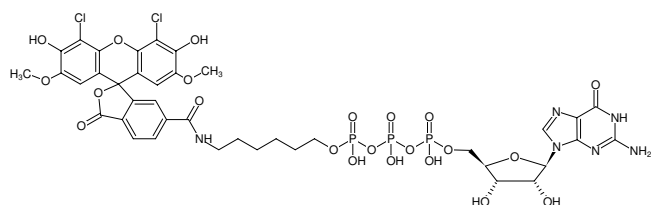




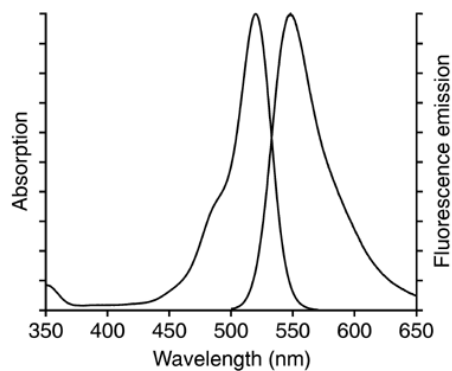
## $\gamma$ -(6-Aminoethyl)-GTP-6-JOE

$\gamma$ -(6-Aminoethyl)-guanosine-5'-triphosphate, labeled with 6-JOE, Triethylammonium salt

Cat. No.	Amount
NU-834-JOE	160 $\mu$ l (1 mM)



Structural formula of  $\gamma$ -(6-Aminoethyl)-GTP-6-JOE



excitation and emission spectrum of 6-JOE

### 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>39</sub>H<sub>41</sub>Cl<sub>2</sub>N<sub>6</sub>O<sub>22</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1108.09 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{exc}$  520 nm,  $\lambda_{em}$  548 nm,  $\epsilon$  75.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)