

# StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells

Catalog no. R7789-120

**Version A** 14 November 2008 *A10855* 

## **Table of Contents**

	Contents and Storage	V
	Additional Products	vi
	Introduction	1
M	ethods	4
	General Information	4
	Thawing and Establishing Cells	6
	Subculturing Cells	8
	Freezing Cells	10
	Differentiation Media	13
	Differentiating StemPro® Alk Phos-expressing Rat MSCs	15
Αį	ppendix	19
	Troubleshooting	19
	Technical Support	21
	Purchaser Notification	22
	References	24

## **Contents and Storage**

#### Shipping

StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells are shipped on dry ice.

## Contents and Storage

Contents and storage conditions for StemPro® Alk Phosexpressing Rat Mesenchymal Stem Cells are listed in the table below. For components of the freezing medium, see page 11.

Product	Amount	Storage
StemPro® Alk Phos-expressing Rat	1 ml	Liquid nitrogen
Mesenchymal Stem Cells (1 × 10 <sup>6</sup> cells/ml in		
freezing medium)		



Handle cells as potentially biohazardous material under at least Biosafety Level 1 (BL-1) containment. This product contains Dimethyl Sulfoxide (DMSO), a hazardous material. Review the Material Safety Data Sheet (MSDS) before handling. Material Safety Data Sheets (MSDSs) are available on our website at www.invitrogen.com/msds.

# Information for European Customers

StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells are genetically modified and carry a chromosomal human Alkaline Phosphatase gene. As a condition of sale, this product must be in accordance with all applicable local legislation and guidelines including EC Directive 90/219/EEC on the contained use of genetically modified organisms.

## **Additional Products**

## Additional Products

The products listed in this section may be used with StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells. For more information, refer to our website (<a href="www.invitrogen.com">www.invitrogen.com</a>) or contact Technical Support (see page 21).

Item	Quantity	Cat. no.
Minimum Essential Medium (MEM) $\alpha$ Medium (1X) with GlutaMAX <sup>TM</sup> -I, ribonucleosides and deoxyribonucleosides	500 ml	32571-036
GlutaMAX™-I Supplement	100 ml	35050-061
Fetal Bovine Serum (FBS), MSC-Qualified	100 ml 500 ml	12662-011 12662-029
StemPro® Adipogenesis Differentiation Kit	100 ml	A10070-01
StemPro® Chondrogenesis Differentiation Kit	100 ml	A10071-01
StemPro® Osteogenesis Differentiation Kit	100 ml	A10072-01
Gentamicin (10 mg/ml)	10 ml	15710-064
Dulbecco's Phosphate Buffered Saline (DPBS), containing no calcium, magnesium, or phenol red	500 ml	14190-144
TrypLE™ Express Dissociation Enzyme without Phenol Red	100 ml 20 × 100 ml	12604-013 12604-039
Antibiotic-Antimycotic (100X), liquid	100 ml	15240-062
Gentamycin Reagent Solution (10 mg/ml), liquid	10 ml	15710-064
Gentamycin Reagent Solution (50 mg/ml), liquid	10 ml	15750-060
Trypan Blue Stain	100 ml	15250-061
LIVE/DEAD® Cell Vitality Assay Kit	1000 assays	L34951
Countess™ Automated Cell Counter (includes 50 Countess™ cell counting chamber slides and 2 ml of Trypan Blue Stain)	1 unit	C10227
ELF® 97 Endogenous Phosphatase Detection Kit	1 kit	E6601
CultureWell™ chambered coverglass (16 wells per coverglass, set of 8)	1 set	C37000

## Introduction

#### Introduction

StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells (MSCs) are produced from bone marrow isolated from transgenic Fischer 344 rats expressing the human placental alkaline phosphatase (hPAP) gene linked to the ubiquitously active ROSA26 (R26) gene promoter (Kisseberth *et al.*, 1999; Mujtaba *et al.*, 2002). The cells were isolated under sterile conditions and cryopreserved from primary cultures. Before cryopreservation, the MSCs are expanded for three passages in  $\alpha$ -MEM medium supplemented with 10% MSC-Qualified FBS and antibiotic/antimycotic solution. The freezing medium consisted of 70%  $\alpha$ -MEM, 20% MSC-Qualified FBS, and 10% DMSO.

Each vial of MSCs contains cells that can differentiate into multiple mature cell phenotypes in vitro, including adipocytes, osteocytes, and chondrocytes (De Ugarte et al., 2003; Meirelles Lda & Nardi, 2003; Pittenger et al., 1999; Wu et al., 2002). In vitro differentiation into non-mesenchymal cell types, such as neuronal and myogenic cells have also been described (Anjos-Afonso et al., 2004; Deng et al., 2001; Han et al., 2002; Han et al., 2004; Moscoso et al., 2005; Phinney et al., 1999; Wakitani et al., 1995). In addition, MSCs are shown to be involved in certain types of cancers (Houghton et al., 2004; Singh et al., 2004), and are known to secrete immunomodulatory, anti-angiogenic, anti-inflammatory, pro-cardiovasculogenic, and pro-arteriogenic factors (Djouad et al., 2003; Gojo et al., 2003; Houghton et al., 2004; Kinnaird et al., 2004; Krampera et al., 2003; Oh et al., 2008; Olivares et al., 2004; Orlic et al., 2001).

StemPro® Alk Phos-expressing Rat MSCs can be used for studies of adult stem cell differentiation, tissue engineering, cell and genetic therapy, and potential future clinical applications. These cells can also be used in transplant studies to track transplanted cells as they differentiate into mature phenotypes.

