

Primary Mouse Embryonic Fibroblast Isolation Kit

Pub. No. MAN0016041

Rev A.0

Pub. Part No. 2162551.0

88279

Number	Description
88279	Pierce Mouse Embryonic Fibroblast (MEF) Isolation Kit , contains sufficient reagents to isolate MEFs from 50 embryonic mouse somatic tissues

Kit Contents:**DMEM for Primary Cell Isolation**, 500mL**MEF Isolation Enzyme (with papain)**, lyophilized, 5 vials**Hanks' Balanced Salt Solution (HBSS without Ca²⁺, Mg²⁺)**, 500mL**Storage:** Upon receipt store kit at 4°C. Product is shipped on ice packs.**Table of Contents**

Introduction	1
Important Product Information	2
Additional Materials Required	2
Material Preparation	2
Procedure for MEF Isolation	2
A. Enzyme Digestion of Mouse Embryonic Tissue	2
B. Cell Yield and Viability Determination	3
C. Plating and Culturing Isolated MEFs	4
Troubleshooting	5
Related Thermo Scientific Products	5
References	6

Introduction

The Thermo Scientific™ Pierce™ Mouse Embryonic Fibroblast (MEF) Isolation Kit provides a reliable and convenient method for the isolation and culture of MEFs (Figure 1). The kit consists of unique tissue-specific dissociation reagents and a validated protocol to ensure a high yield of viable MEFs when used by both experienced and non-experienced users.

MEFs isolated and cultured using the Pierce MEF Isolation Kit can be used as a cell culture model for a diverse range of studies such as gene regulation and stem cell research.^{1, 2, 3, 4} MEFs are used most commonly as feeder layers to maintain mouse embryonic stem cells in an undifferentiated state. With a combination of transcription factors, MEFs can be converted to a pluripotent state or directly reprogrammed to various types of cells such as functional neurons and cardiomyocytes, which could have important implications for studies of development, neurological and cardiac disease modeling, drug discovery, and regenerative medicine.^{3, 5, 6} The Pierce MEF Isolation Kit also can be used to isolate MEFs from different genetically altered mouse models to study growth control and DNA damage response.⁷ MEFs isolated and cultured with the Pierce MEF Isolation Kit can be readily frozen for future use with minimal loss of viability.

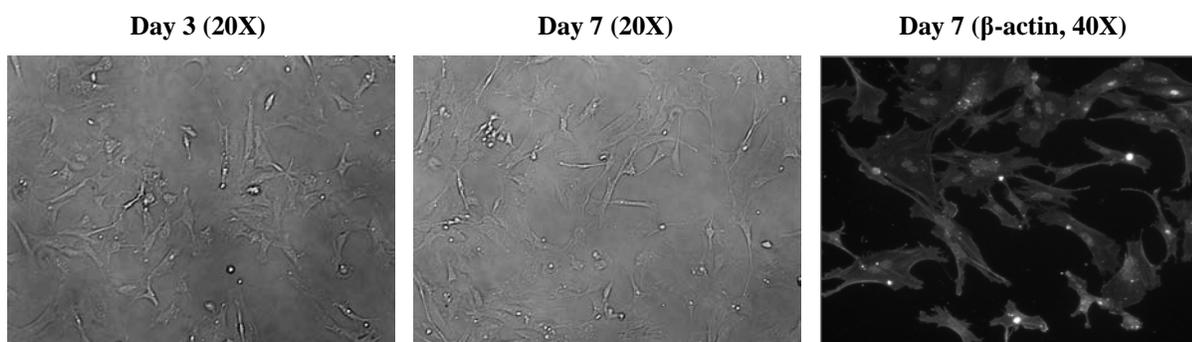


Figure 1. Examples of MEFs in culture following isolation using the Thermo Scientific Pierce Mouse Embryonic Fibroblast Isolation Kit. Phase-contrast and fluorescent images of MEFs after 3 and 7 days in culture. Cultures were plated at a density of 2×10^5 cells per well in a 12-well plate. For the fluorescent image (right panel), MEFs were stained with an antibody specific to β -actin. Images were taken at 20X and 40X magnifications on a Carl Zeiss microscope (AxioVision™ Rel. 4.7).

Important Product Information

- For best cell yield and viability, always isolate MEFs from freshly dissected tissues. The dissection and plating of MEFs should take no longer than two hours.
- Euthanize mice in accordance with the Guide for the Care and Use of Laboratory Animals.⁸
- Perform all tissue digestion, cell manipulations and media handling using sterile technique in a laminar flow cell culture hood to minimize contamination of the isolated MEFs.

Additional Materials Required

- Heat-inactivated fetal bovine serum (FBS) (e.g., Thermo Scientific™ Hyclone™ FBS)
- Penicillin-streptomycin (pen/strep) (e.g., Thermo Scientific™ Hyclone™ Pen/Strep Solution)
- Mouse embryonic tissue freshly dissected from mouse embryos taken from a euthanized mouse (E11-13)
- 37°C heat block or incubator
- Tissue culture incubator at 37°C with humidified, 5% CO₂ atmosphere
- Laminar flow cell culture hood
- Hemocytometer or automated cell counter for counting cells
- Trypan blue stain (e.g., Thermo Scientific™ Hyclone™ Trypan Blue)

Material Preparation

Note: After medium supplementation, medium is stable for approximately one month when stored at 4°C.

