



### EDA-ATPyS-Biotin

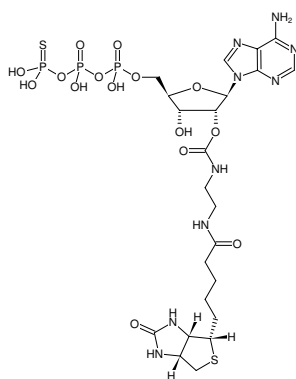
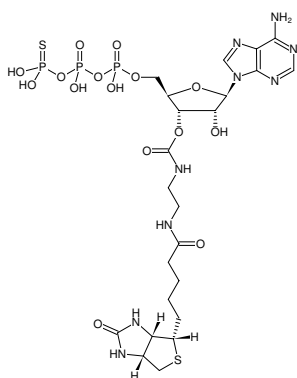
2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-1609-BIO	100 µl (1 mM)

**Please note:** For reasons of stability, please make sure that the pH value of a solution of this product never drops below 7.0.

The limited stability of labeled ATPγS and GTPγS nucleotides makes it impossible to keep stocks of them and to offer them as catalog products that are available immediately.

Therefore, any labeled ATPγS or GTPγS is freshly prepared per order. For this reason the price is higher than for stable nucleotide-dye conjugates which can be synthesized and kept in stock.



Structural formula of EDA-ATPyS-Biotin

**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:** 6 months after date of delivery

**Molecular Formula:** C<sub>23</sub>H<sub>36</sub>N<sub>9</sub>O<sub>15</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 835.64 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC), contains approx. 5 % EDA-ADP-Biotin

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)