



## N<sup>6</sup>-(6-Aminoethyl)-dATP-ATTO-532

N<sup>6</sup>-(6-Aminoethyl)-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 532, Triethylammonium salt

| Cat. No.   | Amount       |
|------------|--------------|
| NU-835-532 | 80 µl (1 mM) |

**For general laboratory use.**

**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>45</sub>H<sub>58</sub>N<sub>9</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1218.04 g/mol (free acid)

**Exact Mass:** 1217.24 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

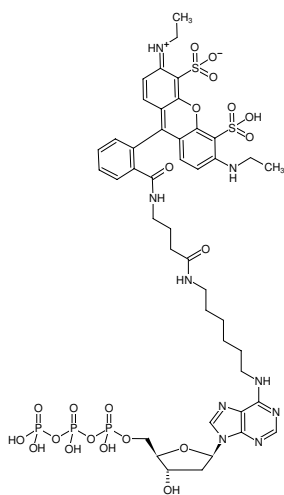
**Form:** solution in water

**Color:** orange

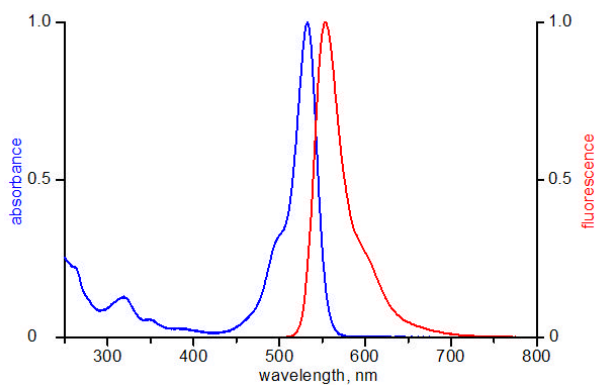
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 532 nm, λ<sub>em</sub> 552 nm, ε 115.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of N<sup>6</sup>-(6-Aminoethyl)-dATP-ATTO-532



excitation and emission spectrum of ATTO 532