We recommend that you use  $\alpha$ -MEM with GlutaMAX<sup>TM</sup>-I and MSC-Qualified FBS (see page vi) for optimal growth and expansion.

## Introduction, continued

### Characteristics of StemPro® Alk Phos-Expressing MSCs

- Are prepared from low-passage (passage 3) adherent rat primary cell cultures
- Express a flow-cytometry cell-surface protein profile positive for CD29, CD73, and CD90 (> 70%), and negative for CD45 (< 10%)</li>
- Stain positive for Alkaline Phosphatase (> 80%)
- Contain cells characteristic of at least tri-potential differentiation that can differentiate into osteogenic, adipogenic, and chondrogenic lineages.

## Isolation and Expansion

StemPro® Alk Phos-expressing Rat MSCs are extracted from the hind leg bones of alkaline phosphatase transgenic Fischer 344 rats through mechanical and enzymatic digestion. Cells are expanded using  $\alpha\textsc{-MEM}$  medium supplemented with 10% MSC-Qualified FBS and antibiotic/antimycotic solution, which supports a cell doubling time of 30  $\pm$  5 hours.

The *in vitro* growth capacity of MSCs has not been definitely established and can vary greatly depending on the culture conditions such as seeding density and growth factors used, but the cells can be expected to expand for at least 30 population doublings before their growth rate decreases significantly (Bruder *et al.*, 1997; Meirelles Lda & Nardi, 2003).

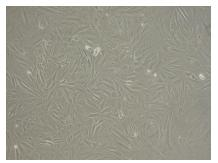
## Differentiation Potential

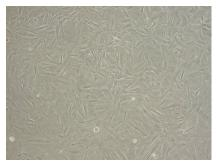
Multiple investigators have demonstrated that MSCs can be differentiated towards multiple mature cell phenotypes. In addition to traditional mesenchymal lineages, MSCs have been differentiated towards cardiomyocytic and neuronal phenotypes using specialized media. The *in vitro* differentiation potential of MSCs has not been definitely established, but long-term culture and high cell density are implicated in the loss of differentiation potential (Meirelles Lda & Nardi, 2003).

## Introduction, continued

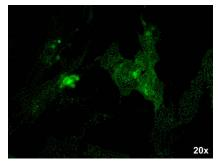
### Alkaline Phosphatase Expression

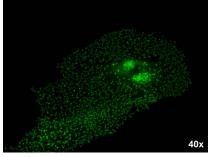
*In vivo* tracking of implanted MSCs in cell and gene therapy protocols is very important as the success of these therapies depends on MSCs' engraftment abilities, especially after systemic infusion (Meirelles Lda & Nardi, 2003). Further, it has been shown that MSCs can fuse with other cells and acquire their characteristics (Spees *et al.*, 2003). The StemPro® Alk Phos-expressing Rat MSCs allow the user to track the implanted cells (Han *et al.*, 2002; Han *et al.*, 2004; Mujtaba *et al.*, 2002) using a simple, fluorescence-based enzymatic assay, where the removal of the phosphate from the substrate provided in the assay kit causes an intense yellow-green fluorescence (ELF® 97 Endogenous Phosphatase Detection Kit, see page vi for ordering information).





Bright field images (10X) of StemPro $^{\circ}$  Alk Phos-expressing Rat MSCs at P4 that have been in culture for 14 days.





Fluorescence images (20X and 40X) of StemPro® Alk Phos-expressing Rat MSCs at P4 that have been in culture for 5 days. Alkaline phosphatase expression is detected using the ELF® 97 Endogenous Phosphatase Detection Kit

## **Methods**

## **General Information**

### General Cell Handling

Follow the general guidelines below to grow and maintain StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells.

- 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.
- Before starting experiments, ensure cells have been established (at least 1 passage), and also have some frozen stocks on hand.
- For differentiation studies and other experiments, we recommend using cells below passage 5.
- For general maintenance of cells, cell confluency should be 60–80%, cell viability should be at least 90%, and the growth rate should be in mid-logarithmic phase prior to subculturing.
- When thawing or subculturing cells, transfer cells into pre-warmed medium.
- Antibiotic-antimycotic containing penicillin, streptomycin, and amphotericin B may be used if required (see page vi for ordering information).



As with other mammalian cell lines, when working with MSCs, handle as potentially biohazardous material under at least Biosafety Level 1 (BL-1) containment. For more information on BL-1 guidelines, refer to *Biosafety in Microbiological and Biomedical Laboratories*, 4<sup>th</sup> ed., published by the Centers for Disease Control, or see the following website: <a href="https://www.cdc.gov/od/ohs/biosfty/bmbl4/bmbl4toc.htm">www.cdc.gov/od/ohs/biosfty/bmbl4/bmbl4toc.htm</a>



It is very important to strictly follow the guidelines for culturing StemPro® Alk Phos-expressing Rat Mesenchymal Stem Cells in this manual to keep them undifferentiated.

## General Information, continued

### Media Requirements

We recommend using Minimum Essential Medium (MEM)  $\alpha$  medium ( $\alpha$ -MEM medium) with GlutaMAX $^{\text{\tiny M}}$ -I and supplemented with 10% MSC-Qualified Fetal Bovine Serum (FBS) for optimal growth and expansion of StemPro $^{\text{\tiny B}}$  Alk Phos-expressing Rat MSCs, and to keep them undifferentiated (see page vi for ordering information).

- Prepare your growth medium prior to use.
- When thawing or subculturing cells, transfer cells into pre-warmed medium at 37°C.
- You may store the complete growth medium in the dark at 4°C for up to four weeks.
- Avoid repeated freeze-thaw cycles of MSC-Qualified FBS.



