

CULTREX[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

3-D Culture Matrix[™] Rat Collagen I

Catalog #: 3447-020-01

Size: 20 ml

Description: 3-D Culture is an innovative approach to modeling the morphological effects of early oncogenesis on three-dimensional microenvironments. When healthy, differentiating cells exhibit a structured, polarized morphology that is critical for cellular formation and function. During carcinoma development, cell cycle controls associated with cellular development, proliferation and death are lost, and as a result, these structures are disrupted. In effect, the morphology of these structures can be used as a measure to study factors in early carcinoma development. In an attempt at standardization, J. Debnath, *et al.* published guidelines for execution of this assay using MCF-10A mammary epithelial cells as a model.¹ To aid in the advancement of this technology, Trevigen has developed the Cultrex[®] 3-D Culture Matrix[™] product line to provide reagents specifically produced for and qualified in 3-D culture studies. The 3-D Culture Matrix[™] Collagen I may be used as a gel on which to grow cells or a media additive alone or in concert with other basement membrane components to study cellular growth and differentiation in three dimensions *in vitro*.

Type I Collagen is the major structural component of extracellular matrices found in connective tissue and internal organs, but is most prevalent in the dermis, tendons, and bone. It is a 300 kDa molecule composed of two alpha₁(I) chains and one alpha₂(I) chain that spontaneously forms a triple helix scaffold at a neutral pH and 37°C. This phenomenon can be exploited to promote cell attachment, growth, differentiation, migration, and tissue morphogenesis during development.

To provide the most standardized Collagen I for use in 3-D cultures, a special process is employed to provide material at a standard concentration of approximately 5 mg/mL. This material is then incorporated in a 3-D culture to validate efficacy.

Specifications:

Concentration: Type I Collagen provided at 5 mg/ml (Sircol Assay).

Source: Rat tail tendons

Storage Buffer: 20 mM Acetic Acid

Storage/Stability: Product is stable for a minimum of 3 months if stored at 4°C. **Do Not Freeze.**

Materials Qualification:

Gelling:

- Type I collagen forms a firm gel at neutral pH and 37°C when diluted to 0.4 mg/ml.

Functional Assays:

- Cell Attachment: Tested for the ability to promote cell attachment and spreading of HT-1080 human fibrosarcoma cells.
- 3-D Culture: Collagen I promotes attachment and growth of murine endothelial SVEC4-10 cells.

© 2009 Trevigen, Inc. All Rights Reserved. Trevigen, and Cultrex are registered trademarks, and DIVAA, 3-D Culture Matrix, and CellSpere are trademarks of Trevigen, Inc. E11/20/09v1

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

Sterility Testing:

- No bacterial or fungal growth detected after incubation at 37°C for 14 days following USP XXIV Chapter 71 sterility test.
- No mycoplasma contamination detected by PCR.
- Endotoxin concentrations \leq 20 EU/ml by LAL assay.

Gelling Procedures:

Note: To prevent contamination maintain aseptic techniques in a laminar flow biological hood throughout this procedure. Working with solutions that are pre-chilled at 4°C, and keeping solutions on ice extends the time that collagen I will remain in solution after neutralization.

Material is qualified at 1 mg/mL, and this is the recommended working concentration.

1. Place the following on ice:
 - a. Type I Collagen (5 mg/ml)
 - b. Sterile 10X PBS
 - c. Sterile, distilled water (dH₂O)
 - d. Sterile 1N NaOH (fresh)
2. Determine the concentration and final volume of Collagen needed for experimentation.
3. Determine the amount of reagents needed so that Collagen I is at the desired concentration in 1X phosphate buffered saline (PBS) neutralized by 1N NaOH:
 - a. Volume of Collagen needed = $\frac{(\text{Final conc. of Collagen}) \times (\text{Total Volume})}{(\text{Initial conc. of Collagen})}$
 - b. Volume of 10X PBS needed = $\frac{\text{Total Volume}}{10}$
 - c. Volume of 1N NaOH needed = (volume of Collagen I) x 0.023 ml
 - d. Volume of dH₂O needed = Total Volume - (sum of volumes from steps A+B+C)
4. In a sterile tube mix the 10X PBS, 1N NaOH and dH₂O.
5. Add the Collagen I to the tube and pipette up and down to mix (do not vortex).
6. Place the Collagen solution into the desired plates or dishes. Solution is stable for up to one hour on ice. Plates may be centrifuged 300 x g for 10 minutes at 4°C to prevent bubbles from forming in the gel.
7. Incubate the plate at 37°C for 1 hour to promote gel formation.

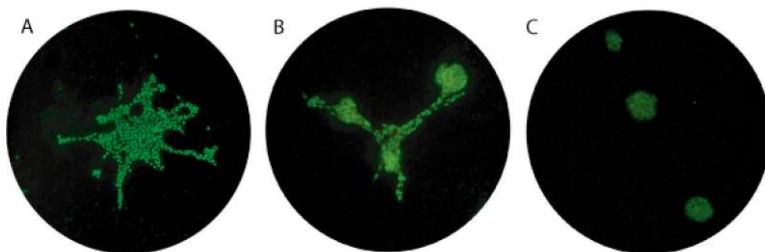
For your cell type, a gelling procedure using 7.5% (w/v) Sodium Bicarbonate for neutralization may be preferred:

