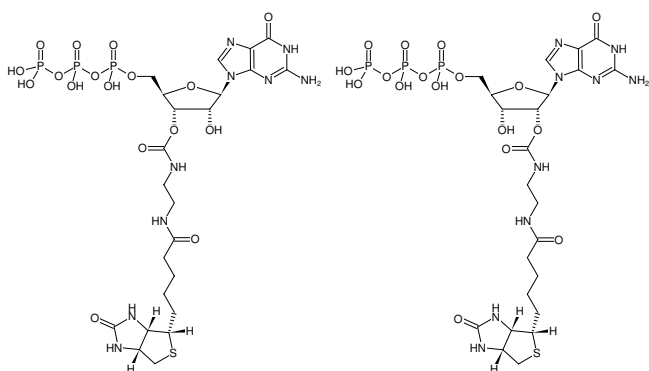
**EDA-GTP-Biotin**

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-820-BIO	50 µl (5 mM)

**Selected References:**

Buettner *et al.* (2014) Site-specific labeling of RNA at internal ribose hydroxyl groups: terbium-assisted deoxyribozymes at work. *J Am Chem Soc.* **136** (22):8131-7.



Structural formula of EDA-GTP-Biotin

**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>23</sub>H<sub>36</sub>N<sub>9</sub>O<sub>17</sub>P<sub>3</sub>S (free acid)**Molecular Weight:** 835.57 g/mol (free acid)**Exact Mass:** 835.12 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 5.0 mM - 5.5 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 252 nm, ε 13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)