


RPMI 1640 Medium

Pub. No. MAN0018935 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](https://www.thermofisher.com/support).

Product description

Roswell Park Memorial Institute (RPMI) 1640 Medium is suitable for culture of a variety of mammalian cells, including HeLa, Jurkat, MCF-7, PC12, PBMC, astrocytes, and carcinomas. RPMI 1640 Medium is unique from other media because it contains the reducing agent glutathione and high concentrations of vitamins. RPMI 1640 Medium contains biotin, vitamin B12, and PABA, which are not found in Eagle's Minimal Essential Medium or Dulbecco's Modified Eagle Medium. In addition, inositol and choline are present in very high concentrations. RPMI 1640 Medium contains no proteins, lipids, or growth factors. Therefore, RPMI 1640 Medium requires supplementation, commonly with 10% fetal bovine serum (FBS). RPMI 1640 Medium uses a sodium bicarbonate buffer system (2.0 g/L) and requires a 5–10% CO₂ environment to maintain physiological pH.


Contents and storage


All quality control testing results are reported on lot-specific Certificate of Analysis available on our website: [thermofisher.com](https://www.thermofisher.com).


Product	Cat. No.	Storage	Shelf life ^[1]
RPMI 1640 (1X) [+] L-Glutamine [-] Phenol Red	11835030 ^[2] 11835055 ^[2] 11835063 ^[3]	2°C to 8°C Protect from light	12 months
RPMI Medium (1X) 1640 [+] L-Glutamine	11875085 ^[4] 11875093 ^[4] 11875101 ^[2] 11875119 ^[4] 11875127 ^[2] 11875135 ^[4] 11875168 ^[2] 11875176 ^[2]	2°C to 8°C Protect from light	12 months
RPMI Medium (1X) 1640 [-] L-Glutamine	21870076 ^[4] 21870084 ^[4] 21870092 ^[4] 21870100 ^[4]	2°C to 8°C Protect from light	12 months
RPMI 1640 W/GLUT (1X)	21875034 ^[3] 21875042 ^[3] 21875059 ^[3]	2°C to 8°C Protect from light	12 months
RPMI 1640 (1X) [+] L-Glutamine [+] HEPES	22400071 ^[2] 22400089 ^[4] 22400097 ^[2] 22400105 ^[4] 22400121 ^[2] 22400197 ^[2]	2°C to 8°C Protect from light	12 months

Product	Cat. No.	Storage	Shelf life ^[1]
RPMI 1640 W/O L-GLUTAMINE	31870017 ^[3] 31870025 ^[3]	2°C to 8°C Protect from light	12 months
RPMI 1640 W/O PHENOL RED	32404014 ^[3]	2°C to 8°C Protect from light	12 months
RPMI 1640 W/25MMHEPES W/OL-GLUT	42401018 ^[3] 42402016 ^[3]	2°C to 8°C Protect from light	12 months
RPMI MEDIUM 1640 W/HEPES	52400017 ^[3] 52400025 ^[3]	2°C to 8°C Protect from light	12 months
RPMI 1640 W/GLUTAMAX-I (1X)	61870010 ^[3] 61870036 ^[2] 61870127 ^[2] 61870143 ^[2] 61870150 ^[2]	2°C to 8°C Protect from light	12 months
RPMI 1640 W/HEPES W/GLUTAMAX-I	72400013 ^[3] 72400021 ^[3]	2°C to 8°C Protect from light	12 months
RPMI (1X) + GlutaMAX -I	72400047 ^[2] 72400120 ^[2] 72400146 ^[2] 72400153 ^[2]	2°C to 8°C Protect from light	12 months

^[1] Shelf life is determined from Date of Manufacture. Do not use beyond the labelled expiration date.

^[2]  Manufacturer: Life Technologies Corporation | 3175 Staley Road | Grand Island, New York 14072 USA

^[3]  Manufacturer: Life Technologies™ Ltd. | 3 Fountain Drive, Inchinnan Business Park | Paisley PA49RF, Scotland, United Kingdom | Tel: +44 (0)141 81416305

^[4]  Dual manufactured.

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

Product	Source
Gentamicin 50 mg/mL	15750078
Gibco Amphotericin B	15290018
Penicillin Streptomycin 10,000 U/mL	15140122
PBS, pH 7.4	10010031
Phytohemagglutinin, M form (PHA-M)	10576015
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.



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



Life Technologies Europe B.V.
Kwartsweg 2, 2665 NN Bleiswijk
The Netherlands



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



Manufacturer:
Dual manufactured products

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. MAN0018935

Revision	Date	Description
2.0	14 March 2022	<ul style="list-style-type: none"> 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 RPMI 1640 Medium.

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.

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