

CULTREX[®] **Instructions**
For Research Use Only. Not For Use In Diagnostic Procedures

Mesenchymal Stem Cell Osteogenic Differentiation Kit

Cat # 5011-024-K

**Reagents for differentiating Mesenchymal
Stem Cells into Osteoblasts**

Mesenchymal Stem Cell Osteogenic Differentiation Kit

Cat# 5011-024-K

Reagents for differentiating Mesenchymal
Stem Cells into Osteoblasts

Table of Contents

	Page
I. Introduction	1
II. Safety Information	1
III. Description of Equipment	1
IV. Materials Required But Not Supplied	2
V. Reagent Preparation	2
VI. Protocol	3
VII. References	10
VIII. Troubleshooting	11
IX. Related Products Available From Trevigen	13

©2009, Trevigen Inc. All rights reserved. Trevigen, Cultrex and PathClear are registered trademarks of Trevigen, Inc. Microsoft and Excel are trademarks of Microsoft Corporation. Corning is a registered trademark of Corning, Inc.

Introduction

Mesenchymal Stem Cells can differentiate into osteogenic cells^{1,2} that will deposit and mineralize extracellular matrix proteins,^{2,3} which is required for the formation of new bone. Measuring the amount of mineralized matrix is an established method for evaluating osteogenic differentiation.²

Mesenchymal stem cells are a self-renewing population of multipotent cells, present in the bone marrow and many other adult tissues, which can differentiate into multiple lineage-specific cells that form bone, fat, cartilage, muscle and tendon.¹⁻³ The differentiation of mesenchymal stem cells into osteogenic cells can be a useful tool for understanding mechanisms involved in bone formation and for identifying potential treatments for osteoporosis, bone repair and genetic disorders such as osteogenesis imperfecta.³

Trevigen's **Mesenchymal Stem Cell Osteogenic Differentiation Kit** contains reagents optimized to direct rat mesenchymal stem cells (RMSC) (cat# 5000-001-01) or human mesenchymal stem cells (HMSC) grown on Cultrex[®] Rat Collagen I⁹ (cat# 3440-100-01) to undergo osteogenic differentiation in a defined growth medium supplemented with ascorbic acid, β-glycerol phosphate, dexamethasone and within 14 days. Osteogenic differentiation is detected by staining with Alizarin Red S⁸. Alizarin red will stain the calcium deposits that occur when differentiated mesenchymal stem cells mineralize the extracellular matrix.⁸

II. Precautions and Limitations

1. For Research Use Only. Not for use in diagnostic procedures.
2. The physical, chemical, and toxicological properties of these products may not yet have been fully investigated; therefore, Trevigen recommends the use of gloves, lab coats, and eye protection while using these chemical reagents. Trevigen assumes no liability for damage resulting from handling or contact with these products.

III. Materials Supplied

<u>Component</u> ¹	<u>Quantity</u>	<u>Storage</u>	<u>Catalog#</u>
Dexamethasone	12 µl	-20°C*	5010-024-01
Ascorbic Acid	1.2 ml	4°C	5011-024-01
β-glycerol Phosphate	1.2 ml	4°C	5011-024-02
Cultrex [®] Rat Collagen I ²	500 µl	4°C	3440-001-01
Alizarin Red S ³	25 ml	RT in dark	5011-024-03
Stain Solubilization Solution ³	35 ml	RT	5011-024-04

1: Reagents are sufficient to differentiate one 24 well plate for up to 21 days

2: Sufficient rat tail collagen type I is provided to coat two 24 well plates

3: Enough reagent is provided to test two 24 well plates.

*Separate from shipping box and store at specified temperature

IV. Materials/Equipment Required But Not Supplied

Equipment

1. 1 - 20 μ l, 20 - 200 μ l, and 200 - 1000 μ l pipettors
2. Laminar flow hood or clean room
3. 37°C CO₂ incubator
4. 37°C Water Bath
5. Hemocytometer or other means to count cells
6. Inverted standard or phase microscope
7. pipette aid
8. Liquid Nitrogen Storage
9. Low speed swinging bucket centrifuge and tubes for cell harvesting
10. Plate Reader

Reagents

1. Mesenchymal Stem Cells (Trevigen's RMSC 5000-001-01 or human equivalent)
2. Qualified MSC Medium (for RMSC: cat# 5000-500-03) or equivalent
3. Cell Harvesting Reagent, trypsin, dispase, etc.
4. Qualified MSC FBS (for RMSC: cat# 5000-050-02)
5. Antibiotic Supplement for Medium (optional)
6. Sterile PBS (Mg²⁺, Ca²⁺ free) or HBSS
7. Trypan blue or equivalent viability stain
8. DMSO
9. 70% Ethanol
10. 10% Formalin

Disposables

1. Cell culture flask, 25 cm², 75 cm², or 185 cm²
2. 24 well Tissue Culture Plates
3. 15 ml tubes
4. 0.22 μ m Filter Unit (optional)
5. 60 ml Syringe
6. 1 - 200 μ l and 200 - 1000 μ l pipette tips
7. 1, 5 and 10 ml serological pipettes
8. gloves
9. Clear 96-well plate
10. Cryovials

V. Reagent Preparation

These procedures should be performed in a biological hood utilizing aseptic technique to prevent contamination.

