

PRODUCT INFORMATION



Fluorescein-12-dGTP

Item No. 36235

Formal Name: 4-((6-((3-(2-amino-7-((2R,4S,5R)-4-hydroxy-5-(((hydroxy((hydroxy(phosphonooxy)phosphoryl)oxy)phosphoryl)oxy)methyl)tetrahydrofuran-2-yl)-4-oxo-4,7-dihydro-3H-pyrrolo[2,3-d]pyrimidin-5-yl)prop-2-yn-1-yl)amino)-6-oxohexyl)carbamoyl)-2-(6-hydroxy-3-oxo-3H-xanthen-9-yl)benzoic acid

Synonym: Fluorescein-12-2'-deoxyguanosine-5'-triphosphate

MF: $C_{41}H_{41}N_6O_{20}P_3$

FW: 1,030.7

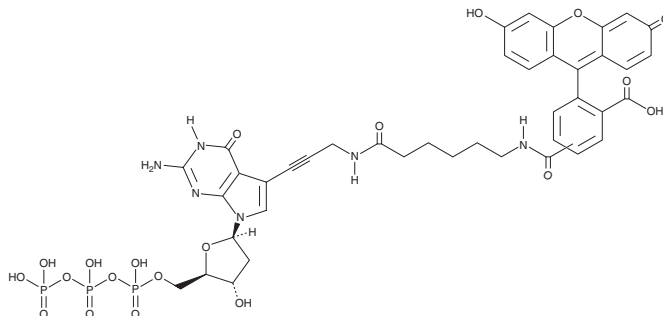
Purity: ≥95%

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-dGTP is supplied as a solution in 1 mM tris buffer (pH 7.5). Fluorescein-12-dGTP is soluble in water.

Description

Fluorescein-12-dGTP is a fluorescently labeled form of the nucleotide 2'-deoxyguanosine-5'-triphosphate (dGTP). It displays excitation/emission maxima of 498/517 nm, respectively. Fluorescein-12-dATP has been used to label DNA in FRET genotyping applications.¹

Reference

1. Duan, X., Yue, W., Liu, L., *et al.* Single-nucleotide polymorphism (SNP) genotyping using cationic conjugated polymers in homogeneous solution. *Nat. Protoc.* **4**(6), 984-991 (2009).

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, 07/30/2026

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