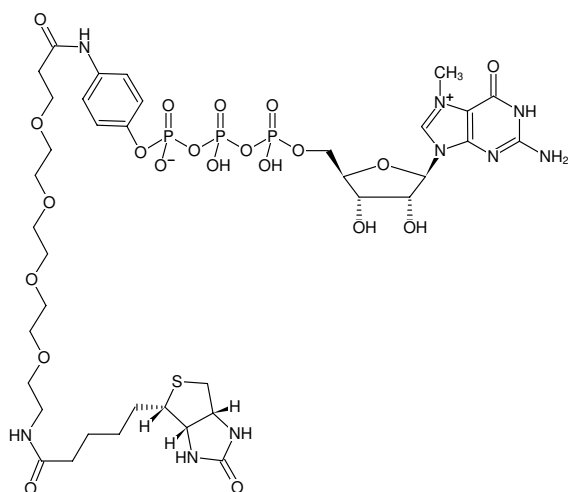



**m<sup>7</sup>GTP-γ-aminophenyl-PEG<sub>4</sub>-Biotin**

 7-Methylguanosine-5'-[γ-(4-aminophenyl)-PEG<sub>4</sub>]-5'-triphosphate-Biotin, Sodium salt

| Cat. No.   | Amount       |
|------------|--------------|
| NU-869-BIO | 50 µl (5 mM) |


 Structural formula of m<sup>7</sup>GTP-γ-aminophenyl-PEG<sub>4</sub>-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:** 12 months after date of delivery

**Molecular Formula:** C<sub>38</sub>H<sub>58</sub>N<sub>9</sub>O<sub>21</sub>P<sub>3</sub>S (free acid)

**Molecular Weight:** 1101.91 g/mol (free acid)

**Exact Mass:** 1101.27 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 5.0 mM - 5.5 mM

**pH:** 7.0 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 260 nm, ε 10.4 L mmol<sup>-1</sup> cm<sup>-1</sup>  
(Tris-HCl pH 7.5)