1. Mesenchymal Complete Growth Medium

For 250 ml of Medium:

Qualified MSC Medium:	225 ml
Qualified FBS:	25 ml

Optional: media can be filter sterilized before use

Store media @ 4°C for one month

Ensure media are at room temperature or 37°C prior to use

2. Mesenchymal Freeze Medium

For 10 ml of Medium:

Qualified MSC Medium:	4 ml
Qualified FBS:	4 ml
DMSO:	2 ml

Mix 1:1 with Growth Medium before use

3. Osteogenic Differentiation Medium*

For 13 ml of Medium

Mesenchymal Growth Medium (From V.I):	12.96 ml
Dexamethasone	1.3 µl
Ascorbic Acid	130 µl
β-Glycerol Phosphate	130 µl

Invert several times to mix.

*Make up Fresh Prior to use

Ensure medium is at 37°C prior to use, addition of cold media can induce "peeling" of cells.

4. Collagen Solution for Coating

To coat one 24 well plate, add 250 µl Cultrex® Collagen I to 12.5 ml of Qualified RMSC Medium (adjust volume as needed accordingly to number of wells)

VI: Protocol

These procedures should be performed in a biological hood utilizing aseptic technique to prevent contamination.

A. Thawing Mesenchymal Stem Cells:

1. Prepare Mesenchymal Complete Growth Medium (Section V.1):
2. Prewarm Complete Growth Medium to 37 °C by placing in 37°C H₂O bath or in Tissue Culture Incubator
3. Immediately before use, remove vial of cryopreserved mesenchymal stem cells (MSC) (e.g. Rat MSC cat# 5000-001-01) from liquid nitrogen freezer.
4. Thaw frozen RMSC quickly in a 37°C H₂O bath.
 - a. Ensure cells are completely thawed before proceeding
 - b. Do not leave cells at 37°C for fast thawing
5. Spray down bottle containing Complete Growth Medium and ampoule containing cells with 70% EtOH before placing in Tissue Culture Hood
6. Aseptically, transfer the thawed cells to a 15 ml conical tube with a 5 ml pipette

7. Add 1 ml of warm Complete Growth Medium to ampoule using a 5 ml pipette, dropwise with gentle swirling of the cells and transfer to a 15 ml conical tube dropwise, gently swirling to mix between drops
8. Add 1 ml of warm Complete Growth Medium to 15 ml conical tube containing cells, gently swirling to mix between drops
Note: Total Volume should be about 3 ml
9. Centrifuge 15 ml conical tube at 250 x g for 5 minutes at room temperature
10. Remove supernatant gently to avoid disturbing cell pellet
11. Resuspend cell pellet in 1 ml of fresh Complete Growth Medium
12. Count cells on hemocytometer (per standard protocol)
13. For Trevigen's RMSC (cat# 5000-001-01): Plate cells at a density of 5.4×10^3 cell/cm² in a T-75 tissue culture treated flask (4.05×10^5 cells) in a total volume of 12-15 ml Complete Growth Medium
Notes: Recommend Corning Tissue Culture Treated Plastic. One vial is sufficient to seed two T-75 or one T-185 flask
14. For other MSC cells, optimal plating density will have to empirically determined and will be cell line specific.
15. Place Tissue Culture Flask/Dish in 5% CO₂ Tissue Culture Incubator at 37°C
16. Change medium on flasks on the next day.