Complete DMEM for Primary Cell Isolation Determine the amount of medium needed (see Table 2 for guidelines). In a sterile bottle, add heat-inactivated FBS (10% final concentration), and pen/strep (1% final concentration) to desired volume of DMEM for Primary Cell Isolation. Pre-warm medium to 37°C before use.

Procedure for MEF Isolation

A. Enzyme Digestion of Mouse Embryonic Tissue

Note: For the medium and buffer removal steps, it is critical to carefully remove the medium/buffer contents without disturbing the tissue. For best results, use a pipette and 1000 μ L tip. Do not aspirate using a vacuum flask.

Note: Equilibrate HBSS to 4°C before procedure.

1. Reconstitute the MEF Isolation Enzyme (with papain) by adding 2.5mL of HBSS to one of the vials. Mix gently for 5 minutes or until completely dissolved. Keep enzyme solution on ice.
Note: 2.5mL of reconstituted MEF Isolation Enzyme (with papain) is sufficient for preparing 10 embryonic tissues.
Note: Reconstituted MEF Isolation Enzyme (with papain) can be stored at -20°C for 6 months and is stable for up to two freeze-thaw cycles. The enzyme solution expires one week following preparation if stored at 4°C. Prolonged storage may result in poor performance.
2. Place freshly dissected embryonic tissue into separate 1.5mL sterile microcentrifuge tubes. Immediately add 500µL ice cold HBSS.
Note: For the best results, use one microcentrifuge tube per one embryonic tissue.
3. Mince each embryonic tissue into 1-3mm³ sized pieces. Wash the minced tissue with 500µL ice cold HBSS twice to remove blood from the tissue.
4. Gently remove HBSS using a pipette to the level of the tissue (450µL). Add 0.2mL reconstituted MEF Isolation Enzyme (with Papain) to each tube. Incubate tubes in a 37°C incubator for 25-30 minutes.
5. Gently remove the MEF Isolation Enzyme (with papain) solution and wash tissue twice with 500µL ice cold HBSS.
6. Add 0.5mL pre-warmed Complete DMEM for Primary Cell Isolation to each tube. Break up the tissue by pipetting up and down 15-20 times using a sterile 1.0mL pipette tip. Avoid air bubbles when pipetting.
Note: Disrupting the tissue by pipetting improves cell yield. However, pipetting too vigorously may result in cell damage.
7. After the tissue is primarily a single-cell suspension, add 1.0mL pre-warmed Complete DMEM for Primary Cell Isolation to each tube to bring the total volume to 1.5mL.
8. Combine individual cell suspensions to determine cell concentration and cell viability.

B. Cell Yield and Viability Determination

1. Mix 25µL single cell suspension obtained in Section A, Step 8, with 25µL 0.4% trypan blue in a 1.5mL microcentrifuge tube.
2. Immediately transfer 10µL trypan blue-stained cell suspension to each of two hemocytometer counting chambers.
3. Count both the total number of cells and the number of stained (blue) cells in the hemocytometer microscopic grid.

$$\text{Cell concentration (cells/mL)} = \text{number of cells} \times \text{dilution factor} \times 10^4$$

Example: If 120 cells in a square, then $120 \times 2 \times 10^4 = 2.4 \times 10^6$ cells/mL.

Cell yield = cell concentration \times total volume of cell suspension, obtained in Section A, Step 8.

$$\text{Viability (\%)} = [(\text{total cells counted} - \text{total stained cells}) / \text{total cells counted}] \times 100\%$$

Typical cell yields and viabilities obtained are shown in Table 1.

4. If using an automated cell counter, determine cell yield and viability according to the manufacturer's instructions.

Table 1. Cell yield and viability from a typical isolation.

Cell Type	Yield (cells/mL)	Viability (Trypan Blue Exclusion)
Mouse MEF (one embryonic tissue in 1.5mL cell suspension)	9.0×10^6	90%

C. Plating and Culturing Isolated MEFs

Note: Determine the desired plating density for the cultured MEFs based on the intended downstream study. A density of 5.0×10^4 cells/cm² is generally recommended for culturing MEFs.

1. Pipette the appropriate cell suspension volume to each well of the culture vessel (see Table 2 for recommendations).

Cell suspension volume/well = [required cell density \times growth area (cm²)]/cell concentration (cells/mL from Section B, Step 3)

Example: For a 12-well Nunc™ plate, a single well is approximately 3.5cm². If the cell count is 2.4×10^6 cells/mL, add 83μL of cell suspension to each well.

