



## EDA-m<sup>7</sup>GDP-DY-480XL

2'/3'-O-(2-Aminoethyl-carbamoyl)-7-methyl-guanosine-5'-diphosphate, labeled with DY 480XL, Triethylammonium salt

| Cat. No.     | Amount       |
|--------------|--------------|
| NU-827-480XL | 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>40</sub>H<sub>51</sub>N<sub>9</sub>O<sub>18</sub>P<sub>2</sub>S (free acid)

**Molecular Weight:** 1039.90 g/mol (free acid)

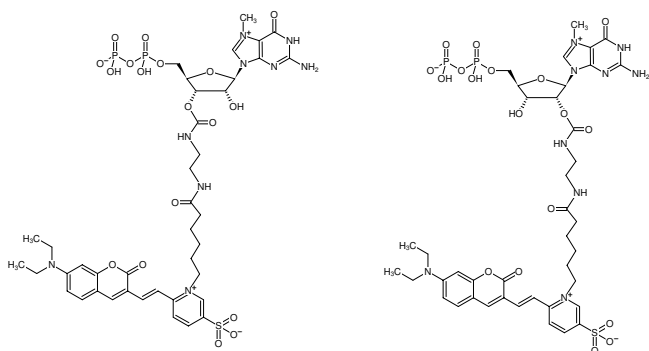
**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

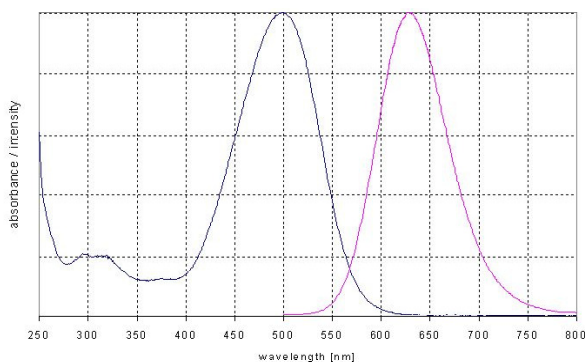
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 500 nm, λ<sub>em</sub> 630 nm, ε 50.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-m<sup>7</sup>GDP-DY-480XL



excitation and emission spectrum of DY 480XL