We have observed that StemPro® Alk Phos-expressing Rat MSCs adhere poorly when plated on media other than  $\alpha$ -MEM medium supplemented with 10% **MSC-Qualified** FBS after their initial thaw. Although they recover and adhere well after their first passage, we suggest that you use the recommended media. If you prefer to culture your MSCs on growth media other than the recommended, we advise you to optimize your growth conditions and treat the your cells gently (*i.e.*, do not vortex, bang the flasks to dislodge the cells, or centrifuge the cells at high speeds).

## **Thawing and Establishing Cells**

#### Introduction

To thaw StemPro® Alk Phos-expressing Rat MSCs and to initiate cell culture, follow the protocol below.

#### Materials Needed

The following materials are required (see page vi for ordering information).

- StemPro® Alk Phos-expressing Rat MSCs, stored in liquid nitrogen
- Ethanol or 70% isopropanol
- α-MEM medium with GlutaMAX<sup>™</sup>-I containing 10% MSC-Qualified FBS plus antibiotic/antimycotic or gentamycin; pre-warmed to 37°C
- Disposable, sterile 15-ml tubes
- 37°C water bath
- 37°C incubator with a humidified atmosphere of 5% CO<sub>2</sub>
- Microcentrifuge
- Tissue-culture treated flasks, plates or dishes
- Hemacytometer, cell counter and Trypan Blue, LIVE/DEAD® Cell Vitality Assay Kit, or the Countess™ Automated Cell Counter



Invitrogen's Countess™ Automated Cell Counter is a benchtop counter designed to measure cell count and viability (live, dead, and total cells) accurately and precisely in less than a minute per sample, using the standard Trypan Blue technique (see page vi for ordering information).

Using the same amount of sample that you currently use with the hemocytometer, the Countess™ Automated Cell Counter takes less than a minute per sample for a typical cell count and is compatible with a wide variety of eukaryotic cells and provides information on cell size.

## Thawing and Establishing Cells, continued

# Thawing Procedure, continued

To thaw and establish StemPro® Alk Phos-expressing Rat MSCs:

- Pre-warm the prepared α-MEM medium with GlutaMAX<sup>™</sup>-I containing 10% MSC-Qualified FBS and antibiotic/antimycotic or gentamycin to 37°C.
- Remove the cells from liquid nitrogen storage, and wipe the cryovial with ethanol or 70% isopropanol before opening. In an aseptic field, briefly twist the cap a quarter turn to relieve pressure and then retighten. Do not expose cells to air before thawing.
- Quickly thaw the vial of cells by swirling it in a 37°C water bath and removing it when the last bit of ice has melted, typically < 2 minutes. Do not submerge the vial completely. Do not thaw the cells for longer than 2 minutes.</li>
- 4. When thawed, immediately transfer cells into a 15-ml sterile tube and add pre-warmed complete  $\alpha$ -MEM medium **dropwise** up to 10 ml.
- 5. Centrifuge cells for 5 minutes at  $300 \times g$ .
- 6. Aspirate supernatant and resuspend cells in 2 ml of complete  $\alpha$ -MEM medium
- 7. Determine the viable cell count using your method of choice, and plate the resuspended cells at a seeding density of 5,000 cells per cm². If necessary, add complete α-MEM medium to the cells to achieve the desired cell concentration and recount the cells.
- 8. Incubate at 37°C, 5% CO<sub>2</sub> and 90% humidity and allow cells to adhere for at least 24 hours.
- 9. The next day, replace the medium with an equal volume of fresh, pre-warmed complete  $\alpha$ -MEM medium.
- 10. Change the medium every 3-4 days.

## **Subculturing Cells**

#### Introduction

Follow the protocol below to culture StemPro® Alk Phosexpressing Rat MSCs. Subculture cells when needed (before colonies start contacting each other), typically every 7–10 days.

#### Materials Needed

The following materials are required (see page vi for ordering information).

- Culture vessels containing StemPro® Alk Phosexpressing Rat MSCs
- Tissue-culture treated flasks, plates or dishes
- α-MEM medium with GlutaMAX<sup>™</sup>-I supplemented with 10% MSC-Qualified FBS and containing antibiotic/antimycotic or gentamycin, pre-warmed to 37°C
- Disposable, sterile 50-ml tubes
- 37°C incubator with humidified atmosphere of 5% CO<sub>2</sub>
- Dulbecco's Phosphate Buffered Saline (DPBS), containing no calcium, magnesium, or phenol red
- TrypLE<sup>™</sup> Express, pre-warmed to 37°C
- Hemacytometer, cell counter and Trypan Blue, LIVE/DEAD® Cell Vitality Assay Kit, or the Countess™ Automated Cell Counter

## Passaging Cells

- 1. Aspirate the complete  $\alpha$ -MEM medium from the cells.
- Rinse the surface of the cell layer with DPBS without Ca<sup>2+</sup> and Mg<sup>2+</sup> (approximately 2 ml DPBS per 10 cm<sup>2</sup> culture surface area) by adding the DPBS to the side of the vessel opposite the attached cell layer, and rocking back and forth several times.
- 3. Aspirate the DPBS and discard.
- To detach the cells, add a sufficient volume of pre-warmed TrypLE<sup>™</sup> Express to cover the cell layer (approximately 0.5 ml/10 cm²).
- 5. Incubate at 37°C for approximately 5–8 minutes.

