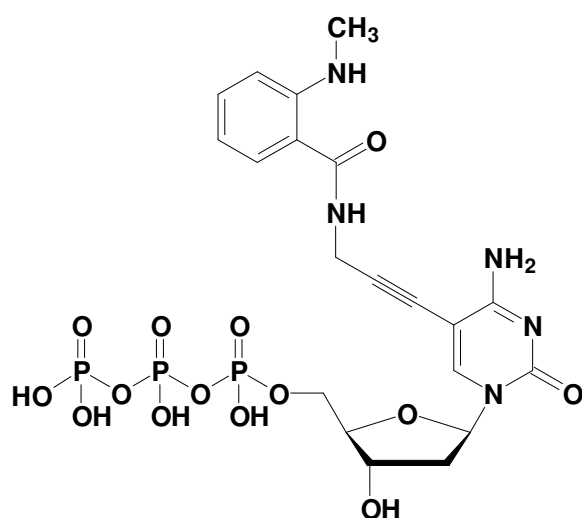


## 5-Propargylamino-dCTP - MANT

5-Propargylamino-2'-deoxy-cytidine-5'-triphosphate, labeled with MANT,  
 Triethylammonium salt

| Cat. No.   | Amount             |
|------------|--------------------|
| NU-809-MNT | 100 $\mu$ l (5 mM) |



structural formula of 5-Propargylamino-dCTP - MANT

**Cat. No.:** NU-809-MNT

**Molecular Formula:** C<sub>20</sub>H<sub>26</sub>N<sub>5</sub>O<sub>14</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 653.37 (free acid)

**Purity:** > 95%, clear solution, 10 mM Tris-HCl, pH 7.5

**Spectroscopic properties:**

$\lambda_{exc}$  335 nm;  $\lambda_{em}$  440 nm;  $\epsilon$  2,700 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!**