B. Growing Mesenchymal Stem Cells:

1. Medium Change (Culture Medium should be changed every 3-4 days)
 - a. Warm Complete Growth Medium to 37 °C by placing in 37°C H₂O bath or in Tissue Culture Incubator
 - b. Spray down bottle containing growth medium with 70% EOH before placing in Tissue Culture Hood
 - c. Remove medium from T-75 flask containing Mesenchymal Stem Cells
 - d. Add 12-15 ml of fresh Complete Growth Medium
 - e. Discard used medium appropriately
2. Passaging Mesenchymal Stem Cells
When MSCs become 70- 80% confluent they are ready to be split. If allowed to over-grow, these cells will lay down a matrix and start to differentiate. As a result, the cells will peel off of the plastic which markedly reduces the ability to passage them.
 - a. Warm Mesenchymal Complete Growth Medium and trypsin solution to 37 °C by placing in 37 °C H₂O bath or in Tissue Culture Incubator
 - b. Spray down bottle containing growth medium, and trypsin bottle with 70% EtOH before placing in Tissue Culture Hood
 - c. Remove medium from T-75 flask containing Mesenchymal Stem Cells
 - d. Gently wash flask with 5-10 ml of sterile 1X PBS (Ca²⁺ and Mg²⁺ free)
 - e. Remove PBS
 - f. Add 3 ml of Trypsin solution to each flask place at 37°C in Tissue Culture Incubator for 3-5 minutes (until cells are no longer attached to plate, this should take no longer than 5 minutes).
 - g. Add 5 ml of warm Complete Growth Medium to flask
 - h. Transfer cells to 15 ml conical tube
 - i. Centrifuge 15 ml conical tube at 200 x g for 3 minutes at room temperature

- j. Remove supernatant gently to avoid disturbing cell pellet
- k. Resuspend cell pellet in 2 ml of fresh Complete Growth Medium
- l. Count cells on hemocytometer (per standard protocol)
- m. For Trevigen's Rat MSC: Passage cells at a density of 5.4×10^3 cell/cm². For a T-75 flask add 4.05×10^5 cells in 12-15 ml of complete growth medium
- n. For other MSC cells, optimal plating density will have to empirically determined and will be cell line specific

Note: One flask of 70-80% confluent cells should be enough to be plated into 1-3 T-75 flasks.

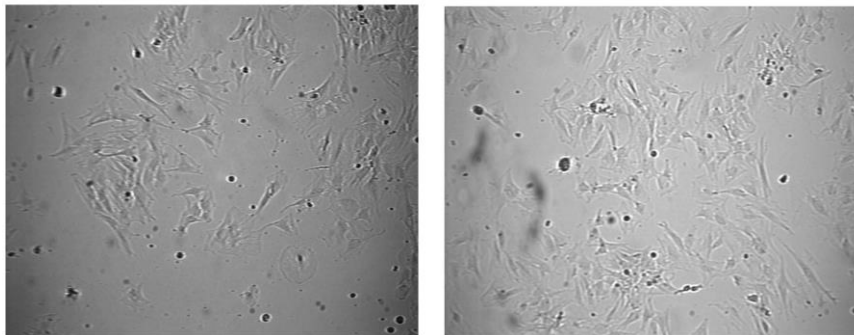


Figure 1: 10X Bright Field Images of Rat Mesenchymal Stem Cells.

C. Freezing Cells

1. Warm Complete Growth Medium and trypsin solution to 37 °C by placing in 37 °C H₂O bath or in Tissue Culture Incubator
2. Make Freeze Medium (see Section V.2), adjust volume according to volume needed (*Will be mixed 1:1 with Complete Growth Medium*)
3. Spray down bottles containing Complete Growth Medium, trypsin, and the Freeze Medium tube with 70% EtOH before placing in Tissue Culture Hood
4. Remove medium from T-75 flask containing Mesenchymal Stem Cells
5. Gently wash flask with 5-10 ml of sterile 1X PBS (Ca²⁺ and Mg²⁺ free)
6. Remove PBS
7. Add 3 ml of Trypsin solution to each flask place at 37 °C in Tissue Culture Incubator for 2-3 minutes (until cells are no longer attached to the plate, which should take no more than 5 minutes).
8. Add 5 ml of Complete Growth Medium to flask
9. Transfer cells to 15 ml conical tube
10. Centrifuge 15 ml conical tube at 200 x g for 3 minutes at room temperature
11. Remove supernatant gently to avoid disturbing cell pellet
12. Resuspend cell pellet in 2 ml of Complete Growth Medium
13. Count cells on hemocytometer (per standard protocol)
14. Dilute cells to a desired concentration for freezing.

Notes: We recommend a concentration of no less than 1×10^6 cell/ml. This concentration is enough to seed one T-75 flask (remember, the cells will be

diluted 1:1 with Freeze Medium). One T-75 flask will provide enough cells for 1-3 vials of 5×10^6 cells/vial.

