

PRODUCT INFORMATION



Fluorescein-12-dUTP

Item No. 36232

Formal Name: 2'-deoxy-5-[3-[[[3',6'-dihydroxy-3-oxospiro[isobenzofuran-1(3H),9'-[9H]xanthen]-5(or 6)-yl]carbonyl]amino]-1-oxohexyl]amino]-1-propen-1-yl]-uridine 5'-(tetrahydrogen triphosphate), trisodium salt

Synonym: FITC-12-dUTP

MF: C₃₉H₃₈N₄O₂₁P₃ • 3Na

FW: 1,060.6

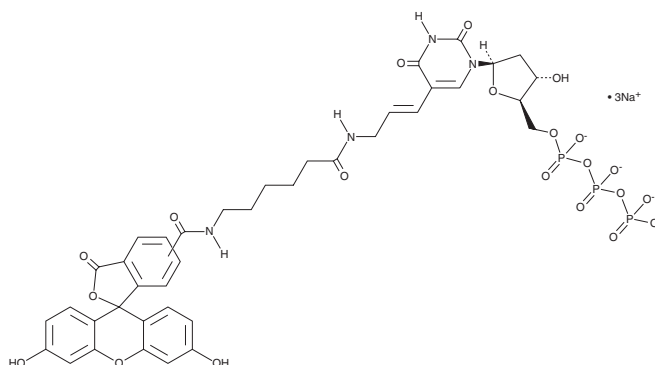
Purity: ≥95%

Ex./Em. Max: 498/517 nm

Supplied as: A solution in 1 mM tris buffer (pH 7.5)

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Fluorescein-12-dUTP is supplied as a solution in 1 mM tris buffer (pH 7.5). To change the solvent, simply evaporate the 1 mM tris buffer (pH 7.5) under a gentle stream of nitrogen and immediately add the solvent of choice. A stock solution may be made by dissolving the fluorescein-12-dUTP in the solvent of choice. Fluorescein-12-dUTP is soluble in water.

Description

Fluorescein-12-dUTP is a fluorescently tagged form of deoxy-UTP that can be used to generate fluorescently labeled DNA via reverse transcription, nick translation, random primed labeling, or PCR. It displays ex/em maxima of 498/517 nm, respectively. Fluorescein-12-dUTP has been used to detect nucleic acids *in situ*.¹

As a starting point for labeling, utilize approximately 0.05 µg/µl DNA, 25-50 U/ml DNA polymerase, and 10-100 µM fluorescein-12-dUTP in 10 mM Tris-HCl, pH 7.5, containing 1 mM EDTA, 5 mM NaCl, and 0.1 mM DTT, as well as dATP, dCTP, dGTP, and dTTP at 1 mM each in a total volume of 20 µl.

Reference

1. Koch, J., Mogensen, J., Pedersen, S., *et al.* Fast one-step procedure for the detection of nucleic acids *in situ* by primer-induced sequence-specific labeling with fluorescein-12-dUTP. *Cytogenet. Cell Genet.* 60(1), 1-3 (1992).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/18/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM