

# CTS™ Opti-MEM™ I

Catalog Numbers A4124801, A4124802

Pub. No. MAN0018638 Rev. A.0

 **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

Gibco™ CTS™ Opti-MEM™ I is a Minimal Essential Medium (MEM) containing insulin, transferrin, hypoxanthine, thymidine, and trace elements. CTS™ Opti-MEM™ I is recommended for use with cationic lipid transfection reagents.

## Contents and storage

Contents	Cat. No.	Amount	Storage	Shelf Life <sup>[1]</sup>
CTS™ Opti-MEM™ I	A4124801	100 mL	2°C to 8°C	18 months
	A4124802	500 mL	Protect from light	

<sup>[1]</sup> Shelf Life duration is determined from Date of Manufacture.

## Safety information

Human origin materials are non-reactive (donor level) for anti-HIV 1 & 2, anti-HCV, and HBsAg. Handle in accordance with established biosafety practices.

## Procedural guidelines

- We recommend CTS™ Opti-MEM™ I for use as a complexation buffer.
- For detailed instructions on using CTS™ Opti-MEM™ I for transfection, refer to CTS™ LV-MAX™ *Lentiviral Production System User Guide* (MAN0018450) at [thermofisher.com](http://thermofisher.com).

## Related products

Unless otherwise indicated, all materials are available through [thermofisher.com](http://thermofisher.com).

Item	Source
CTS™ Viral Production Cells	A3152801
CTS™ LV-MAX™ Production Medium	A4124001
CTS™ LV-MAX™ Transfection Kit	A4132601
CTS™ Opti-MEM™ I	A4124801

## Explanation of symbols

				
Temperature Limitation	Manufacturer	Batch code	Use By	Catalog Number
				
Caution, consult accompanying documents	Consult instructions for use	Keep away from light	Sterilized using aseptic processing techniques	European Community

## 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).



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.

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