15. Add an equal volume of Freeze Medium to the cells, mix gently
16. Aliquot 1 ml of cells into labeled cryovials
17. Place on ice of 15-30 minutes
18. Transfer cells to -80°C freezer and incubate overnight
19. Transfer to liquid nitrogen freezer for long term storage

Note: Vapor phase is recommended to ensure viability

D. Osteogenic Differentiation of Mesenchymal Stem Cells

When MSCs become 70- 80% confluent they are ready to be split. If allowed to over-grow, these cells will lay down a matrix and start to differentiate, thereby allowing them to peel off which markedly reduces the ability to passage them.

Day before plating of the cells, coat 24-well plate with Cultrex[®] Collagen I (Rat)

a: Add 500 μl of Collagen Solution for Coating (sect V.4) to each well

b: Place plate at 37°C CO_2 incubator overnight

Note: If not working in a 24 well plate, adjust coating volume accordingly.

1. Plating Cells

- a. Warm Complete Growth Medium and trypsin solution to 37°C by placing in 37°C H_2O bath or in Tissue Culture Incubator
- b. Spray down bottle containing Complete Growth Medium, and trypsin bottle with 70% EtOH before placing in hood
- c. Remove medium from T-75 flask containing Mesenchymal Stem Cells
- d. Gently wash flask with 5-10 ml of sterile 1X PBS (Ca^{2+} and Mg^{2+} free)
- e. Remove PBS
- f. Add 3 ml of Trypsin to each flask and place at 37°C in Tissue Culture Incubator for 3-5 minutes (until cells are no longer attached to plate, which should take no more than 5 minutes).
- g. Add 5 ml of warm Complete Growth Medium to flask
- h. Transfer cells to 15 ml conical tube
- i. Centrifuge 15 ml conical tube at 200 x g for 3 minutes at room temperature
- j. Remove supernatant gently to avoid disturbing cell pellet
- k. Resuspend cell pellet in 2 ml of fresh Complete Growth Medium
- l. Count cells on hemocytometer (per standard protocol)
- m. Remove medium from Collagen coated 24-well plate
- n. For Trevigen's Rat MSC: Plate cells at 5,000 cells/ cm^2 . For a 24-well plate add 10,000 cells/well in 0.5 ml of complete growth medium per well.
- o. For HMSC [e.g. Zen-Bio, Inc cat# ACS-F], Plate cells at 30,000 cells/ cm^2 . For a 24-well plate add 60,000 cells/well in 0.5 ml of complete growth medium per well.
- p. For other MSC, initial plating concentration will have to be optimized for endogenous matrix deposition by bone marrow derived MSC.
Note: If not using 24-well plate, adjust medium volumes accordingly. Components are sufficient to differentiate one 24-well plate using the stated medium volumes.
- q. Incubate overnight at 37°C

2. Differentiation

- a. Observe the plates under phase microscope to ensure that the mesenchymal stem cells have attached to the plate.
- b. Medium needs to be changed every 3-4 days for 4 days.
- c. Procedure:

- i. Day 1

1. Label wells for Differentiated or Undifferentiated control groups
2. For Differentiation: Remove Complete Growth Medium and add 0.5 ml of freshly prepared Osteogenic Differentiation Medium (sect. V.3) to each well
3. For Undifferentiated controls: Remove Complete Growth Medium and add 0.5 ml of fresh Complete Growth Medium (V.1) to control wells

- ii. Days 4, 7, 10, 13 (if extending culture to 21 days continue changing medium every 3-4 days)

1. For Differentiation: Remove spent Differentiation Medium and add 0.5 ml of freshly prepared Osteogenic Differentiation Medium (section V.3) to each well
2. For Undifferentiated controls: Remove growth medium and add 0.5 ml of Complete Growth Medium (V.1) to control wells

Note: Confluent cultures undergoing osteogenic differentiation tend to lift off or "peel." Osteogenic differentiation occurs within 7 days of growth in osteogenic medium. Peeling can begin to occur within 12 days in culture. Osteogenic differentiation induction can be extended up to 21 days depending on cell number and reduced growth rate.

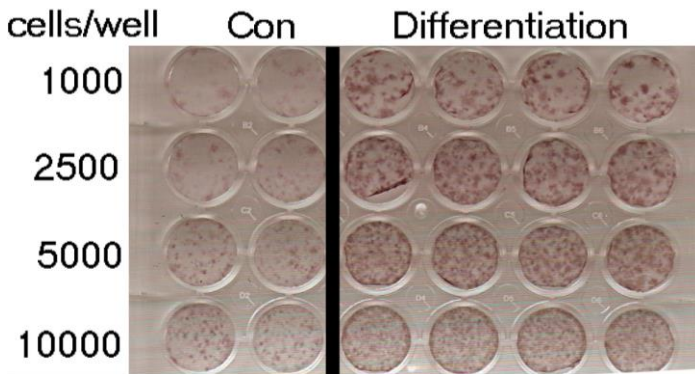


Figure 2: Rat MSCs [Trevigen cat# 5000-001-01] grown on Rat Tail Collagen I-Coated wells for 14 days in Osteogenic Differentiation Medium, fixed and stained with Alizarin Red S

