



The following is StemCell Technologies' recommended procedure to isolate mesenchymal progenitor cells from mouse compact bone prior to EasySep® enrichment.

1. Clean a mortar and pestle with 70% isopropanol. Remove the isopropanol from the mortar and pestle, and allow to air dry in a sterile biohazard safety cabinet for 30 minutes. Rinse mortar and pestle with sterile Phosphate Buffered Saline (PBS) just prior to use.
2. Wipe scissors and forceps with an alcohol wipe and air dry. Ensure they are completely dry as residual alcohol may reduce marrow cell viability.
3. After sacrificing the mouse, wet the pelt thoroughly with 70% isopropanol, then clip and peel it back to expose the hind limbs. Using sterile sharp scissors (to avoid splitting of the bone), cut the knee joint in the center and remove ligaments and excess tissue.
4. Remove the femur and tibia by severing them from the mouse at the hip and ankle respectively.
5. Using a scalpel, scrape bones thoroughly to remove muscle, and cut to remove epiphyses. Ensure that the bones are cleaned thoroughly with no remaining muscle tissue attached.
6. Place clean bones in the mortar containing 10 mL PBS with 2% FBS (Catalog #07905) and 1 mM EDTA. The solution of PBS with 2% FBS and 1 mM EDTA is now referred to as 'Buffer'.
7. Crush bones gently with pestle, using only enough force to crack open the bones. Agitate gently to free bone marrow (BM) from bone fragments and pipette Buffer off. Buffer containing BM can be filtered through a 70 µm cell strainer (Falcon Catalog #352350) and used for other applications (for example, culture expansion or to perform the CFU-F assay).
8. Add 10 mL fresh Buffer and repeat agitation and removal of BM. Repeat wash step an additional 4 times (for a total of 6 washes) or until the majority of the BM has been removed (bone fragments turn white in color).

A loss of cell viability and excess debris will be generated when bones are harshly ground. It is important to only use gentle pressure to crack open the bones.

9. Transfer the bone fragments to a 100 mm dish. Add 2 mL of 0.25% Collagenase Type I in PBS containing 20% FBS (Catalog #07902). Ensure bones are completely covered in solution. Let sit for 3 - 5 minutes.

This step softens the bone, allowing it to be chopped more easily.

10. Using a scalpel, chop the remaining bone fragments into fine pieces (1 – 2 mm fragments).

Proper bone fragmentation is required to release sufficient amounts of cells for cell separation.

11. Transfer the bone fragments and collagenase solution to a 50 mL polypropylene tube and add further 0.25% Collagenase Type I (Catalog #07902) to a final volume of 2 mL per mouse used, or a minimum of 10 mL.
12. Seal lid with Parafilm® and place tube in a shaking 37°C waterbath at maximum speed for 45 minutes. If using a bacterial culture shaker, set speed to ~200 rpm.
13. After 45 minutes, remove the tube from the shaker and add Buffer (refer to Step 6) to a final volume of 30 mL. Collect supernatant and filter through a 70 µm cell strainer (Falcon Catalog #352350). Wash bone fragments by mixing with an additional 10 mL of Buffer and allowing fragments to settle for 3 - 4 minutes. Filter the wash through the 70 µm strainer, combining with the previously collected cells (for a final volume of 40 mL).
14. Centrifuge at 300 x g (~1200 rpm) for 10 minutes at room temperature(15-25°C) with the brake on. Remove the supernatant and resuspend the cell pellet in ~200 - 500 µL of medium (note that small particles and debris may be visible in the cell suspension):
 - a. For cell culture, resuspend cells in Complete MesenCult® Medium (Mouse) (Catalog #05511).
 - b. For cell separation experiments, resuspend cells in the recommended medium described in the Mouse Mesenchymal Progenitor Enrichment Kit for Compact Bone Product Information Sheet (Catalog #29122, provided).
Transfer the cells into a smaller tube (e.g. Falcon 5 mL polystyrene round-bottom tubes (BD Catalog #352058).
15. Place cells on ice until ready for use.
16. Remove a small aliquot of cells and dilute 1/20 to 1/100 in 3% Acetic Acid with Methylene Blue (Catalog #07060). Count nucleated cells using a hemacytometer.
17. Expected cell recovery: 1.5 – 3.5 x 10⁶ cells per mouse (2 femurs and 2 tibias). If the cell yield is >5 x 10⁶ cells/mouse, this is an indication that the marrow was not depleted sufficiently.

CFU-F assay and culture expansion protocols are described on the reverse side.

