

Mouse (ICR) Inactivated Embryonic Fibroblasts

Description

Mouse (ICR) Inactivated Embryonic Fibroblasts are used as feeder layers for culturing embryonic stem cells (ESCs), including mouse and human ESCs, in their undifferentiated state. The fibroblasts were isolated from ICR mouse embryos under sterile conditions, expanded for up to two passages, and then mitotically inactivated by γ -irradiation. The growth-arrested feeder layer supports the ESC culture by providing nutrients, growth factors, and matrix components, and it enables ESCs to survive and proliferate more readily in culture.

Product	Catalog No.	Amount	Storage
Mouse (ICR) Inactivated Embryonic Fibroblasts	A24903	1 mL ($\geq 2 \times 10^6$ viable cells/mL)	Liquid nitrogen (vapor phase)

Product Use

For Research Use Only. Not for use in diagnostic procedures.

Important Information

Follow the guidelines below to use inactivated mouse embryonic fibroblasts (MEFs) as feeder layers to culture mouse and human embryonic stem cells (ESCs).

- **All solutions and equipment that come in contact with the cells must be sterile.** Always use proper aseptic technique and work in a laminar flow hood.
- Make sure to prepare MEF feeder layer at least one day before culturing ESCs.
- After thawing, transfer MEFs into pre-warmed medium.
- Plate MEFs on culture vessels coated with 1× Attachment Factor (AF) solution.
- Use MEF dishes or plates within 3–4 days of preparation.
- Before starting experiments, ensure that ESCs have been established (at least 1 passage), and also have some frozen ESC stocks on hand.

Safety Information

Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Culture Conditions

Media: Dulbecco's Modified Eagle medium (DMEM) containing 4.5 g/L glucose (Cat. no. 10569), and supplemented with 10% FBS, ES-cell qualified (Cat. no. 16141).

Culture Type: Adherent

Recommended Substrate: Attachment Factor (Cat. no. S-006-100), which is a sterile 1× solution containing 0.1 % gelatin.

Temperature Range: 36°C to 38°C

Incubator Atmosphere: Humidified atmosphere of 5% CO₂.

Coat Culture Vessels with Attachment Factor

1. Cover the whole surface of each culture vessel with Attachment Factor (AF) solution and incubate the vessels for 30 minutes at 37°C or for 2 hours at room temperature. Refer to the recommended coating volumes in the table, next page.
2. Using sterile technique in a laminar flow culture hood, completely remove the AF solution from the culture vessel by aspiration.

Note: It is not necessary to wash the culture surface before adding cells or medium.
3. You can use the coated vessels immediately or store them at room temperature for up to 24 hours.

Prepare MEF Medium

1. Prepare 500 mL of MEF medium by mixing the following components (pre-warmed in a 37°C, 5% CO₂ incubator):

D-MEM	450 mL
FBS	50 mL

Thaw MEFs

1. Remove the cryovial containing inactivated MEFs from the liquid nitrogen storage tank.
2. Briefly roll the vial between hands to remove frost, and swirl it gently in a 37°C water bath.
3. When only a small ice crystal remains in the vial, remove it from water bath. Spray the outside of the vial with 70% ethanol before placing it in the cell culture hood.
4. Pipet the thawed cells gently into a 50-mL conical tube.
5. Add 10 mL of pre-warmed MEF medium **dropwise** to the cells while gently swirling the conical tube. Gently mix by pipetting up and down.

Note: Adding the medium slowly helps the cells to avoid osmotic shock.
6. Transfer entire cell suspension to a 15-mL conical tube and centrifuge at 200 × g for 5 minutes.
7. Aspirate the supernatant and resuspend the cell pellet in an appropriate volume of pre-warmed MEF medium.
8. Use an appropriate volume of the cell suspension to determine the viable cell number using your method of choice.