Procedure continued on next page

## Subculturing Cells, continued

# Passaging Cells, continued

Procedure continued from previous page

- 6. Observe the cells under a microscope. If the cells are less than 90% detached, continue incubating and observe within 2 minutes for complete detachment of the cells. Tap the vessel gently to expedite cell detachment.
- 7. When ≥ 90% of the cells have detached, tilt the vessel for a minimal length of time to allow the cells to drain. Add the equivalent of 2 volumes (twice the volume used for TrypLE™ Express) of pre-warmed complete α-MEM medium. Disperse the medium by pipetting over the cell layer surface several times.
- 8. Transfer the cells to a 50-ml conical tube and centrifuge at  $300 \times g$  for 5 minutes at room temperature. Aspirate and discard the medium
- 9. Resuspend the cell pellet in a minimal volume of pre-warmed complete  $\alpha$ -MEM medium and remove a sample for counting.
- 10. Determine the total number of cells and percent viability using your method of choice. If necessary, add complete  $\alpha$ -MEM medium to the cells to achieve the desired cell concentration and recount the cells.
- Determine the total number of vessels to inoculate by using the following equation:
   Number of vessels = Number of viable cells ÷ (growth
  - area of vessel in  $cm^2 \times 5,000$  cells per  $cm^2$  recommended seeding density)
- 12. Add complete  $\alpha$ -MEM medium to each vessel so that the final culture volume is 0.2–0.5 ml per cm<sup>2</sup>.
- 13. Add the appropriate volume of cells to each vessel and incubate at 37°C, 5% CO<sub>2</sub> and 90% humidity.
- 14. 3–4 days after seeding, completely remove the medium. Replace with an equal volume of complete  $\alpha$ -MEM medium.

## **Freezing Cells**

#### Introduction

Guidelines and procedures for preparing freezing medium and freezing cells are provided in this section.

#### Materials Needed

The following materials are required (see page vi for ordering information).

- Culture vessels containing StemPro<sup>®</sup> Alk Phosexpressing Rat MSCs
- α-MEM medium
- Fetal Bovine Serum, MSC-Qualified
- DMSO (use a bottle set aside for cell culture; open only in a laminar flow hood)
- Disposable, sterile 15-ml conical tubes.
- DPBS, containing no calcium, magnesium, or phenol red
- TrypLE<sup>™</sup> Express
- Hemacytometer, cell counter and Trypan Blue, LIVE/DEAD® Cell Vitality Assay Kit, or the Countess™ Automated Cell Counter
- Sterile freezing vials

#### Guidelines

When freezing MSCs, we recommend the following:

- Freeze cells at a density of  $1-2 \times 10^6$  viable cells/ml.
- Use a freezing medium composed of final concentrations of 20% MSC-Qualified FBS and 10% DMSO.
- Bring the cells into freezing medium in two steps, as described in this section.

## Freezing Cells, continued

### Preparing Freezing Media

Prepare Freezing Medium A and B immediately before use. You will need enough of each freezing medium to resuspend cells at a density of  $1-2 \times 10^6$  cells/ml (see the freezing procedure below).

 In a sterile 15-ml tube, mix together the following reagents for every 1 ml of Freezing Medium A needed:

 $\alpha$ -MEM medium 0.6 ml FBS, MSC-Qualified 0.4 ml

2. In another sterile 15-ml tube, mix together the following reagents for every 1 ml of **Freezing Medium B** needed:

 $\alpha$ -MEM medium 0.8 ml DMSO 0.2 ml

Place tube with Freezing Medium B on ice until use (leave Freezing Medium A at room temperature).

**Note:** Discard any remaining freezing medium after use.

## Freezing Cells Procedure

- 1. Aspirate complete  $\alpha$ -MEM medium from the flask, well, or dish.
- Rinse the surface with DPBS without Ca<sup>2+</sup> and Mg<sup>2+</sup> (approximately 2 ml DPBS per 10 cm<sup>2</sup> culture surface area) by adding the DPBS to the side of the vessel opposite the attached cell layer and rocking back and forth several times.
- 3. Aspirate the DPBS and discard.
- 4. To detach the cells, add a sufficient volume of pre-warmed TrypLE<sup>™</sup> Express to cover the cell layer (approximately 0.5 ml/10 cm²).
- 5. Incubate at 37°C for approximately 5–8 minutes.
- 6. Observe the cells under a microscope. If the cells are less than 90% detached, continue incubating and observe within 2 minutes for complete detachment of the cells. Gently tap the vessel to expedite cell detachment.

Procedure continued on next page

## Freezing Cells, continued

#### Freezing Cells Procedure, continued

Procedure continued from previous page

- 7. When ≥90% of the cells have detached, tilt the vessels on end for a minimal length of time to allow the cells to drain. Add the equivalent of 2 volumes (twice the volume used for the TrypLE™ Express) of pre-warmed complete α-MEM medium to each vessel. Disperse the medium by pipetting over the cell layer surface several times.
- 8. Transfer the cells to a 15-ml conical tube and centrifuge at  $300 \times g$  for 5 minutes at room temperature. Aspirate the medium used for washing the cells (step 7).
- 9. Resuspend the cell pellet in a minimal volume of pre-warmed complete  $\alpha$ -MEM medium and remove a sample for counting.
- Determine the total number of cells using your method of choice.
- 11. Gently aspirate media from the vessel and resuspend the cells to a concentration of  $4\times10^6$  cells/ml in Freezing Medium A.
- 12. Add the same volume of Freezing Medium B to cells in a **dropwise** manner.
- 13. Aliquot 1 ml to each freezing vial and store at -80°C overnight in an isopropanol chamber.
- 14. The next day, transfer the frozen vials to a liquid nitrogen tank for long-term storage.

**Note:** You may check the viability and recovery of frozen cells 24 hours after storing cryovials in liquid nitrogen by following the procedure outlined in **Thawing and Establishing Cells**, page 6.

