

Fluorescein-12-dUTP, molecular biology grade

Fluorescein-6-carboxaminocaproyl-[5-(3-aminoallyl)-2'-deoxyuridine-5'-triphosphate]

Catalog Number R0101

Pub. No. MAN0013126 Rev. B.00



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

Fluorescein-12-dUTP (fluorescein-5(6)-carboxaminocaproyl-[5-(3-aminoallyl)-2'-deoxyuridine-5'-triphosphate]) is supplied as 1 mM aqueous solution. The nucleotide is designed for enzymatic non-radioactive labeling of DNA during cDNA synthesis, PCR, nick-translation, random primed labeling, or primer extension. The nucleotide can be enzymatically incorporated into DNA with Reverse Transcriptases, *Taq* DNA Polymerase, phi29 DNA Polymerase, Klenow Fragment, *exo-*, Klenow Fragment and DNA Polymerase I.

General characteristics

λ_{\max} =495 nm, ϵ =70.0 $\times 10^3$ M⁻¹cm⁻¹ (pH 9.0).

Excitation maximum at 495 nm;

Emission maximum at 520 nm (pH 9.0).

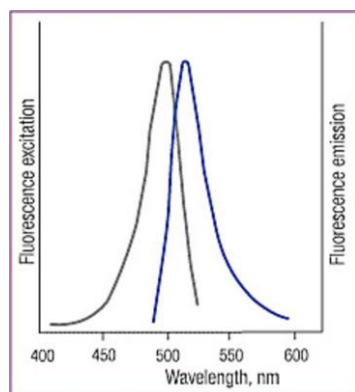


Fig.1. Normalized excitation-emission spectra of Fluorescein-12-dUTP.

Contents and storage

Contents	Formula	Molecular weight	Amount	Storage
Fluorescein-12-dUTP	C ₃₉ H ₄₁ N ₄ O ₂₁ P ₃	994.7	25 μ L, 25 nmol, 1 mM	-25 °C to -15 °C in the dark

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.