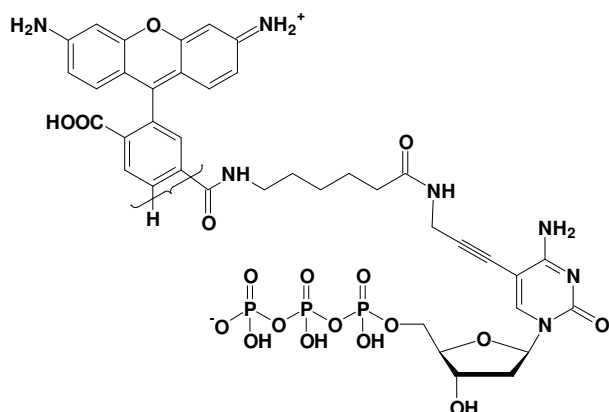


Rhodamine-12-dCTP

Rho-green-X-dCTP, RhG-dCTP, R110-dCTP

5/6-Rhodamine-X-5-propargylamino-2'-deoxy-cytidine-5'-triphosphate,
Triethylammonium salt

Cat. No.	Amount
NU-809-RHOX	30 μ l / (1 mM)



Cat. No.: NU-809-RHOX

Molecular Formula: C₃₉H₄₂N₇O₁₈P₃ (free acid)

Molecular Weight: 989.71 (free acid)

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

Spectroscopic properties:

λ_{exc} 505 nm; λ_{em} 530 nm; ϵ 85,000 cm⁻¹ M⁻¹

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:

PCR Incorporation^[1-3]

DNA-sequencing^[4, 5]

GISH^[6]

SNP-analysis^[7]

Selected References:

[1] Kaur *et al.* (2008) Genes controlling plant growth habit in *Leymus* (Triticeae): maize barren stalk1 (ba1), rice lax panicle, and wheat tiller inhibition (tin3) genes as possible candidates. *Funct Integr Genomics*. **8** (4):375.

[2] Brakmann *et al.* (2004) Optimal enzymes for single-molecule sequencing. *Curr Pharm Biotechnol*. **5** (1):119.

[3] Tasara *et al.* (2003) Incorporation of reporter molecule-labeled nucleotides by DNA polymerases. II. High-density labeling of natural DNA. *Nucleic Acids Res*. **31** (10):2636.

[4] Ramanathan *et al.* (2005) High density polymerase-mediated incorporation of fluorochrome-labeled nucleotides. *Anal. Biochem*. **337**:1.

[5] Seo *et al.* (2005) Four-color DNA sequencing by synthesis of a chip using photocleavable fluorescent nucleotides. *PNAS USA* **102**:5926.

[6] Park *et al.* (2007) Genomic changes in bilateral breast cancer by comparative genomic hybridisation. *Clin. Exp. Med*. **7**:1.

[7] Rockenbauer *et al.* (2005) SNP genotyping using microsphere-linked PNA and flow cytometric detection. *Cytometry, Part A* **64A**:80.