

EpiLife™ Medium

Catalog Number MEPI500CA

Pub. No. MAN0001626 Rev. 3.0

 **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

Gibco™ EpiLife™ Medium is a sterile, liquid tissue culture medium prepared with 60 µM calcium chloride for the growth of normal human epidermal keratinocytes and normal human corneal epithelial cells. This medium is intended for use as one component in a complete culture environment. EpiLife™ Medium is a basal medium containing essential and non-essential amino acids, vitamins, other organic compounds, trace minerals, and inorganic salts. EpiLife™ Medium does not contain antibiotics, antimycotics, hormones, growth factors, or proteins. This medium is HEPES and bicarbonate buffered and is designed for use in an incubator with an atmosphere of 5% CO₂/95% air. To support plating and long term proliferation of human keratinocytes or human corneal epithelial cells, EpiLife™ Medium must be supplemented with selected hormones and growth factors.

Content and storage

Table 1 EpiLife™ Medium (Cat. No. [MEPI500CA](#))

Contents	Amount	Storage ^[1,2]
EpiLife™ Medium	500 mL	4°C; Protect from light

^[1] EpiLife™ Medium is shipped at ambient temperature. Upon receipt, the medium should be stored at 4° C and should not be frozen. Protect from light. Several components of this tissue culture medium are light-labile, and we recommend that the medium not be exposed to light for lengthy periods of time. If the medium is warmed prior to use, do not exceed 37°C.

^[2] When stored in the dark at 4°C, the product is stable until the expiration date on the label.

Intended use

EpiLife™ Medium is intended for use in the routine culture of human epidermal keratinocytes or human corneal epithelial cells. Additional applications for use may include primary isolation of keratinocytes or corneal epithelial cells. See “Storage and stability of supplemented EpiLife™ Medium” on page 2 for more information.

Prepare supplemented EpiLife™ Medium

Depending on your application, obtain one of the growth supplement products. See “Storage and stability of supplemented EpiLife™ Medium” on page 2. Growth supplement are sold separately. To supplement one bottle of EpiLife™ Medium with the appropriate growth supplement, we recommend the following protocol:

1. Thaw the chosen growth supplement according to the instructions provided with that product. Make sure that the caps of all of the bottles are tight.
2. Gently swirl the bottle(s) of supplement. Avoid splashing the supplement into the cap of the bottle or causing the supplement to foam.
3. Obtain one bottle of EpiLife™ Medium with calcium. Wipe the outside of the containers with a disinfecting solution such as 70% ethanol or isopropanol and move the containers to a laminar flow culture hood.
4. To add the EpiLife™ Defined Growth Supplement (EDGS), Human Keratinocyte Growth Supplement (HKGS) or Human Corneal Growth Supplement (HCGS), use sterile technique in a laminar flow culture hood and transfer the entire contents of the bottle of supplement to the bottle of EpiLife™ Medium. To add the Human Keratinocyte Growth Supplement Kit (HKGS Kit), use sterile technique and transfer the desired amount of each component of the HKGS Kit to the bottle of medium.

- To add the EDGS, HKGS or HCGS, use sterile technique in a laminar flow culture hood and transfer the entire contents of the bottle of supplement to the bottle of medium. To add the HKGS Kit, use sterile technique and transfer the desired amount of each component of the HKGS Kit to the bottle of medium.

Note: Addition of less than the entire amount of any component may affect the performance of the supplemented

medium. If antibiotics/antimycotics are desired, add the antibiotic/antimycotic solution included in HKGS Kit using the same technique as above.

- Tightly cap the bottle of supplemented medium and swirl the contents to ensure a homogeneous solution. Avoid causing the medium to foam.

Storage and stability of supplemented EpiLife™ Medium

Once EpiLife™ Medium has been supplemented with EDGS, HKGS, HKGS Kit, or HCGS, the supplemented medium should be stored at 4°C and should not be frozen. When stored at 4°C, the supplemented medium is stable for 1 month.

EpiLife™ Extended-Lifespan Culture Systems			
	For keratinocytes:		For corneal epithelial cells:
	Defined	Undefined	Undefined
Basal Medium	EpiLife™ Medium (MEPI500CA) or EpiLife™ CF Kit (MEPICF500) or EpiLife™ CF/PRF Kit (MEPICFPRF500)	EpiLife™ Medium (MEPI500CA) or EpiLife™ CF Kit (MEPICF500) or EpiLife™ CF/PRF Kit (MEPICFPRF500)	EpiLife™ Medium (MEPI500CA) or EpiLife™ CF Kit (MEPICF500) or EpiLife™ CF/PRF Kit (MEPICFPRF500)
Growth supplement	EDGS ^[1] (S0125)	HKGS (S0015) or HKGS Kit (S001K)	HCGS (S0095)
Coating Matrix Kit	Coating Matrix Kit (R-011-K)	N/A	N/A
Subculture reagent	Trypsin/EDTA Solution (R001100)	Trypsin/EDTA Solution (R001100)	Trypsin/EDTA Solution (R001100)
Subculture reagent	Defined Trypsin Inhibitor (R007100)	Trypsin Neutralizer (R002100)	Trypsin Neutralizer (R002100)

^[1] For optimal cell growth, EDGS should be used in conjunction with Coating Matrix Kit and Defined Trypsin Inhibitor.

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 Corporation | 3175 Staley Road | Grand Island, New York 14072 USA
For descriptions of symbols on product labels or product documents, go to 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: 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.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.