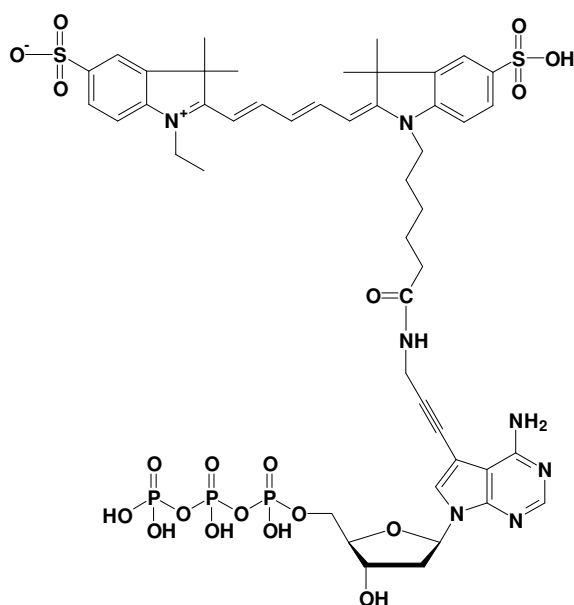


## 7-Propargylamino-7-deaza-dATP - Cy5

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, labeled with Cy5,  
 Triethylammonium salt

Cat. No.	Amount
NU-1611-CY5	60 $\mu$ l / (0.5 mM)



structural formula of 7-Propargylamino-7-deaza-dATP - Cy5

**Cat. No.:** NU-1611-CY5

**Molecular Formula:** C<sub>47</sub>H<sub>56</sub>N<sub>7</sub>O<sub>19</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1182.05 (free acid)

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

**Spectroscopic properties:**

$\lambda_{exc}$  649 nm;  $\lambda_{em}$  670 nm;  $\epsilon$  2,5 x 10<sup>5</sup> 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!**