

# Bio-11-UTP

Catalog Numbers AM8450, AM8451

Pub. No. MAN0013581 Rev. B



**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](http://thermofisher.com/support).

## Product description

Biotinylated UTP is intended for use as a substrate in *in vitro* transcription reactions with a variety of RNA polymerases, including T7 RNA polymerase, SP6 RNA polymerase, or T3 RNA polymerase. Biotinylated RNA probes may be used in place of radioactively labeled RNA probes in many applications, including *in situ* hybridization and chromosome mapping.

Biotinylated probes can be detected using a variety of methods based on streptavidin binding. Biotin-bound streptavidin can be detected directly, using fluorescently labeled streptavidin, or indirectly, using a streptavidin-conjugated enzyme assay or appropriately labeled or conjugated anti-streptavidin antibody.

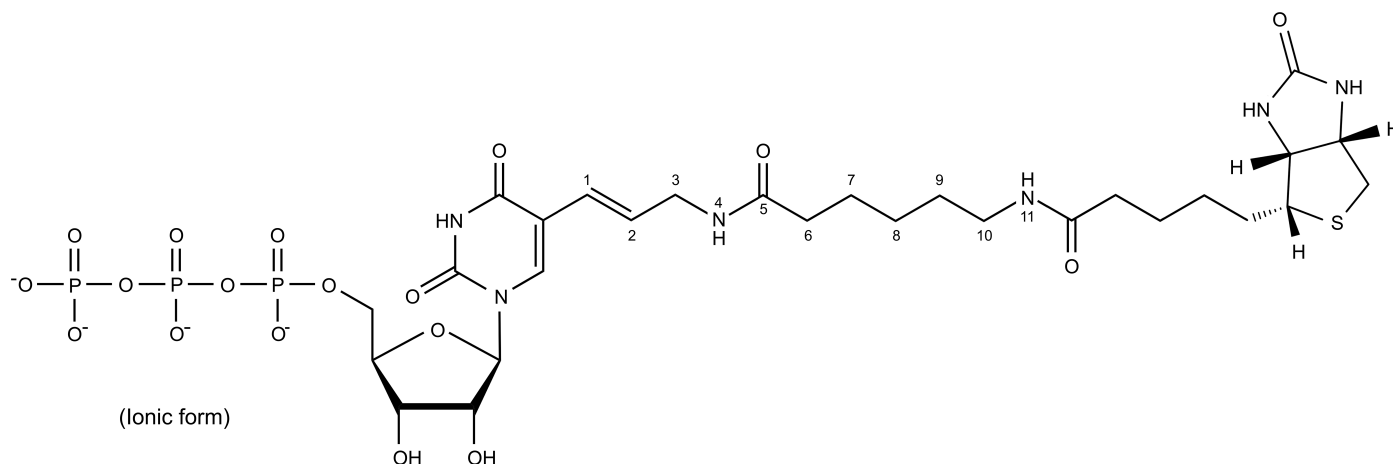


Figure 1 Structure

## Contents and storage

	Cat. No. <a href="#">AM8450</a>	Cat. No. <a href="#">AM8451</a>
Product Description	Bio-11-UTP (Biotin-11-Uridine-5'-triphosphate)	
Amount	250 nmol	2250 nmol
Volume	25 µL	30 µL
Concentration	10 mM	75 mM
Storage <sup>[1]</sup>	Store at or below –70°C. Avoid multiple freeze-thaw cycles.	
Storage Buffer	20 mM Tris-HCl pH 7.5	

<sup>[1]</sup> Dividing the product into single-use aliquots after the first use is recommended.

Specifications

Product specifications	
Molecular weight	944.62 (acid form)
UV absorption maximum	Extinction coefficient = 7,100 M <sup>-1</sup> cm <sup>-1</sup> at 289 nm
Purity	95% by HPLC at 289 nm Column: Ion exchange Buffer A: 0.01 M Na <sub>2</sub> HPO <sub>4</sub> pH 6.6 Buffer B: 0.7 M Na <sub>2</sub> HPO <sub>4</sub> pH 6.6 0–100% B over 35 minutes, 1.0 mL/minute

Limited product warranty

Life Technologies Corporation and its affiliates 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](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://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](http://thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN0013581 B

Revision	Date	Description
B	12 September 2024	Branding update.
A	9 June 2010	New manual for Bio-11-UTP.

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 these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2010-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.