

## 7-Propargylamino-7-deaza-dATP - DY 751

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with DY 751,  
 Triethylammonium salt

Cat. No.	Amount
NU-1611-751	10 $\mu$ l / (1 mM)

**Cat. No.:** NU-1611-751

**Molecular Formula:** C<sub>57</sub>H<sub>71</sub>N<sub>6</sub>O<sub>20</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1317.26 (free acid)

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

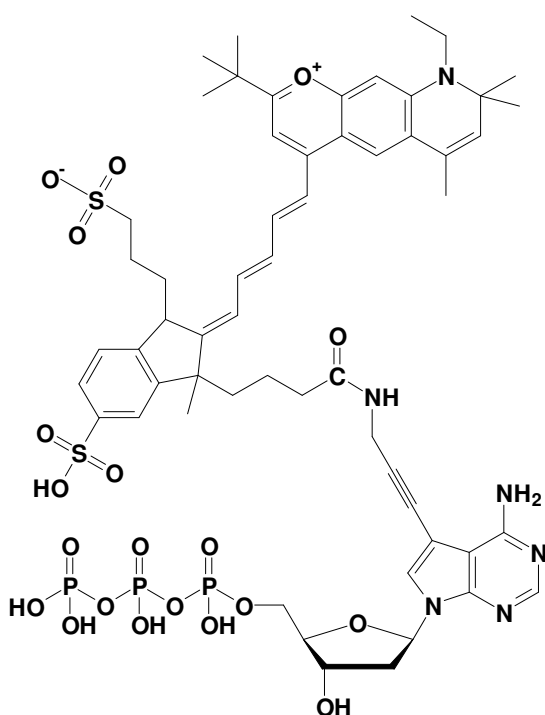
**Spectroscopic properties:**

$\lambda_{exc}$  751 nm;  $\lambda_{em}$  779 nm;  $\epsilon$  220,000 cm<sup>-1</sup> M<sup>-1</sup>

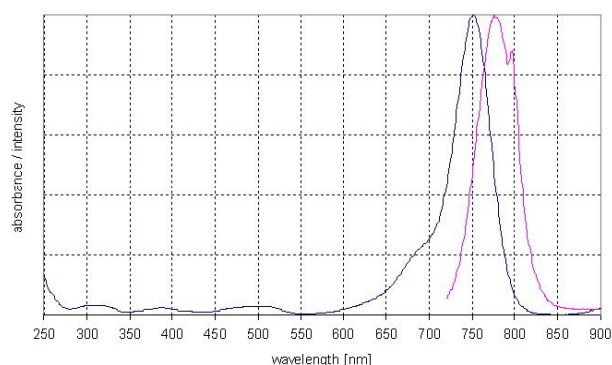
**Storage conditions:**

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!**



structural formula of 7-Propargylamino-7-deaza-dATP - DY 751



excitation and emission spectrum of DY 751