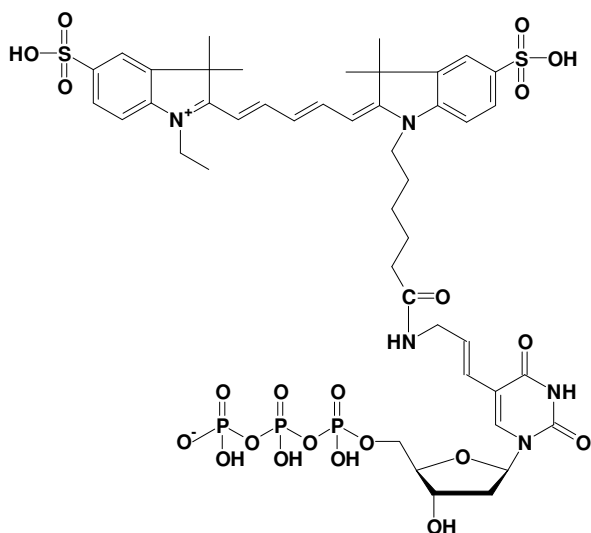


## Aminoallyl-dUTP - Cy5

5-(3-Aminoallyl)-2'-deoxy-uridine-5'-triphosphate, labeled with Cy5, Triethylammonium salt

Cat. No.	Amount
NU-803-CY5	30 $\mu$ l (1 mM)



structural formula of Aminoallyl-dUTP - CY5

**Cat. No.:** NU-803-CY5

**Molecular Formula:** C<sub>45</sub>H<sub>58</sub>N<sub>5</sub>O<sub>21</sub>P<sub>3</sub>S<sub>2</sub> (free acid)

**Molecular Weight:** 1162.01 (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!**