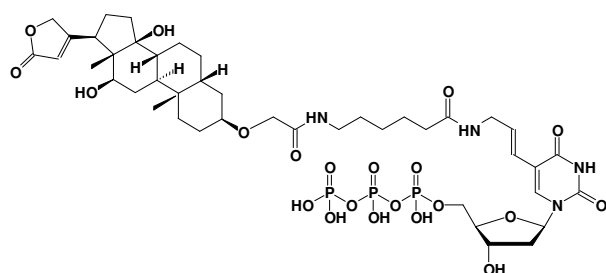


Digoxigenin-11-dUTP

Digoxigenin-X-5-aminoallyl-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

| Cat. No. | Amount |
|--------------|--------------------|
| NU-803-DIGXS | 25 μ l / 1mM |
| NU-803-DIGXL | 125 μ l / 1 mM |



Cat. No.: NU-803-DIGX

Molecular Formula: C₄₃H₆₅N₄O₂₁P₃ (free acid)

Molecular Weight: 1066.91 (free acid)

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

Spectroscopic properties: λ_{\max} 290 nm;
 ϵ 8780

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:

Nick translation^[1]

PCR^[1, 2]

Selected References:

[1] McCreery (1997) Digoxigenin labeling. *Mol Biotechnol.* **7(2)**:121.

[2] Jackson (1991) Detection of Shiga Toxin-Producing *Shigella dysenteriae* Type 1 and *Escherichia coli* by Using Polymerase Chain Reaction with Incorporation of Digoxigenin-11-dUTP. *J Clin Microbiol.* **29(9)**:1910.