Table 2. Recommended seeding densities for common culture vessels.*

Nunc Culture Dishes/ Chamber Slides	Well Diameter (mm)	Approximate Growth Area (cm ²)	Media Volume (mL)	Total Number of Cells Required to Seed Each Well (5×10^4 cells/cm ²)
75cm ² flask	NA	75	15	4×10^6
35mm dish	35	9.0	2.0	5×10^5
6-well plate	35	9.6	2.0	5×10^5
12-well plate	22	3.5	0.8	2×10^5
24-well plate	16	1.8	0.5	1×10^5
48-well plate	11	1.1	0.3	5×10^4
96-well plate	4.3	0.14	0.1	1×10^4
4-well chamber slide	NA	1.8	0.5	1×10^5
8-well chamber slide	NA	0.8	0.2	4×10^4

*Additional medium may be required for long-term or less dense cultures.

2. Add Complete DMEM for Primary Cell Isolation to each well to bring the total volume to the recommended level.

Example: For the 12-well Nunc plate example in Step 1, add 717μL Complete DMEM for Primary Cell Isolation to each well to bring the total volume to 0.8mL.

3. Incubate the dishes/chamber slides at 37°C in a 5% CO₂ incubator for 24 hours.
4. After 24 hours, replace the medium with an equivalent volume of fresh Complete DMEM for Primary Cell Isolation.
5. Place the dishes/chambers back into the 37°C, 5% CO₂ incubator.
6. Continue to feed the cultures by replacing the culture medium with fresh Complete DMEM for Primary Cell Isolation every 48-72 hours.

Troubleshooting

Problem	Possible Cause	Solution
Low yield/low viability	Insufficient dissociation	Use freshly reconstituted MEF Isolation Enzyme
		Confirm that the volume of MEF Isolation Enzyme used was the quantity recommended
		Follow recommended incubation time
		Check the expiration date on the product. Do not use after expiration date
	Tissue was not disrupted with pipette	Pipette tissue to aid in dissociation of MEFs
	Over-digestion	Do not exceed an enzyme incubation time of 35 minutes in Section A, Step 4
Ensure that MEF Isolation Enzyme (with papain) is reconstituted to the recommended concentration		
	Disruption by pipette was too vigorous	Do not over-pipette when disrupting digested tissue
Slow cell growth	Medium and/or supplement stored incorrectly, or expired medium and/or supplements	Confirm that the product components were stored at the proper temperature
Cell density too high or too low	Inaccurate total cell numbers seeded in the culture vessels	Use Table 2 to calculate the total cell numbers required for each type of cell culture vessel

Related Thermo Scientific Products

88290	Mouse Embryonic Fibroblast Isolation Enzyme (with papain)
88284	Hanks' Balanced Salt Solution (HBSS, without Ca ²⁺ , Mg ²⁺), 500mL
88287	DMEM for Primary Cell Isolation
88281	Pierce Cardiomyocyte Isolation Kit
88288	Cardiomyocyte Isolation Enzyme 1 (with papain)
88289	Cardiomyocyte Isolation Enzyme 2 (with thermolysin)
88280	Pierce Neuron Isolation Kit
88285	Neuronal Isolation Enzyme (with papain)
88286	Neuronal Culture Media Supplement
88283	Neuronal Culture Medium, 500mL
87790	Subcellular Protein Fractionation Kit for Tissue
78510	T-PER™ Tissue Protein Extraction Reagent

References

1. Wang, S., *et al.* (2012). Inhibition of AMPK α by doxorubicin accentuates genotoxic stress and cell death in mouse embryonic fibroblasts and cardiomyocytes: role of p53 and SIRT1. *J Biol Chem* **287**(11):8001-12.
2. Blackman, B.E., *et al.* (2009). PDE4D and PDE4B function in distinct subcellular compartments in mouse embryonic fibroblasts. *J Biol Chem* **286** (14):12590-601.
3. Vierbuchen, T., *et al.* (2010). Direct conversion of fibroblasts to functional neurons by defined factors. *Nature* **463**:1035-41.
4. Imamura, M., *et al.* (2006). Induction of pluripotent stem cells from mouse embryonic and adult fibroblast cultures by defined factors. *Cell* **126**(4):663-76.
5. Smith, A.W., *et al.* (2013). Direct reprogramming of mouse fibroblasts to cardiomyocyte-like cells using Yamanaka factors on engineered polyethylene glycol (PEG) hydrogels. *Biomaterials* **34**(28):6559-71.
6. Efe, J.A., *et al.* (2011). Conversion of mouse fibroblasts into cardiomyocytes using a direct reprogramming strategy. *Nat Cell Biol* **13**(3):215-22.
7. Lengner, C.J., *et al.* (2004). Primary mouse embryonic fibroblasts: a model of mesenchymal cartilage formation. *J Cell Physiol* **200**(3):327-33.
8. National Research Council of the National Academies (2011). The guide for the care and use of laboratory animals: *The National Academies Press*. Eighth Edition.

Tytuł artykułu

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2016 Thermo Fisher Scientific Inc. All rights reserved. AxioVision is a trademark of Carl Zeiss Corporation. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.