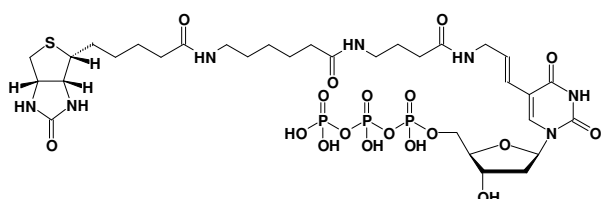


Biotin-16-dUTP

γ -[N-(Biotin-6-amino-hexanoyl-6-aminobutanoyl)]-5-(3-aminoallyl)-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-BIO16	200 μ l / 5 mM



Cat. No.: NU-803-BIO16

Molecular Formula: C₃₂H₅₂N₇O₁₈P₃S (free acid)

Molecular Weight: 947.78 (free acid)

Purity: > 98%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{\max} 240 nm; ϵ 10,700 M⁻¹cm⁻¹ Tris HCl 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 12 months after date of delivery.

For research use only!

Applications:

Comparative genomic hybridization^[1,2]

FISH^[2]

Nick-translation^[3, 4]

Selected References:

[1] Mattfeldt *et al.* (2002) Chromosomal regions in prostatic carcinomas studied by comparative genomic hybridization, hierarchical cluster analysis and self-organizing feature maps. *Anal. Cell. Pathology* **24**:167.

[2] Klink *et al.* (2010) Glioblastomas with oligodendroglial component - common origin of the different histological parts and genetic subclassifications. *Anal. Cell. Pathology/Cell. Oncology* **33**:37.

[3] Hermesen *et al.* (2005) Chromosomal changes in relation to clinical outcome in larynx and pharynx squamous cell carcinoma. *Cell. Oncology* **27**:191.

[4] Ghadimi *et al.* (2006) Distinct chromosomal profiles in metastasizing and non-metastasizing colorectal carcinomas. *Cellular Oncology* **28**:273.