## **Differentiation Media**

#### Introduction

One critical hallmark of MSCs is their ability to differentiate into three or more mature cell types. Traditional and modern bioassays are used to demonstrate the multipotency of MSCs to differentiate along the osteogenic, adipogenic, and chondrogenic lineages. This section provides guidelines for preparing media that are used for inducing StemPro® Alk Phos-expressing Rat MSCs to differentiate into osteogenic, adipogenic and chondrogenic cell types.

#### Mesenchymal Stem Cell Basal Medium

MSC basal medium is used a cell attachment medium and as a negative control during differentiation experiments. It consists of  $\alpha\textsc{-MEM}$  medium with GlutaMAX $^{\!\scriptscriptstyle{\text{\tiny M}}}\textsc{-I}$  containing 10% MSC-Qualified FBS and 5  $\mu\textsc{I}/\text{ml}$  gentamicin (see page vi).

Component	Final Conc.	For 500 ml
α-MEM medium with GlutaMAX™-I	1X	450 ml
FBS, MSC-Qualified	10%	50 ml
Gentamicin (10 mg/ml)	5 μg/ml	250 µl

# Osteogenic Differentiation Medium

To prepare osteogenic differentiation (OD) medium, combine the following in a sterile flask. Although you may use the StemPro® Osteocyte/Chondrocyte Differentiation Basal Media, differentiation appears to be more efficient with  $\alpha\text{-MEM}$  as the basal media. Store the OD medium at 4°C in the dark up to four weeks.

Component	Final Conc.	For 100 ml
α-MEM medium with GlutaMAX™-I	1X	90 ml
StemPro® Osteogenesis Supplement	1X	10 ml
Gentamicin (10 mg/ml)	5 µg/ml	50 µl

## **Differentiation Media**, continued

### Adipogenic Differentiation Medium

To prepare adipogenic differentiation (AD) medium, combine the following in a sterile flask. Although you may use the StemPro® Adipocyte Differentiation Basal Media, differentiation appears to be more efficient with  $\alpha$ -MEM as the basal media. Store the AD medium at 4°C in the dark up to four weeks.

Component	Final Conc.	For 100 ml
$\alpha$ -MEM medium with GlutaMAX $^{\text{\tiny TM}}$ -I	1X	90 ml
StemPro® Adipogenesis Supplement	1X	10 ml
Gentamicin (10 mg/ml)	5 µg/ml	50 µl

# Chondrogenic Differentiation Medium

To prepare chondrogenic differentiation (CD) medium, combine the following in a sterile flask. Although you may use the StemPro® Osteocyte/Chondrocyte Differentiation Basal Media, differentiation appears to be more efficient with  $\alpha\textsc{-MEM}$  as the basal media. Store the CD medium at 4°C in the dark up to four weeks.

Component	Final Conc.	For 100 ml
α-MEM medium with GlutaMAX™-I	1X	90 ml
StemPro® Chondrogenesis Supplement	1X	10 ml
Gentamicin (10 mg/ml)	5 µg/ml	50 µl

## Differentiating StemPro® Alk Phos-expressing Rat MSCs

#### Introduction

This section provides guidelines and instructions for inducing StemPro® Alk Phos-expressing Rat MSCs to differentiate into osteogenic, adipogenic, and chondrogenic cell types.

#### Materials Needed

The following materials are required (see page vi for ordering information).

- Culture vessels containing your MSCs
- Tissue-culture treated flasks, plates, or dishes
- MSC Basal Medium, prewarmed to 37°C (see page 13)
- Appropriate Differentiation Medium, pre-warmed to 37°C (see pages 13–14)
- Dulbecco's Phosphate Buffered Saline (DPBS), containing no calcium, magnesium, or phenol red
- Disposable, sterile 50-ml tubes
- 37°C incubator with humidified atmosphere of 5% CO<sub>2</sub>
- TrypLE<sup>™</sup> Express, pre-warmed to 37°C
- Hemacytometer, cell counter and Trypan Blue, LIVE/DEAD<sup>®</sup> Cell Vitality Assay Kit, or the Countess<sup>™</sup> Automated Cell Counter

## Harvesting MSCs

Follow the protocol below to harvest your StemPro® Alk Phosexpressing Rat MSCs for differentiation experiments. We recommend that you expand your cells to  $\leq 70\%$  confluency in a tissue-culture treated T-225 flask, and prepare the appropriate differentiation medium ahead of time.

- 1. Aspirate complete  $\alpha$ -MEM medium from the flask and rinse the surface with DPBS without Ca<sup>2+</sup> and Mg<sup>2+</sup> (approximately 2 ml DPBS per 10 cm<sup>2</sup> culture surface area) by adding the DPBS to the side of the vessel opposite the attached cell layer and rocking back and forth several times.
- 2. Aspirate the DPBS and discard.
- To detach the cells, add a sufficient volume of pre-warmed TrypLE<sup>™</sup> Express to cover the cell layer (approx. 0.5 ml/10 cm²).
- 4. Incubate at 37°C for approximately 5–8 minutes.

Procedure continued on next page

# Differentiating StemPro® Alk Phos-expressing Rat MSCs, continued

# Harvesting MSCs, continued

Procedure continued from previous page

- 5. Observe the cells under a microscope. If the cells are less than 90% detached, continue incubating and observe within 2 minutes for complete detachment of the cells. Gently tap the vessel to expedite cell detachment.
- 6. Spin for 5 minutes at  $300 \times g$  at room temperature. While the cells are spinning, perform a viable cell count using your method of choice; note total cell number. Calculate required amount of MSC basal medium to obtain the appropriate seeding concentration (see below).
- Resuspend cells in the appropriate amount of MSC basal medium.
- 8. Dispense cell solution according to differentiation condition being tested (see protocols below).

