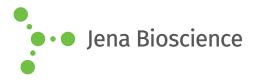
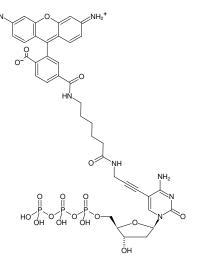
# **DATA SHEET**



## Rhodamine-12-dCTP

Rhodamine-X-5-propargylamino-dCTP Rho-green-X-dCTP, RhG-dCTP, R110-dCTP 5/6-Rhodamine-X-5-propargylamino-2'-deoxycytidine-5'-triphosphate, Triethylammonmium salt

Cat. No.	Amount
NU-809-RHOX-S	10 μl (1 mM)
NU-809-RHOX-L	5 x 10 μl (1 mM)



Structural formula of Rhodamine-12-dCTP

### 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>39</sub>H<sub>42</sub>N<sub>7</sub>O<sub>18</sub>P<sub>3</sub> (free acid)

Molecular Weight: 989.71 g/mol (free acid)

Exact Mass: 989.18 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

Form: solution in water

Color: orange

Concentration: 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{exc}$  505 nm,  $\lambda_{em}$  530 nm,  $\epsilon$  85.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Jena Bioscience GmbH Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com

### Applications:

Incorporation into DNA/cDNA by

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

#### **Description**:

Rhodamine-12-dCTP is recommended for direct enzymatic labeling of DNA/cDNA by Nick Translation. It is incorporated as substitute for its natural counterpart dCTP. 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 cytidine.

Recommended Rhodamine-12-dCTP/dCTP ratio for Nick Translation: 30-50% Rhodamine-12-dCTP/ 70-50% dCTP

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