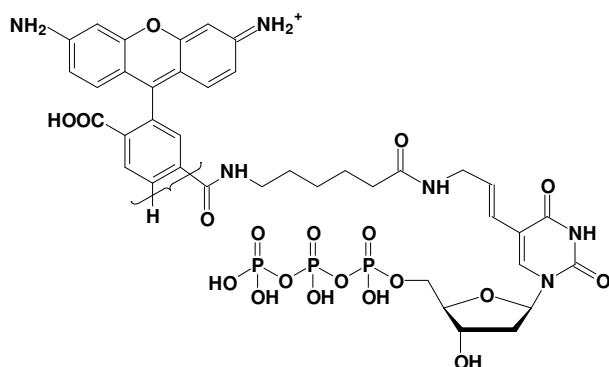


Rhodamine-12-dUTP

Rho-green-X-dUTP, RhG-dUTP, R110-dUTP

5/6-Rhodamine-X-5-Aminoallyl-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-RHOX	30 μ l / 1 mM



Selected References:

- [1] Zhong *et al.* (2010) Dual-color, Break-apart FISH Assay on Paraffin-embedded Tissues as an Adjunct to Diagnosis of Xp11 Translocation Renal Cell Carcinoma and Alveolar Soft Part Sarcoma. *Am. J. Surg. Pathol.* **34 (6)**:757.
- [2] 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.
- [3] Anderson *et al.* (2005) Incorporation of reporter-labeled nucleotides by DNA polymerases. *BioTechniques* **38 (2)**:257.

Cat. No.: NU-803-RHOX

Molecular Formula: C₃₉H₄₄N₆O₁₉P₃ (free acid)

Molecular Weight: 993.72 (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:

FISH^[1]

PCR Incorporation^[2-3]