#### Osteogenic Differentiation Protocol

Follow the protocol below to differentiate your StemPro® Alk Phos-expressing Rat MSCs into an osteogenic phenotype.

- Seed the MSCs into culture vessels at 1.9 × 10<sup>4</sup> cells/cm<sup>2</sup>.
   For classical stain differentiation assays, seed into a
   12-well plate. For gene-expression profile studies, seed into a T-75 flask. For immunocytochemistry studies, seed into a 16-well CultureWell™ chambered coverglass or 96-well plate.
- 2. To six wells of a 12-well plate, add 1 ml of cell solution per well and let attach in the 37°C, 5% CO<sub>2</sub> incubator for a minimum of two hours.
- 3. Replace three wells with MSC basal medium as negative controls, and other three wells with fresh OD medium. Let culture at 37°C with 5% CO<sub>2</sub>.
- 4. Refeed cultures every 2–3 days with media prepared at initiation of differentiation. MSCs will continue to expand as they differentiate under osteogenic conditions.
- 5. After specific periods of cultivation, osteogenic cultures can be processed for alkaline phosphatase staining (7–14 days) or Alizarin Red S staining (>21 days), gene expression analysis, or protein detection. For long term culture (>21 days), we recommend that you reduce the seeding density by half  $(9.5 \times 10^3 \text{ cells/cm}^2)$  to prevent overgrowth and cell detachment.

# Differentiating StemPro® Alk Phos-expressing Rat MSCs, continued

### Adipogenic Differentiation Protocol

Follow the protocol below to differentiate your StemPro® Alk Phos-expressing Rat MSCs into an adipogenic phenotype.

- Seed the MSCs into culture vessels at 7.6 × 10<sup>4</sup> cells/cm<sup>2</sup>.
   For classical stain differentiation assays, seed into a
   12-well plate. For gene-expression profile studies, seed into a T-75 flask. For immunocytochemistry studies, seed into a 16-well CultureWell™ chambered coverglass or 96-well plate.
- 2. To six wells of a 12-well plate, add 1 ml of cell solution per well, and let attach in the 37°C, 5% CO<sub>2</sub> incubator for a minimum of two hours.
- Replace three wells with MSC basal medium as negative controls, and other three wells with fresh AD medium. Let culture at 37°C and 5% CO<sub>2</sub>.
- Refeed cultures every 3–4 days with media prepared at initiation of differentiation. MSCs will continue to undergo limited expansion as they differentiate under adipogenic conditions.
- 5. After specific periods of cultivation, adipogenic cultures can be processed for Oil Red O or LipidTOX™ staining (beginning at 7–14 days), gene expression analysis, or protein detection.

# Differentiating StemPro® Alk Phos-expressing Rat MSCs, continued

### Chondrogenic Differentiation Protocol

Follow the protocol below to differentiate your StemPro® Alk Phos-expressing Rat MSCs into a chondrogenic phenotype.

- Detach cells using TrypLE™ Express and perform a cell count as described in Harvesting MSCs, pages 15–16 (through Step 6).
- 2. Resuspend the cells in MSC basal medium to a concentration of  $8 \times 10^6$  cells/ml.
- 3. To six wells in a 12-well tissue-culture dish, spot 10  $\mu$ l of cells per well.
- 4. Incubate for two hours at  $37^{\circ}$ C, 5% CO<sub>2</sub> and 90% humidity.

**Note**: If this step is not performed under high humidity conditions, the spots may dehydrate and the formation of chondrogenic pellets inhibited.

- To three of the spotted wells, add 1 ml of MSC basal medium as a negative control. To the other three wells, add 1 ml of CD medium.
- 6. Incubate at 37°C, 5% CO<sub>2</sub>, and 90% humidity. Refeed cultures every 2–3 days with same media, prepared at the initiation of differentiation.
- Check for chondrogenesis after a set period of cultivation. You may perform alcian blue staining on the pellets (to detect glycosaminoglycans) after 14 days, or paraffin section of pellets for collagen 2a immunohistological staining after ~21 days.



## **Appendix**

## **Troubleshooting**

## Culturing Cells

The table below lists some potential problems and solutions that help you troubleshoot your cell culture problems.

Problem	Cause	Solution
No viable cells after thawing stock	Stock not stored correctly	Order new stock and store in liquid nitrogen. Keep in liquid nitrogen until thawing.
	Home-made stock not viable	Freeze cells at a density of $1-2 \times 10^6$ viable cells/ml.
		Use low-passage cells to make your own stocks.
		Follow procedures in <b>Freezing Cells</b> (page 10) exactly. Slow freezing and fast thawing is the key. Add Freezing Medium B drop wise manner (slowly). At time of thawing, thaw quickly and do not expose vial to the air but quickly change from nitrogen tank to 37°C water bath.
		Obtain new StemPro® Alk Phos-expressing Rat MSCs.
	Thawing medium not correct	Use pre-warmed complete $\alpha$ -MEM medium, prepared as described on page 5. Be sure to use MSC-Qualified FBS.
	Cells too diluted	Generally we recommend thawing one vial at a density of 5,000 cells per cm <sup>2</sup> .
	Cell not handled gently.	StemPro® Alk Phos-expressing Rat MSCs are fragile; treat your cells gently, do not vortex, bang the flasks to dislodge the cells, or centrifuge the cells at high speeds.
Cells grow slowly	Growth medium not correct	Use prewarmed complete $\alpha$ -MEM medium.
	Cells too old	Use healthy MSCs, under passage 5; do not overgrow.
Cells differentiated	Culture conditions not correct	Thaw and culture fresh vial of new StemPro® Alk Phos-expressing Rat MSCs. Follow thawing instructions (page 6) and subculture procedures (page 8) exactly.
	Cells too old	MSCs above passage 5 may become differentiated.

