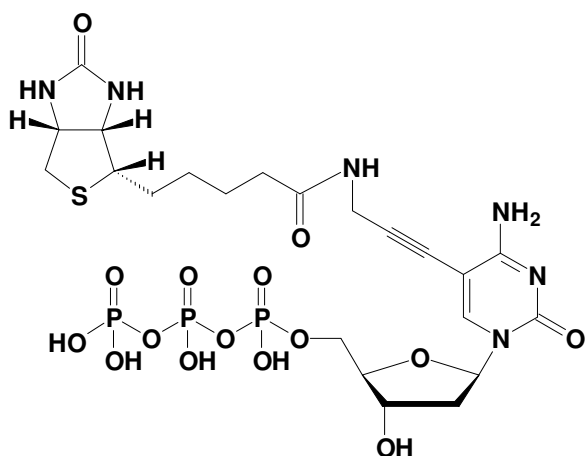


Biotin-4-dCTP

5-Propargylamino-2'-deoxy-cytidine-5'-triphosphate - Biotin, Triethylammonium salt

Cat. No.	Amount
NU-809-BIO	200 μ l / 5 mM



Cat. No.: NU-809-BIO

Molecular Formula: $C_{22}H_{33}N_6O_{15}P_3S$ (free acid)

Molecular Weight: 746.51 (free acid)

Purity: > 98%, 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!