



## EDA-m<sup>7</sup>GDP - ATTO-Thio12

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO Thio12

Cat. No.	Amount
NU-827-THIO12	40 µl (1 mM)

**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>43</sub>H<sub>52</sub>N<sub>10</sub>O<sub>14</sub>P<sub>2</sub>S (free acid)

**Molecular Weight:** 1026.94 g/mol (free acid)

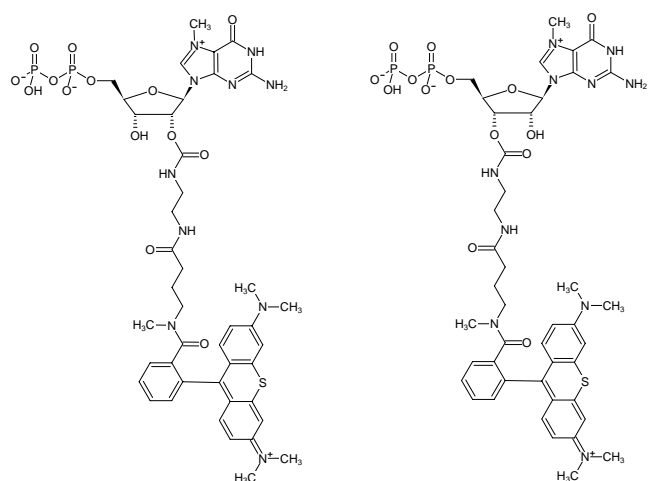
**Purity:** ≥ 95 % (HPLC)

**Form:** clear aqueous solution, pH 7.5 ± 0.5

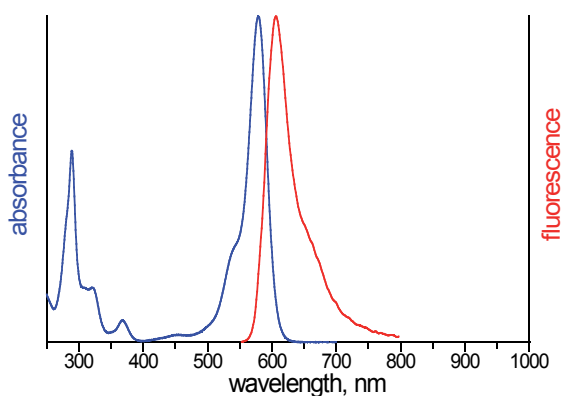
**Concentration:** 1 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 579 nm; λ<sub>em</sub> 609 nm; ε 110.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-m<sup>7</sup>GDP - ATTO-Thio12



excitation and emission spectrum of ATTO Thio12