## Troubleshooting, continued

### Culturing Cells, continued

The table below lists some potential problems and solutions that help you troubleshoot your cell culture problems.

Problem	Cause	Solution
Cells not adherent after initial thaw	Used serum other than MSC- Qualified FBS	Be sure to prepare your culture medium using <b>MSC-Qualified</b> FBS (see page vi for ordering information).
Cannot detect expression of alkaline phosphatase	Assay system not sensitive enough	Use ELF® 97 Endogenous Phosphatase Detection Kit (see page vi for ordering information).

### Differentiating Cells

The table below lists some potential problems and solutions that help you troubleshoot your cell culture problems.

Problem	Cause	Solution
Cells fail to differentiate	Used StemPro® Osteocyte/Chon- drocyte or Adipocyte Differentiation Basal Media	Although you may use the StemPro® Osteocyte/Chondrocyte or Adipocyte Differentiation Basal Media for your differentiation studies, we have observed that differentiation is more efficient with $\alpha$ -MEM as the basal media. Repeat your differentiation studies using $\alpha$ -MEM as the basal media
	Initial spotting step not performed under high humidity (if differentiating into chondrocytes)	If this step is not performed under high humidity conditions, the spots may dehydrate and the formation of chondrogenic plates inhibited. Repeat the initial spotting step at 37°C, 5% CO <sub>2</sub> , and 90% humidity.
Cells have overgrown the culture plates and have detached	Initial seeding density too high	For long term culture (>21 days), we recommend that you seed at a lower cell density of $3 \times 10^3$ cells/cm <sup>2</sup> to prevent overgrowth and cell detachment.

## **Technical Support**

#### Web Resources



Visit the Invitrogen website at <u>www.invitrogen.com</u> for:

- Technical resources, including manuals, vector maps and sequences, application notes, MSDSs, FAQs, formulations, citations, handbooks, etc.
- Complete Technical Support contact information
- Access to the Invitrogen Online Catalog
- Additional product information and special offers

#### **Contact Us**

For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our website (<a href="www.invitrogen.com">www.invitrogen.com</a>).

### Corporate Headquarters:

Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 USA Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288

Fax: 1 760 602 6500

E-mail:

tech\_support@invitrogen.com

#### Japanese Headquarters:

Invitrogen Japan LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Tel: 81 3 5730 6509 Fax: 81 3 5730 6519

E-mail: jpinfo@invitrogen.com

#### **European Headquarters:**

Invitrogen Ltd Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF, UK Tel: 44 (0) 141 814 6100 Tech Fax: 44 (0) 141 814 6117 E-mail:

E-mail: eurotech@invitrogen.com

### Material Safety Data Sheets (MSDSs)

MSDSs (Material Safety Data Sheets) are available on our website at <a href="https://www.invitrogen.com/msds">www.invitrogen.com/msds</a>.

### Certificate of Analysis

The Certificate of Analysis provides detailed quality control information for each product. Certificates of Analysis are available on our website. Go to <a href="https://www.invitrogen.com/support">www.invitrogen.com/support</a> and search for the Certificate

of Analysis by product lot number, which is printed on the box.

## **Purchaser Notification**

## Limited Warranty

Invitrogen is committed to providing our customers with highquality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, contact our Technical Support Representatives.

Invitrogen warrants that all of its products will perform according to specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Support representatives.

Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

## Purchaser Notification, continued

Limited Use Label License No. 5: Invitrogen Technology

The purchase of this product conveys to the buyer the nontransferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For products that are subject to multiple limited use label licenses, the most restrictive terms apply. Invitrogen Corporation will not assert a claim against the buyer of infringement of patents owned or controlled by Invitrogen Corporation which cover this product based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Licensing Department, Invitrogen Corporation, 5791 Van Allen Way, Carlsbad, California 92008. Phone (760) 603-7200. Fax (760) 602-6500. Email: outlicensing@invitrogen.com.

## References

- Anjos-Afonso, F., Siapati, E. K., and Bonnet, D. (2004) In vivo contribution of murine mesenchymal stem cells into multiple cell-types under minimal damage conditions. J Cell Sci 117, 5655-5664
- Bruder, S. P., Jaiswal, N., and Haynesworth, S. E. (1997) Growth kinetics, self-renewal, and the osteogenic potential of purified human mesenchymal stem cells during extensive subcultivation and following cryopreservation. J Cell Biochem *64*, 278-294
- De Ugarte, D. A., Morizono, K., Elbarbary, A., Alfonso, Z., Zuk, P. A., Zhu, M., Dragoo, J. L., Ashjian, P., Thomas, B., Benhaim, P., Chen, I., Fraser, J., and Hedrick, M. H. (2003) Comparison of multi-lineage cells from human adipose tissue and bone marrow. Cells Tissues Organs 174, 101-109
- Deng, W., Obrocka, M., Fischer, I., and Prockop, D. J. (2001) In vitro differentiation of human marrow stromal cells into early progenitors of neural cells by conditions that increase intracellular cyclic AMP. Biochem Biophys Res Commun 282, 148-152
- Djouad, F., Plence, P., Bony, C., Tropel, P., Apparailly, F., Sany, J., Noel, D., and Jorgensen, C. (2003) Immunosuppressive effect of mesenchymal stem cells favors tumor growth in allogeneic animals. Blood 102, 3837-3844
- Gojo, S., Gojo, N., Takeda, Y., Mori, T., Abe, H., Kyo, S., Hata, J., and Umezawa, A. (2003) In vivo cardiovasculogenesis by direct injection of isolated adult mesenchymal stem cells. Exp Cell Res 288, 51-59
- Han, S. S., Kang, D. Y., Mujtaba, T., Rao, M. S., and Fischer, I. (2002) Grafted lineage-restricted precursors differentiate exclusively into neurons in the adult spinal cord. Exp Neurol *177*, 360-375
- Han, S. S., Liu, Y., Tyler-Polsz, C., Rao, M. S., and Fischer, I. (2004)

