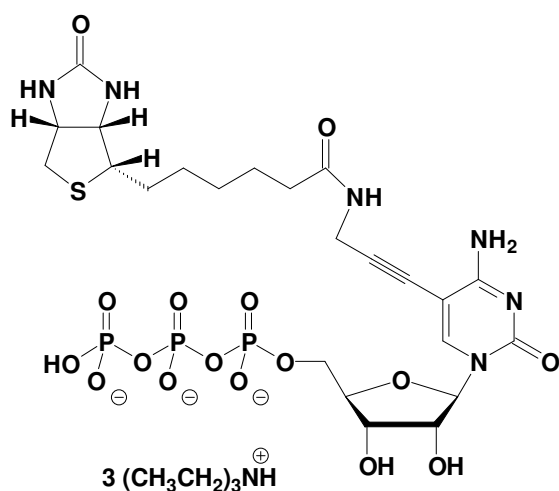


5-Propargylamino-CTP - Biotin

5-Propargylamino-cytidine-5'-triphosphate - Biotin

Cat. No.	Amount
NU-831-BIO	25 μ l / 5 mM



Cat. No.: NU-831-BIO

Molecular Formula: $\text{C}_{23}\text{H}_{32}\text{N}_6\text{O}_{16}\text{P}_3\text{S}$ (Anion)

Molecular Weight: 773.51 (Anion)

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

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!