Plate MEFs

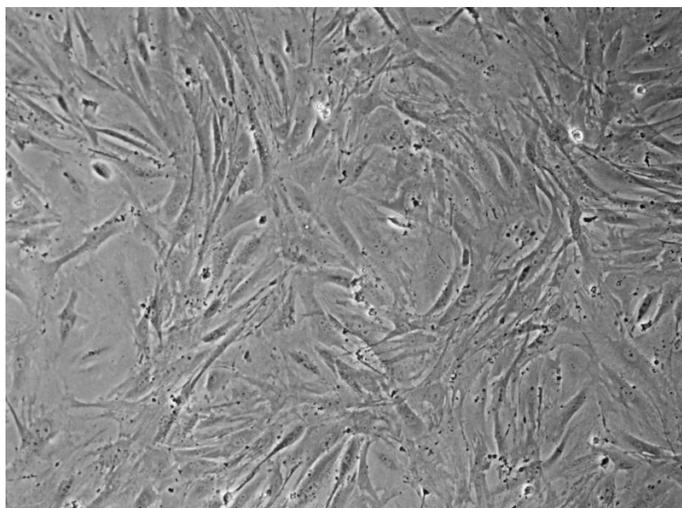
1. Aspirate the gelatin solution from the AF-coated culture vessels.
2. Add the appropriate amount of MEF medium into each culture vessel (refer to the table, next page).
3. Into each of these culture vessels, add the appropriate amount of MEF suspension (refer to the table, next page).

Note: The appropriate cell density should be optimized for the specific application. We recommend 3.0×10^4 MEFs/cm² as a good starting point.
4. Move the culture vessels in several quick back-and-forth and side-to-side motions to disperse the cells across the surface of the vessels.
5. Incubate the cells in a 37°C incubator with a humidified atmosphere of 5% CO₂.
6. Use the MEF culture vessels within 3–4 days after plating.

Vessel Size	AF Coating Volume	Volume of Media	Number of MEFs
96-well plate	0.1 mL	0.1 mL	1.0×10^4 /well
24-well plate	0.3 mL	0.5 mL	6.0×10^4 /well
12-well plate	0.5 mL	1 mL	1.1×10^5 /well
6-well plate	1 mL	2 mL	2.9×10^5 /well
60-mm dish	3 mL	5 mL	5.9×10^5
100-mm dish	9 mL	10 mL	1.8×10^6
25-cm ² flask	3 mL	5 mL	7.5×10^5
75-cm ² flask	9 mL	15 mL	2.3×10^6

Expected Results

The bright field image below shows Mouse (ICR) Inactivated Embryonic Fibroblasts plated at the recommended density on culture dishes coated with 1× AF solution. The image was taken with a 10× objective.



Related Products

Product	Cat. no.
Dulbecco's Modified Eagle Medium (D-MEM), high glucose	A10569
Dulbecco's PBS (DPBS) without Calcium and Magnesium	14190
Fetal Bovine Serum (FBS), ES-Cell Qualified	A16141
Attachment Factor	S-006-100

Explanation of Symbols and Warnings

The symbols present on the product label are explained below:

Temperature limitation	Use by	Batch code	Catalog number	Manufacturer

References

- Evans, M., Kaufman, M. (1981) Establishment in culture of pluripotent cells from mouse embryos. *Nature* 292, 154–156.
- Thomson, J.A., Itskovitz-Eldor, J., Shapiro, S.S., Waknitz, M.A., Swiergiel, J.J., Marshall, V.S., Jones, J.M. (1998) Embryonic stem cell lines derived from human blastocysts. *Science* 282, 1145–1147.
- Williams, R.L., Hilton, D.J., Pease, S., Willson, T.A., Stewart, C.L., Gearing, D.P., Wagner, E.F., Metcalf, D., Nicola, N.A., Gough, N.M. (1988) Myeloid leukaemia inhibitory factor maintains the developmental potential of embryonic stem cells. *Nature* 336, 684–687.

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 found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

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.

Made in USA

7335 Executive Way Frederick, MD 21704

For additional technical information such as Safety Data Sheets (SDS), Certificates of Analysis, visit www.lifetechnologies.com/support. For further assistance, email techsupport@lifetech.com

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

www.lifetechnologies.com

11 June 2014

life
technologies™