  Transplantation of glial-restricted precursor cells into the adult spinal cord: survival, glial-specific differentiation, and preferential migration in white matter. Glia 45, 1-16
- Houghton, J., Stoicov, C., Nomura, S., Rogers, A. B., Carlson, J., Li, H., Cai, X., Fox, J. G., Goldenring, J. R., and Wang, T. C. (2004) Gastric cancer originating from bone marrow-derived cells. Science *306*, 1568-1571
- Kinnaird, T., Stabile, E., Burnett, M. S., Shou, M., Lee, C. W., Barr, S., Fuchs, S., and Epstein, S. E. (2004) Local delivery of marrow-derived stromal cells augments collateral perfusion through paracrine mechanisms. Circulation 109, 1543-1549
- Kisseberth, W. C., Brettingen, N. T., Lohse, J. K., and Sandgren, E. P. (1999) Ubiquitous expression of marker transgenes in mice and rats. Dev Biol 214, 128-138

## References, continued

- Krampera, M., Glennie, S., Dyson, J., Scott, D., Laylor, R., Simpson, E., and Dazzi, F. (2003) Bone marrow mesenchymal stem cells inhibit the response of naive and memory antigen-specific T cells to their cognate peptide. Blood *101*, 3722-3729
- Meirelles Lda, S., and Nardi, N. B. (2003) Murine marrow-derived mesenchymal stem cell: isolation, in vitro expansion, and characterization. Br J Haematol 123, 702-711
- Moscoso, I., Centeno, A., Lopez, E., Rodriguez-Barbosa, J. I., Santamarina, I., Filgueira, P., Sanchez, M. J., Dominguez-Perles, R., Penuelas-Rivas, G., and Domenech, N. (2005) Differentiation "in vitro" of primary and immortalized porcine mesenchymal stem cells into cardiomyocytes for cell transplantation. Transplant Proc 37, 481-482
- Mujtaba, T., Han, S. S., Fischer, I., Sandgren, E. P., and Rao, M. S. (2002) Stable expression of the alkaline phosphatase marker gene by neural cells in culture and after transplantation into the CNS using cells derived from a transgenic rat. Exp Neurol *174*, 48-57
- Oh, J. Y., Kim, M. K., Shin, M. S., Lee, H. J., Ko, J. H., Wee, W. R., and Lee, J. H. (2008) The anti-inflammatory and anti-angiogenic role of mesenchymal stem cells in corneal wound healing following chemical injury. Stem Cells 26, 1047-1055
- Olivares, E. L., Ribeiro, V. P., Werneck de Castro, J. P., Ribeiro, K. C., Mattos, E. C., Goldenberg, R. C., Mill, J. G., Dohmann, H. F., dos Santos, R. R., de Carvalho, A. C., and Masuda, M. O. (2004) Bone marrow stromal cells improve cardiac performance in healed infarcted rat hearts. Am J Physiol Heart Circ Physiol 287, H464-470
- Orlic, D., Kajstura, J., Chimenti, S., Jakoniuk, I., Anderson, S. M., Li, B., Pickel, J., McKay, R., Nadal-Ginard, B., Bodine, D. M., Leri, A., and Anversa, P. (2001) Bone marrow cells regenerate infarcted myocardium. Nature 410, 701-705
- Phinney, D. G., Kopen, G., Isaacson, R. L., and Prockop, D. J. (1999) Plastic adherent stromal cells from the bone marrow of commonly used strains of inbred mice: variations in yield, growth, and differentiation. J Cell Biochem 72, 570-585
- Pittenger, M. F., Mackay, A. M., Beck, S. C., Jaiswal, R. K., Douglas, R., Mosca, J. D., Moorman, M. A., Simonetti, D. W., Craig, S., and Marshak, D. R. (1999) Multilineage potential of adult human mesenchymal stem cells. Science 284, 143-147
- Singh, S. K., Hawkins, C., Clarke, I. D., Squire, J. A., Bayani, J., Hide, T., Henkelman, R. M., Cusimano, M. D., and Dirks, P. B. (2004) Identification of human brain tumour initiating cells. Nature 432, 396-401

## References, continued

- Spees, J. L., Olson, S. D., Ylostalo, J., Lynch, P. J., Smith, J., Perry, A., Peister, A., Wang, M. Y., and Prockop, D. J. (2003) Differentiation, cell fusion, and nuclear fusion during ex vivo repair of epithelium by human adult stem cells from bone marrow stroma. Proc Natl Acad Sci U S A 100, 2397-2402
- Wakitani, S., Saito, T., and Caplan, A. I. (1995) Myogenic cells derived from rat bone marrow mesenchymal stem cells exposed to 5-azacytidine. Muscle Nerve 18, 1417-1426
- Wu, Y. Y., Mujtaba, T., Han, S. S., Fischer, I., and Rao, M. S. (2002) Isolation of a glial-restricted tripotential cell line from embryonic spinal cord cultures. Glia 38, 65-79

©2008 Invitrogen Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

## **Notes**

## **Notes**



#### **Corporate Headquarters**

Invitrogen Corporation 5791 Van Allen Way Carlsbad, CA 92008 T: 1 760 603 7200 F: 1 760 602 6500 E: tech\_support@invitrogen.com

For country-specific contact information, visit our web site at www.invitrogen.com