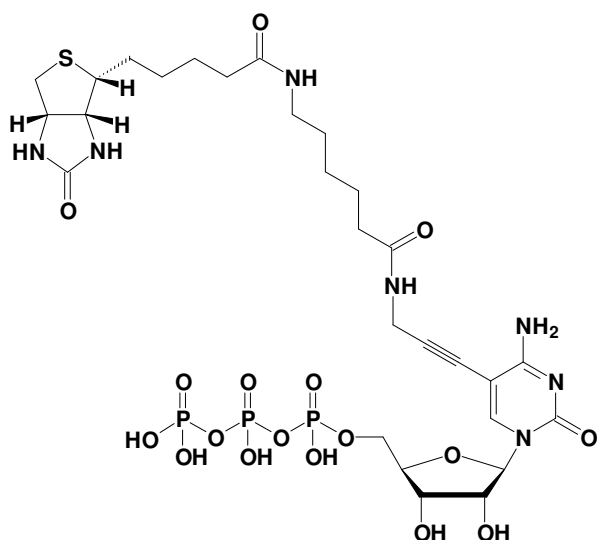


Biotin-11-CTP

γ -[N-(Biotin-6-amino-hexanoyl)]-5-propargylamino-cytidine-5'-triphosphate,
 Triethylammonium salt

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



Applications:

Microarray based hybridization detection
 [1, 2]

Ribonuclease protection assay^[3]

Selected References:

[1] Li *et al.* (2006) Identification of estrogen-responsive genes in parenchyma and fat pad of the bovine mammary gland by microarray analysis. *Physiol. Genomics* **27**:42.

[2] Jelinsky *et al.* (1999) Global response of *Saccharomyces cerevisiae* to alkylating agent. *Proc. Natl. Acad. Sci. USA* **96**: 1486.

[3] Kenrick *et al.* (1995) A homogeneous method to quantify mRNA levels: a hybridization of Rnase protection and scintillation proximity assay technologies. *Oxford University Press*:2947.

Cat. No.: NU-831-BIOX

Molecular Formula: C₂₈H₄₄N₇O₁₇P₃S (free acid)

Molecular Weight: 875.67 (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 12 months after date of delivery.

For research use only!