- d. Evaluation of Osteogenic Differentiation: Alizarin Red S staining (Day 4):
 - i. Remove the media from each well
 - ii. Gently wash wells with 1 ml of 1X PBS
 - iii. Remove PBS
 - iv. Add 0.5 ml of 10% Formalin in PBS to each well
 - v. Incubate plate at room temperature for 30 minutes
 - vi. Remove 10% Formalin
 - vii. Wash each well with 1 ml of ddH₂O
 - viii. Add 0.5 ml of ddH₂O to each well and incubate plate for 5 minutes
 - ix. Remove ddH₂O
 - x. Add 0.5 ml of Alizarin Red S to each well

- xi. Incubate plate for 20 minutes at room temperature with gentle shaking (shaking optional)
- xii. Remove Alizarin Red S solution
- xiii. Wash wells twice with 1 ml of ddH₂O
- xiv. Remove ddH₂O and let the wells air dry.

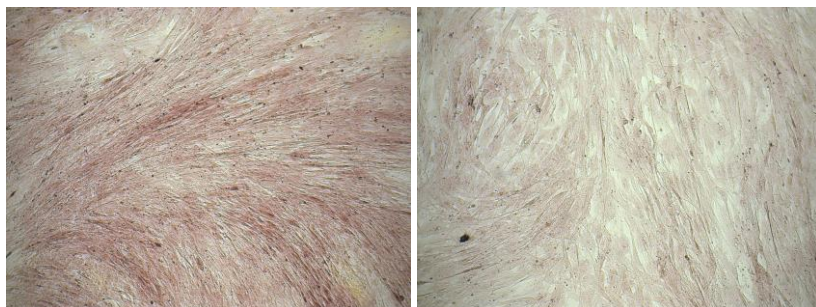


Figure 3: HMSC [Zen-Bio, Inc, cat# ACS-F] grown on Rat Tail Collagen I-Coated wells for 14 days in Osteogenic Differentiation Medium (left), grown in control medium (right), fixed and stained with Alizarin Red S

e. Quantification of Alizarin Red S Stain

i. Standard Curve

- 1. Prepare Samples according to the following table. Use 1.7 ml tubes

	Standard (concentration of Alizarin Red S stain)	Alizarin Red Stain	Solubilization Solution
A	1 mg/ml	50 µl Stock Solution	950 µl
B	500 µg/ml	500 µl of A	500 µl
C	250 µg/ml	500 µl of B	500 µl
D	100 µg/ml	400 µl of C	600 µl
E	50 µg/ml	500 µl of D	500 µl
F	25 µg/ml	500 µl of E	500 µl
G	10 µg/ml	400 µl of F	600 µl
H	5 µg/ml	500 µl of G	500 µl
I	0 µg/ml	0 µl	500 µl

- 2. Transfer 150 µl of standard solutions B-I to a clear 96-well plate in triplicate.

ii. Extraction of Alizarin Red Stain from wells

- 1. Add 500 µl of Stain Solubilization Solution to each well of the 24 well plate
- 2. Incubate plate at room temperature for 10 min.
- 3. Transfer 150 µl of sample to the 96-well clear plate in triplicate

iii. Read absorbance of standards and extracted Alizarin Red Stain on a plate reader at 405 nm

iv. Calculate the Alizarin Red S present in each sample relative to the standard curve:

1. Plot the standards with Absorbance at A_{405} on the Y-axis and concentration of Alizarin Red S standard on the X-axis, if using Excel, use the X-Y graph feature. Be sure to subtract the $0 \mu\text{g/ml}$ (I) value from all standard (and sample) values.
2. Insert a trendline and determine the equation of the standard curve line. Your trendline should have an R^2 value ≥ 0.975
3. Take sample readings at A_{405} (use $150 \mu\text{l}$ per triplicate sample). Using the equation from the standard curve, calculate the amount of Alizarin Red S present in the samples.
4. Plot the amount of Alizarin Red S present in the samples and controls on a bar graph.
5. The amount of Alizarin Red S stain present in a sample is relative to the amount of differentiation occurring in the culture. By day 14 of the assay, cells will be confluent so the amount of stain is representative of the differentiation state; not the number of cells present.

