

Aminoallyl-dUTP, molecular biology grade

5-(3-aminoallyl)-2'-deoxyuridine-5'-triphosphate, sodium salt

Catalog Number R1101

Pub. No. MAN0012968 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.

Product description

Aminoallyl-dUTP (5-[3-aminoallyl]-2'-deoxyuridine-5'-triphosphate) is supplied as 50 mM aqueous solution titrated to pH 7.0-7.5 with NaOH. The nucleotide is designed for enzymatic indirect non-radioactive labeling of DNA during cDNA synthesis, PCR, nick-translation, random-primed labeling, or primer extension. Aminoallyl-dUTP can be enzymatically incorporated into DNA with Reverse Transcriptases, *Taq* DNA Polymerase, phi29 DNA Polymerase, Klenow Fragment, Klenow Fragment, exo- and DNA Polymerase I. The resulting amine-containing DNA can be subsequently labeled with any amine-reactive fluorescent dye, biotin or hapten.

General characteristics

λ_{\max} =240 nm, ϵ = 11.9×10^3 M⁻¹cm⁻¹ (pH 7.0);

λ_{\max} =290 nm, ϵ = 7.8×10^3 M⁻¹cm⁻¹ (pH 7.0).

Contents and storage

Contents	Formula	Molecular weight	Amount	Storage
Aminoallyl-dUTP	C ₁₂ H ₁₇ N ₃ O ₁₄ P ₃ Na ₃	589.2 (acid form: 523.2)	50 µL, 2.5 µmol, 50 mM	-25 °C to -15 °C

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.

For Research Use Only. Not for use in diagnostic procedures.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

ThermoFisher
SCIENTIFIC