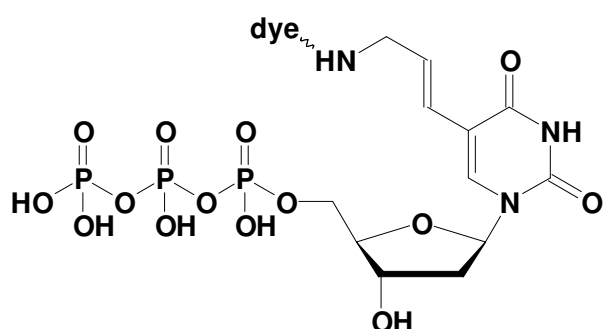


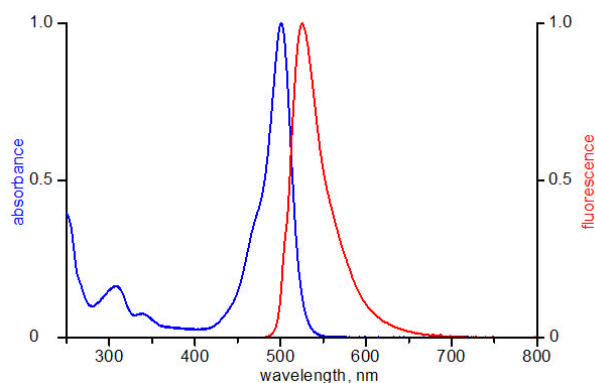
## Aminoallyl-dUTP - ATTO 488

5-(3-Aminoallyl)-2'-deoxy-uridine-5'-triphosphate, labeled with ATTO 488,  
 Triethylammonium salt

Cat. No.	Amount
NU-803-488	20 $\mu$ l / 1 mM



structural formula of Aminoallyl-dUTP



excitation and emission spectrum of ATTO 488

**Cat. No.:** NU-803-488

**Molecular Formula:** C<sub>12</sub>H<sub>20</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> - ATTO 488 (free acid)

**Molecular Weight:** 1094.22 (free acid)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  501 nm;  $\lambda_{em}$  523 nm;  $\epsilon$  90,000 cm<sup>-1</sup> M<sup>-1</sup>

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**

**Selected References:**

Vitali *et al.* (2010) Long nuclear-retained non-coding RNAs and allele-specific higher-order chromatin organization at imprinted snoRNA gene arrays. *J. Cell Sci.* **123** (1):70.