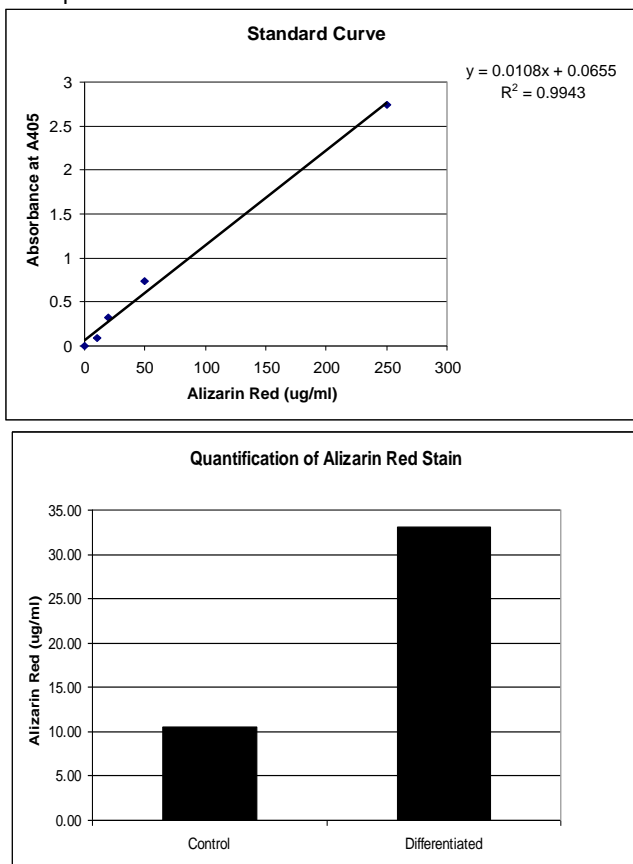


Figure 5: Quantitation of Alizarin Red Staining of Rat Mesenchymal Stem Cells [Trevigen, Inc cat# 5000-001-01] cultured in Osteogenic Differentiation Medium for 14 days

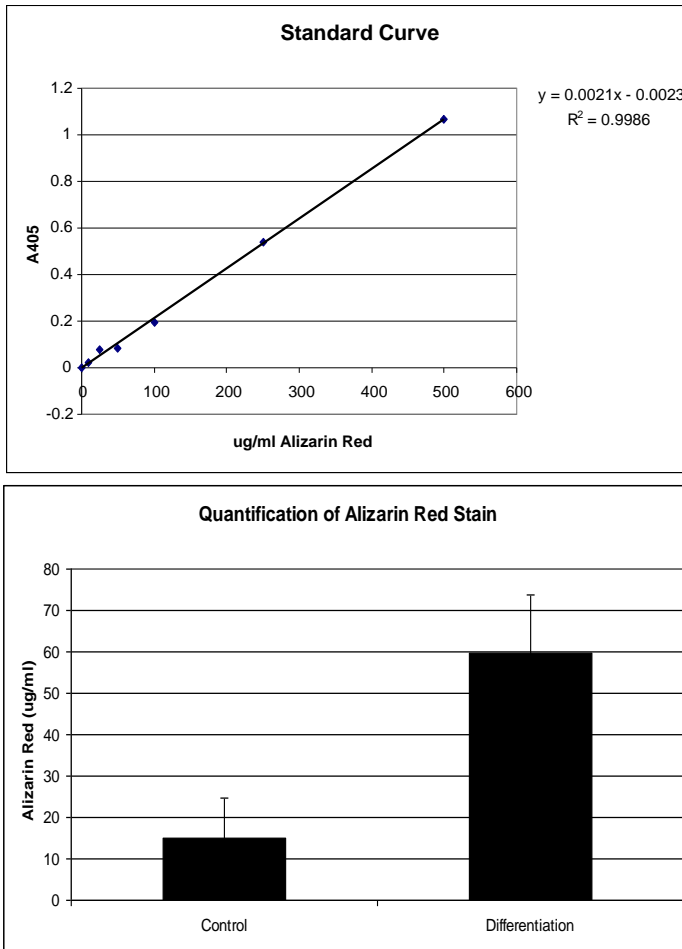


Figure 6: Quantitation of Alizarin Red Staining of Human Mesenchymal Stem Cells [Zen-Bio, Inc cat# ACS-F] cultured in Osteogenic Differentiation Medium for 14 days

VII. References

1. Satija NK, Gurudutta GU, Sharma S, Afrin F, Gupta P, Verma K, Singh VK, Tripathi RP. 2007 *Mesenchymal Stem Cells: Molecular Targets for Tissue Engineering*. *Stem Cells and Development* 16:7-23.
2. Jones JL, Scotting P, Sottile V. 2007. *Adult Mesenchymal stem cell: Differentiation potential and therapeutic applications*. *J Postgrad Med* 53(2): 121-127.
3. Li F, Wang, X, Niyibizi C. 2007. *Distribution of Single-Cell Expanded Marrow Derived Progenitors in a Developing Mouse Model of Osteogenesis Imperfecta Following Systemic Treatment*. *Stem Cells*. 25:3183-3193.
4. Phinney DG, Prockop DJ. 2007. Concise Review: Mesenchymal Stem/Multipotent Stromal Cells: The State of Transdifferentiation and Models of Tissue Repair-Current Views. *Stem Cells* 25: 2896-2902.

