# Leibovitz's L-15 Medium

Pub. No. MAN0018899 Rev. 2.0

**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.

# **Product description**

Leibovitz's L-15 Medium can be used to culture a variety of cells in suspension or adherent culture. It supports the growth of HEP-2 monkey kidney cells and primary explants of embryonic and adult human tissue. L-15 is buffered by phosphates and free base amino acids instead of sodium bicarbonate. This medium is designed for supporting cell growth in environments without CO<sub>2</sub> equilibration.

## Contents and storage

All quality control testing results are reported on lot-specific Certificate of Analysis available on our website: thermofisher.com.

Product	Cat. No.	Storage	Shelf life <sup>[1]</sup>
L-15 (LEIBOVITZ) MEDIUM	11415049 <sup>[2]</sup>	2°C to 8°C Protect from light	12 months
LEIBOVITZ's (1X) L 15 MED [+] L-Glutamine	11415064 <sup>[3]</sup>		
LEIBOVITZ's (1X) L 15 MED [+] L-Glutamine	11415114 <sup>[3]</sup>		
L-15 LEIBOVITZ W/GLUTAMAX-I	31415029 <sup>[2]</sup>		

<sup>[1]</sup> Shelf life is determined from Date of Manufacture. Do not use beyond the labeled expiration date.

🛛 🖬 Manufacturer: Life Technologies<sup>™</sup> Ltd. | 3 Fountain Drive, Inchinnan Business Park | Paisley PA49RF, Scotland, United Kingdom | Tel: +44 (0)141 81416305

[3] Manufacturer: Life Technologies Corporation | 3175 Staley Road | Grand Island, New York 14072 USA

# Guidelines for use

Ready-to-use commercial cell culture media undergoes strict quality control to ensure sterility, but may become contaminated while handling. Follow these guidelines for sterile handling to avoid contamination.

- Always wipe your gloved hands and work area with 70% ethanol.
- Wipe the outside of the containers, flasks, plates, and dishes with 70% ethanol before placing them in the cell culture hood.
- Use sterile pipette tips and pipettes to work with liquids, and use each pipette tip only once to avoid cross-contamination. Do not unwrap sterile pipettes until they are ready to be used. Keep pipettes and tips within the clean work area.
- Do not talk while performing sterile procedures and perform your cell culture as efficiently and carefully as possible to minimize contamination.

# Quality control

Standard evaluations for cell culture media are pH, osmolality, endotoxins and sterility testing for liquid products. All quality control testing results are reported on lot specific Certificate of Analysis available on our website: thermofisher.com.

# **Related products**

Unless otherwise indicated, all materials are available through thermofisher.com.

Item	Source
Gentamicin (50 mg/mL)	15750078
Amphotericin B	15290018
Penicillin-Streptomycin (10,000 U/mL)	15140122
PBS, pH 7.4	10010031
Fetal Bovine Serum (FBS)	16000044

# Labeling symbols

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



# 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.

Life Technologies Europe B.V.

The Netherlands

Kwartsweg 2, 2665 NN Bleiswijk

EC REP



Manufacturer: Life Technologies Corporation | 3175 Staley Road | Grand Island, New York 14072 USA

### Manufacturer:

Life Technologies<sup>™</sup> Ltd. | 3 Fountain Drive, Inchinnan Business Park | Paisley PA49RF, Scotland, United Kingdom | Tel: +44 (0)141 81416305

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.

#### Revision history: Pub. No. MAN0018899

Revision	Date	Description	
2.0 14 March 2022		The regulatory label was changed to "For Research Use or Further Manufacturing. Not for diagnostic use or direct administration into humans or animals."	
	The content required by previous IVD label was deleted or updated.		
	The document type was changed to Product Information Sheet to reflect changed scope of the document.		
1.0	12 November 2019	New document for Leibovitz's L-15 Medium.	

Important Licensing Information: This product 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.

@2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.