

# TheraPure™ CTP, 200 mM Tris Solution

Catalog Number R00045T1

Pub. No. MAN0030281 Rev. A.00



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

## Contents and storage

Contents	Amount	Storage
TheraPure™ CTP, 200 mM, Tris Solution	200 µL	-25 °C to -15 °C

Thermo Scientific™ TheraPure™ CTP (cytidine 5'-triphosphate) is a superior quality nucleotide with exceptional purity of over 99%, supplied as a 200 mM Tris solution. It serves as a core building block for mRNA synthesis during development of mRNA therapeutics and vaccines.

TheraPure™ products are designed to support mRNA production and are recommended for process optimization and early development of mRNA vaccines and therapeutics: proof of concept, research, and preclinical studies.

## Features

- Animal-origin free (AOF) raw materials used in manufacturing process
- β lactam-free manufacturing technology
- Validated analytical tests methods
- Suitable for translational research
- Scalable: available as OEM product as TheraPure™ and TheraPure™ GMP.

Revision history: Pub. No. MAN0030281

Revision	Date	Description
A.00	2024-01-24	New

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](http://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](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).

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.

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

[thermofisher.com/support](http://thermofisher.com/support) | [thermofisher.com/askaquestion](http://thermofisher.com/askaquestion)

[thermofisher.com](http://thermofisher.com)

24 January 2024