5. Kolf CM, Cho E, Tuan RS. 2007. Biology of Adult Mesenchymal Stem Cells: Regulation of Niche, Self-Renewal and Differentiation. *Arthritis Research and Therapy* 9:204
6. Javazon EH, Colter DC, Schwarz EJ, Prockop DJ. 2001. Rat Marrow Stromal Cells are More Sensitive to Plating Density and Expand More Rapidly from Single-Cell-Derived Colonies than Human Marrow Stromal Cells. *Stem Cells* 19:219-225.
7. Li Yi, McIntosh K, Chen J, Zhang C, Gao, Q, Borneman J, Ragniski K, Mitchell J, Shen L, Zhang J, Lu D, Chopp M. Allogeneic bone marrow stromal cells promote glial-axonal remodeling without immunologic sensitization after stroke in rats. *Experimental Neurology* 198:313-325.
8. Reger RL, Tucker AH, Wolfe MR. 2008. Differentiation and Characterization of Human MSCs. In Prockop DJ, Phinney DG, Bunnell BA (eds) *Methods in Molecular Biology Mesenchymal Stem Cells: Methods and Protocols* vol 449. Humana Press, Totawa, NJ, USA pp 93-107.
9. Salasznyk RM, Williams WA, Boskey A, Batorsky A, Plopper GE. 2004. *Adhesion to Vitronectin and Collagen I Promotes Osteogenic Differentiation of Human Mesenchymal Stem Cells*. J. Biomedicine and Biotechnology 2004: 24-34.

VIII. Troubleshooting

PROBLEM	CAUSE	ACTION
Poor Viability from initial freeze	To rough in thawing of cells	Ensure medium is added slowly to re-equilibrate the MSC from freeze medium Ensure cells were removed from freeze medium immediately after vial has been thawed Ensure Vial of cells was thawed at 37 °C Fresh media was pre-warmed to 37 °C
Poor proliferation	Medium and Fetal Bovine Serum not optimized for support of MSC growth	For RMSC: Use Qualified RMSC Medium and FBS from Trevigen For other MSC: Try different lots of medium and fetal bovine serum or purchase qualified medium and serum appropriate for cell origin.
	Tissue Culture Labware not ideal for MSC	Use Corning or Nunc Treated Labware
	Frequency of Medium Change	Ensure medium is changed every 3-4 days Ensure pH of fresh medium has not changed.
	CO ₂ Incubator not humidified	Add sterile water to CO ₂ incubator per manufacturer's instructions
	No gas exchange is allowed by flask	Ensure cap is loosened to allow air gas or use vented flask
Cells were allowed to become over confluent and lay down matrix		Extend time in trypsin
		Titrate cells to remove as many cells as possible from matrix Remove visual matrix aggregated from tube before spinning (will reduce cell recovery) Pass cell suspension through cell strainer (will reduce cell recovery)

PROBLEM	CAUSE	ACTION
Cells were clumpy after passaging, limited recovery of single cells.	Cells were allowed to become over confluent and lay down matrix	Extend time in trypsin Titrate cells to remove as many cells as possible from matrix Remove visual matrix aggregated from tube before spinning (will reduce cell recovery) Pass cell suspension through cell strainer (will reduce cell recovery)
Contamination of Cells	Contaminated Medium Improper Aseptic Technique Hood is working improperly Contaminated CO ₂ Incubator.	To prevent contamination, filter medium through a 0.22 µm filter before use <i>Never use contaminated medium once cloudy or after microorganisms are visible under the microscope.</i> Spray down hands, reagents and hood with 70% ethanol before opening any flasks Check to make sure blower is on and functioning Ensure hood is currently certified Wipe down hood with 70% ethanol Ensure CO ₂ incubator is free of microbial growth
Poor Cell Recovery from flask (for cell growth)	Too high seeding density	Passage cells at lower confluency
No Osteogenic Differentiation or low levels of differentiation	Serum Medium was not prepared properly Change basal medium type Too long between medium changes Cell were plated too sparse	Cells were grown in serum that did not support undifferentiated growth, for RMSC use Trevigen's Qualified RMSC FBS. For other MSC try different fetal bovine serum Ensure the supplements were added to medium in proper amounts Ensure basal medium contains calcium Ensure medium is being changed every 3-4 days Plate cells at a higher density
Osteogenic differentiation present in control medium	Cells plated too high density	Plate cells at lower density
Cells peeling from wells	Cells over-confluent	Plate cells at a lower concentration
High background in control wells for during quantification	Presence of non-specific stain in the well	Increase number of washes

