



## EDA-ATP-ATTO-594

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-triphosphate, labeled with ATTO 594, Triethylammonium salt

Cat. No.	Amount
NU-808-594	40 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>54</sub>H<sub>67</sub>N<sub>10</sub>O<sub>23</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1381.22 g/mol (free acid)

**Exact Mass:** 1380.30 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

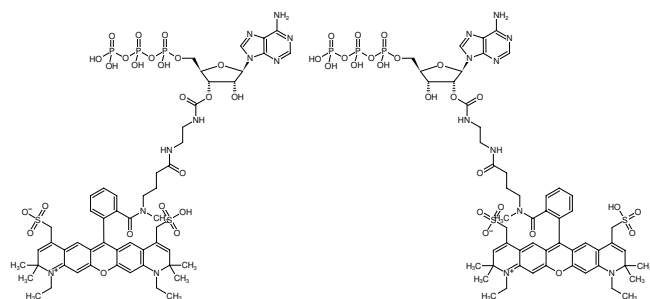
**Form:** solution in water

**Color:** red-violet

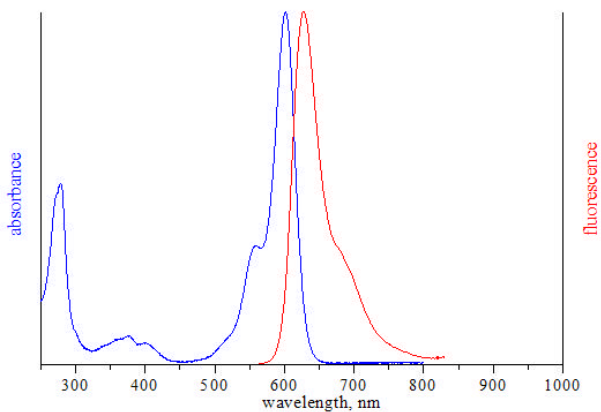
**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 603 nm, λ<sub>em</sub> 626 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of EDA-ATP-ATTO-594



excitation and emission spectrum of ATTO 594