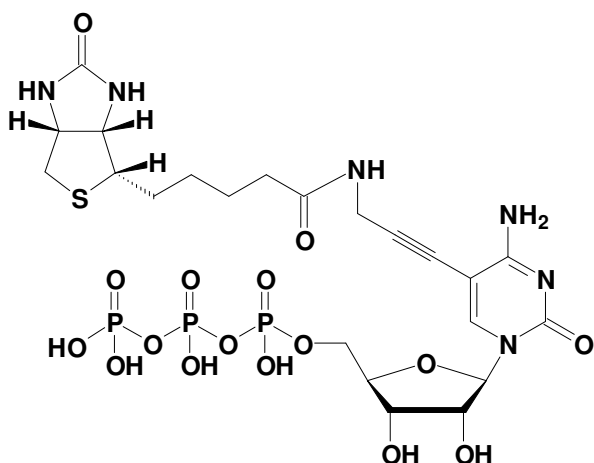


Biotin-4-CTP

5-Propargylamino-cytidine-5'-triphosphate - Biotin, Triethylammonium salt

| Cat. No. | Amount |
|------------|-------------------|
| NU-831-BIO | 60 μ l / 5 mM |



Cat. No.: NU-831-BIO

Molecular Formula: C₂₂H₃₃N₆O₁₆P₃S (free acid)

Molecular Weight: 762.51 (free acid)

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

Spectroscopic properties: λ_{\max} 294 nm;
 ϵ 9300

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 6 months after date of delivery.

For research use only!