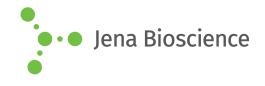
# **DATA SHEET**





## **DEAC-dUTP**

Aminoallyl-dUTP - DEAC Diethylaminocoumarin-5-dUTP

Diethylaminocoumarin-(5-aminoallyl)-2'-deoxyuridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-803-DEAC	25 μl (1 mM)

Structural formula of DEAC-dUTP

#### For general laboratory use.

**Shipping:** shipped on gel packs **Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient

temperature possible.

**Shelf Life:** 12 months after date of delivery **Molecular Formula:** C<sub>26</sub>H<sub>33</sub>N<sub>4</sub>O<sub>17</sub>P<sub>3</sub> (free acid) **Molecular Weight:** 766.48 g/mol (free acid)

Exact Mass: 766.11 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

Form: filtered solution (30 kDa) in 10 mM Tris-HCl

Color: yellow

Concentration: 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{exc}$  426 nm,  $\lambda_{em}$  480 nm,

 $\epsilon$  57.0 L mmol  $^{\text{-1}}$  cm  $^{\text{-1}}$  (Tris-HCl pH 7.5)

#### **Applications:**

Incorporation into DNA/cDNA by

- Nick Translation with DNAse I/DNA Polymerase I in-house data

### **Description:**

DEAC-dUTP is recommended for direct enzymatic labeling of DNA/cDNA by Nick Translation. It is incorporated as substitute for its natural counterpart dTTP. The resulting Dye-labeled DNA/cDNA probes are ideally suited for fluorescence hybridization applications such as FISH or microarray-based gene expression profiling.Optimal substrate properties and thus labeling efficiency is ensured by an optimized linker attached to the C5 position of uridine.

Recommended DEAC-dUTP/dTTP ratio for Nick Translation: 30-50% DEAC-dUTP/ 70-50% dTTP

Please note: Protect the Dye-labeled dUTP from exposure to light and carry out experimental procedures in low light conditions. The optimal final concentration of the Dye-labeled dUTP may very depending on the application and assay conditions. For optimal product yields and high incorporation rates an individual optimization of the Dye-labeled-dUTP/dTTP ratio is recommended.