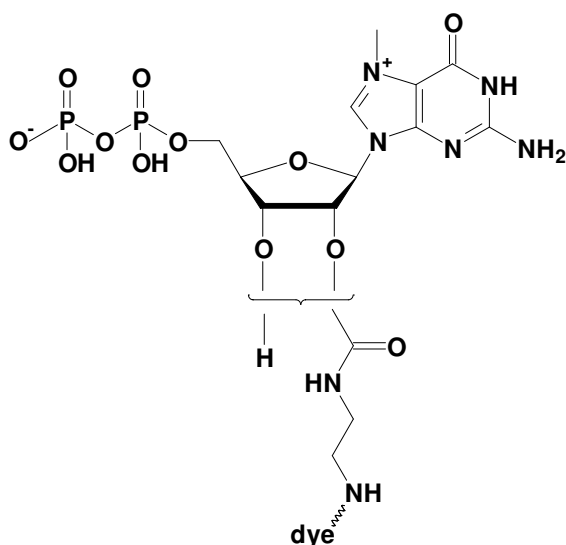


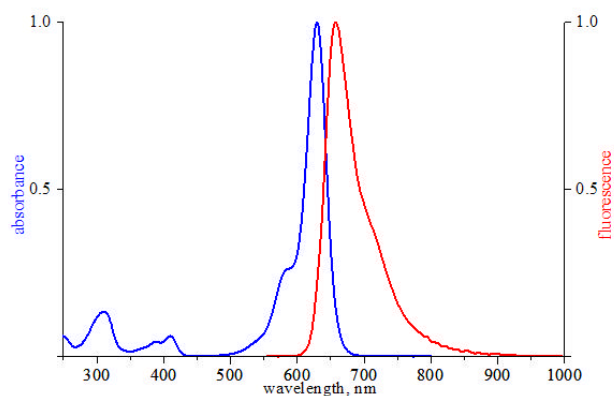
## EDA-m<sup>7</sup>GDP - ATTO 633

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with ATTO 633, Triethylammonium salt

Cat. No.	Amount
NU-827-633	40 $\mu$ l (1 mM)



structural formula of EDA-m<sup>7</sup>GDP



excitation and emission spectrum of ATTO 633

**Cat. No.:** NU-827-633

**Molecular Formula:** C<sub>14</sub>H<sub>24</sub>N<sub>7</sub>O<sub>12</sub>P<sub>2</sub> - ATTO 633 (free acid)

**Molecular Weight:** 1077.33 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  629 nm;  $\lambda_{em}$  657 nm;  $\epsilon$  130,000 cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**