1. Place the following on ice:
 - a. Type I Collagen (5 mg/ml)
 - b. Sterile 10X PBS
 - c. Sterile, distilled water (dH₂O)
 - d. 7.5% Sodium Bicarbonate, sterile
2. Determine the concentration and final volume of Collagen needed for experimentation.
3. Determine the amount of reagents needed so that Collagen I is at the desired concentration in 1X phosphate buffered saline (PBS), neutralized by 7.5% sodium bicarbonate.
 - a. Volume of Collagen needed = $\frac{(\text{Final conc. of Collagen}) \times (\text{Total Volume})}{(\text{Initial conc. of Collagen I})}$
 - b. Volume of 10X PBS needed = $\frac{\text{Total Volume}}{10}$
 - c. Volume of 7.5% sodium bicarbonate needed = (Volume of Collagen I, step a) x 0.0125 ml
 - d. Volume of dH₂O needed = Total Volume - (sum of volumes from steps A+B+C)

4. In a sterile tube mix the 10X PBS, and dH₂O and 7.5% sodium bicarbonate.
5. Add the Collagen I to the tube and pipette up and down to mix (do not vortex).
6. Place the neutralized Collagen I solution into the desired plates or dishes. This solution is stable for up to 1 hour on ice. Plates may be centrifuged 300 x g for 10 minutes at 4°C to prevent bubbles from forming in the gel.
7. Incubate the plate at 37 °C for 1 hour to promote gel formation.

High Concentration Collagen gel method:

1. Place Collagen I (5 mg/ml), 7.5% sodium bicarbonate solution, sterile tube and cell culture plate on ice.
2. Add necessary amount of Collagen I into sterile tube.
3. Add 5 µl of 7.5% sodium bicarbonate per 0.1 ml of Collagen I (5 mg/ml)
4. Pipette Collagen I up and down to mix. (Do not vortex.)
5. Place neutralized collagen into a cell culture plate. Plate may be centrifuged for 300 x g for 10 minutes at 4 °C to prevent bubbles from forming in the gel.
6. Incubate the plate at 37 °C for 1 hour to promote gel formation.



Mammary epithelial cells, MCF-10A cultured on 3-D Culture Matrix™ Collagen I are induced to differentiate with the addition of 3-D Culture Matrix™ Laminin-1 at: a) 0 mg/mL, b) 1 mg/mL, and c) 2 mg/mL.

References:

1. Debnath J, Muthuswamy SK, Brugge JS. Morphogenesis and oncogenesis of MCF-10A mammary epithelial acini grown in three-dimensional basement membrane cultures. *Methods*, 2003 pp. 256-268.
2. Chen, S., R. Revoltella, S. Papini, M. Michelini, W. Fitzgerald, J. Zimmerberg, and L. Margolis. 2003. Multilineage differentiation of rhesus monkey embryonic stem cells in three-dimensional culture systems. *Stem Cells*. **21**:281-295.
3. Kokenyesi, R., K. Murray, A. Benshushan, E. Huntley, and M. Kao. 2003. Invasion of interstitial matrix by a novel cell line from primary peritoneal carcinosarcoma, and by established ovarian carcinoma cell lines: role of cell-matrix adhesion molecules, proteinases and E-cadherin expression. *Gynecol Oncol*. **89**:60-72.
4. Kutznetsova, N., S. Chi, and S. Leikin. 1998. Sugars and polyols inhibit fibrillogenesis of type I collagen by disrupting hydrogen-bonded water bridges between the helices. *Biochem*. **37**:11888-11895.
5. Kutznetsova, N., and S. Leikin. 1999. Does the triple helical domain of type I collagen encode molecular recognition and fiber assembly while telopeptides serve as catalytic domains. *J. Bio. Chem*. **274**:36083-36088.
6. Leikin, S., D. Rau, and V. Parsegian. 1994. Direct measurement of forces between self-assembled proteins: Temperature-dependent exponential forces between collagen triple helices. *Proc. Natl. Acad. Sci. USA*. **91**:276-280.
7. Leikina, E., M. Meritts, N. Kuznetsova, and S. Leikin. 2002. Type I collagen is thermally unstable at body temperature. *Proc. Natl. Acad. Sci. USA*. **99**:1314-1318.

8. O' Shaughnessy, T., H. Lin, and W. Ma. 2003. Functional synapse formation among rat cortical neurons grown on three-dimensional collagen gels. *Neuroscience Letters*. **340**:169 - 172.
9. Park, D., D. Choi, H. Ryu, H. Kwon, H. Joo, and C. Min. 2003. A well-defined in vitro three-dimensional culture of human endometrium and its applicability to endometrial cancer invasion. *Cancer Letters*. **195**:185-192.
10. Ritty, T., and J. Herzog. 2003. Tendon cells produce gelatinases in response to type I collagen attachment. *J. Ortho. Res.* **21**:442-450.
11. Van Oostveldt, K., M. Paape, and C. Burvemich. 2002. Apoptosis of bovine neutrophils following diapedesis through a monolayer of endothelial and mammary epithelial cells. *J. Dairy Sci. Ass.* **85**:139-147.

Related Products:

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 Factors	5 ml
3433-005-01	Cultrex® BME; no Phenol Red; Reduced Growth Factors	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
3442-050-01	Cultrex® Bovine Collagen I	50 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® Ploy-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

*New Zealand Herd-Derived



Rat Collagen I

Catalog #: 3447-020-01

Storage: 4 °C

(Do Not Freeze)

TREVIGEN®

1-800-873-8443