StemCell Technologies

In North America
Tel: 1.604.877.0713
Fax: 1.604.877.0704
Toll Free Tel: 1.800.667.0322
Toll Free Fax: 1.800.567.2899
e-mail: info@stemcell.com
www.stemcell.com

In the United Kingdom
Tel: +44.(0).20.7691.3561
Fax: +33.(0).4.76.18.99.63
Toll Free within United Kingdom:
Tel: 0800.731.27.14
Fax: 0800.731.27.13
e-mail: info@stemcellgb.com

In Europe
Tel: +33.(0).4.76.04.75.30
Fax: +33.(0).4.76.18.99.63
e-mail: info@stemcellfrance.com

November 2007

FOR RESEARCH USE ONLY

#29134

CFU-F Assay for Compact Bone Cells

1. Mouse CFU-F assays initiated from compact bone cells should be initiated at 3 different cell concentrations, to ensure that there are sufficient colony numbers to accurately assess the progenitor cell content of the mouse bone cells.
 - a. For cells from compact bone prepared according to the protocol described on the reverse side and grown in a hypoxic environment (5% O₂/10% CO₂/85% N₂ in a Hypoxia Chamber, Catalog #27310):
 - In 2 mL Complete MesenCult® Medium (Mouse) (Catalog #05511) plate 1,000, 5,000 and 10,000 cells/cm² in duplicate in a 6-well plate (e.g. plate 10,000, 50,000 and 100,000 cells/well).
 - b. For cells from compact bone prepared according to the protocol described on the reverse side and grown in a normal environment (20% O₂/5% CO₂):
 - In 2 mL Complete MesenCult® Medium (Mouse) (Catalog #05511) plate 10,000, 20,000, and 40,000 cells/cm² in duplicate in a 6-well plate (e.g. plate 100,000, 200,000 and 400,000 cells/well).
 - c. For enriched compact bone cells (e.g. enriched for mesenchymal progenitor cells using the EasySep® Mesenchymal Progenitor Enrichment Kit Catalog #19771) grown in a hypoxic environment (5% O₂/10% CO₂/85% N₂ in a Hypoxia Chamber, Catalog #27310):
 - In 2 mL Complete MesenCult® Medium (Mouse) (Catalog #05511) plate 50 - 250 cells/cm² in duplicate in a 6-well plate (e.g. plate 500 - 2,500 cells/well).
 - c. For enriched compact bone cells (e.g. enriched for mesenchymal progenitor cells using the EasySep® Mesenchymal Progenitor Enrichment Kit Catalog #19771) grown in a normal environment (20% O₂/5% CO₂):
 - In 2 mL Complete MesenCult® Medium (Mouse) (Catalog #05511) plate 200 - 1000 cells/cm² in duplicate in a 6-well plate (e.g. plate 2,000 - 10,000 cells/well).
2. Culture cells for 10 - 13 days. No medium change is required.

Cells from compact bone grown in the CFU-F assay occasionally undergo spontaneous differentiation to form adipocytic or chondrogenic cells. For additional information, please contact StemCell Technologies Technical Support (techsupport@stemcell.com).

Expansion Protocol for Cells Isolated from Compact Bone

1. Harvest and process mouse compact bone cells as described in the protocol on the reverse side.
2. For cells grown in a hypoxic environment (5% O₂/10% CO₂/85% N₂ in a Hypoxia Chamber, Catalog #27310):
 - In 2 mL Complete MesenCult® Medium (Mouse) (Catalog #05511) plate 2 - 5 x 10⁵ cells in 1 well of a 6-well plate, or
 - In 10 mL Complete MesenCult® Medium (Mouse) plate 6 x 10⁵ - 1.2 x 10⁶ cells in a T-25 cm² tissue culture flask.

For growth at normal oxygen levels, more cells must be plated. We recommend starting with 5 - 10X more cells, and optimizing for your own experimental conditions.

3. For enriched compact bone cells (e.g. enriched for mesenchymal progenitors using the EasySep® Mouse Mesenchymal Progenitor Enrichment Kit, Catalog #19771) grown in a hypoxic environment (5% O₂/10% CO₂/85% N₂ in a Hypoxia Chamber, Catalog #27310):
 - In 2 mL Complete MesenCult® Medium (Mouse) (Catalog #05511) plate 1 - 5 x 10⁴ cells in 1 well of a 6-well plate, or
 - In 10 mL Complete MesenCult® Medium (Mouse) plate 4 x 10⁴ - 1.2 x 10⁵ cells in a T-25 cm² tissue culture flask.

For growth at normal oxygen levels, more cells must be plated. We recommend starting with 5 - 10X more cells, and optimizing for your own experimental conditions.

4. Grow cells at 37°C in desired atmosphere for 8 - 14 days until an adherent cell layer has formed. After 8 days, if the color of the medium has become orange, a half-medium change can be performed.

For growth at normal oxygen levels, grow cells for 10 - 21 days.

5. Observe mesenchymal cells under the microscope after 7 days to determine confluency. Once the cells have reached 80% confluency, they are ready to be passaged. 80% confluency may be reached in 7 - 16 days.
6. Passage according to the protocol provided in Section 5.1 of the MesenCult® Technical Manual (Mouse; Catalog #28374), Steps 5 - 8.

StemCell Technologies

In North America

Tel: 1.604.877.0713
 Fax: 1.604.877.0704
 Toll Free Tel: 1.800.667.0322
 Toll Free Fax: 1.800.567.2899
 e-mail: info@stemcell.com
www.stemcell.com

In the United Kingdom

Tel: +44.(0).20.7691.3561
 Fax: +33.(0).4.76.18.99.63
 Toll Free within United Kingdom:
 Tel: 0800.731.27.14
 Fax: 0800.731.27.13
 e-mail: info@stemcellgb.com

In Europe

Tel: +33.(0).4.76.04.75.30
 Fax: +33.(0).4.76.18.99.63
 e-mail: info@stemcellfrance.com

November 2007

FOR RESEARCH USE ONLY

#29134