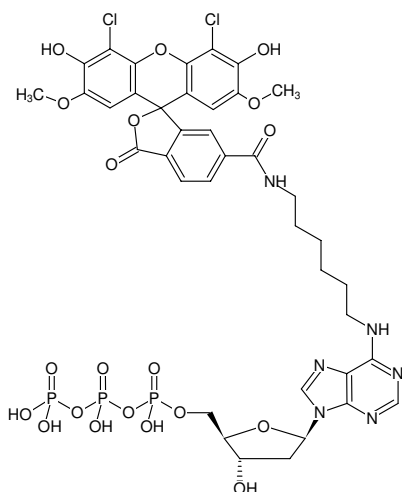




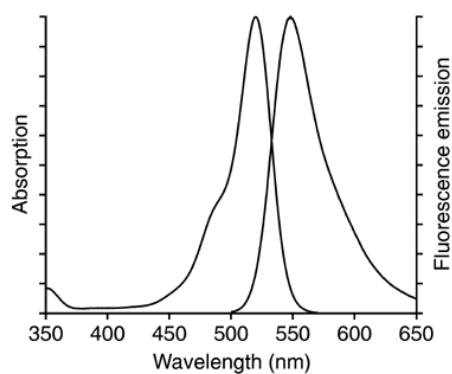
### N<sup>6</sup>-(6-Aminohexyl)-dATP-6-JOE

N<sup>6</sup>-(6-Aminohexyl)-2'-deoxyadenosine-5'-triphosphate, labeled with 6-JOE, Triethylammonium salt

Cat. No.	Amount
NU-835-JOE	160 µl (1 mM)



Structural formula of N<sup>6</sup>-(6-Aminohexyl)-dATP-6-JOE



excitation and emission spectrum of 6-JOE

#### For research use only!

**Shipping:** shipped on gel packs

**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>20</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1076.10 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> 520 nm, λ<sub>em</sub> 548 nm, ε 75.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)