IX. Related Products Available From Trevigen

Contact Trevigen for details of our unique product line for studying DNA damage and repair. All of Trevigen's kits include highly qualified enzymes, substrates.

Differentiation:

Catalog #	Description	Size
5000-001-K	Cultrex® Rat Mesenchymal Stem Cell Starter Kit	1 kit
5000-001-R	Cultrex® Rat Mesenchymal Stem Cell Replenisher Kit	1 kit
5000-001-01	Cultrex® Rat Mesenchymal Stem Cells	1 vial
5010-024-K	Cultrex® Adipogenic Differentiation Kit	24 samples

3D Culture Kits:

Catalog #	Description	Size
3445-096-K	Cultrex® 3D Culture BME Cell Proliferation Assay Kit	96 tests
3446-096-K	Cultrex® 3D Culture Laminin I Cell Proliferation Assay	96 tests
3447-096-K	Cultrex® 3D Culture 96 Well Collagen I Cell Prolif Assay	96 tests
3448-020-K	Cultrex® 3D Culture Cell Harvesting Kit	96 tests

Invasion/Migration Kits:

Catalog#	Description	Size
3455-024-K	Cultrex® 24 Well BME Cell Invasion Assay	24 inserts
3460-024-K	CultreCoat® 24 Well BME-Coated Cell Invasion Assay	24 inserts
3465-096-K	Cultrex® 96 Well Cell Migration Assay	96 samples
3465-024-K	Cultrex® 24 Well Cell Migration Assay	12 samples
3455-096-K	Cultrex® 96 well BME Cell Invasion Assay	96 samples
3456-096-K	Cultrex® 96 well Laminin I Cell Invasion Assay	96 samples
3457-096-K	Cultrex® Collagen I Cell Invasion Assay	96 samples
3458-096-K	Cultrex® Collagen IV Cell Invasion Assay	96 samples

Accessories:

Catalog#	Description	Size
3415-001-02	Cultrex® Human BME, PathClear®	1 ml
3432-005-02	Cultrex® BME, PathClear®	5 ml
3432-005-01	Cultrex® BME without Phenol Red	5 ml
3431-005-01	Cultrex® BME with Phenol Red; Reduced Growth Factor	5 ml
3433-005-01	Cultrex® BME; no Phenol Red; Reduced Growth Factor	5 ml
3430-005-02	Cultrex® BME with Phenol Red, PathClear®	5 ml
3431-005-02	Cultrex® BME with Phenol Red, Reduced Growth Factor PathClear®	5 ml
3400-010-01	Cultrex® Mouse Laminin I	1 mg
3440-100-01	Cultrex® Rat Collagen I	100 mg
3410-010-01	Cultrex® Mouse Collagen IV	1 mg
3420-001-01	Cultrex® Human Fibronectin, PathClear®	1 mg
3416-001-01	Cultrex® Bovine Fibronectin, NZHD*	1 mg
3421-001-01	Cultrex® Human Vitronectin, PathClear®	50 µg
3417-001-01	Cultrex® Bovine Vitronectin, NZHD	50 µg
3438-100-01	Cultrex® Poly-L-Lysine	100 ml
3439-100-01	Cultrex® Poly-D-Lysine	100 ml
3445-048-01	Cultrex® 3-D Culture Matrix™ BME	15 ml
3446-005-01	Cultrex® 3-D Culture Matrix™ Laminin I	5 ml
3447-020-01	Cultrex® 3-D Culture Matrix™ Collagen I	100 mg
3430-005-01	Cultrex® BME with Phenol Red	5 ml
3437-100-K	Cultrex® Cell Staining Kit	100 ml
3450-048-05	CellSpere™	15 ml

*New Zealand Herd Derived

The product accompanying this document is intended for research use only and is not intended for diagnostic purposes or for use in humans.

Trevigen, Inc.

8405 Helgerman Ct. Gaithersburg, MD 20877

Tel: 1-800-873-8443 • 301-216-2800

Fax: 301-560-4973

e-mail: info@trevigen.com

www.trevigen.com



Please
Recycle