
Product Manual

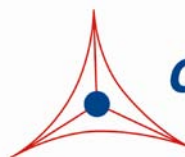
CytoSelect™ 24- Well Cell Contraction Assay Kit (Floating Matrix Model)

Catalog Number

CBA- 5020

24 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

Wound healing is comprised of three processes: epithelialization, connective tissue deposition, and contraction. The contraction process is believed to be mediated by specialized fibroblasts called myofibroblasts. Three-dimensional collagen gels have been widely used to study fibroblast contraction, integrin signaling, cell apoptosis and cytoskeleton reorganization. Since three-dimensional matrix adhesions differ in structure, localization, and function from two-dimensional adhesions; and therefore, three-dimensional cell-matrix interactions may be more relevant biologically.

The signaling mechanisms used by fibroblasts to regulate collagen matrix contraction depend on whether the cells are mechanically loaded or unloaded at the time that contraction is initiated as well as on the growth factor used to initiate contraction. For instance, stimulation of fibroblasts by lysophosphatidic acid (LPA) but not by platelet-derived growth factor (PDGF) causes robust force generation in restrained matrices, whereas LPA and PDGF stimulate floating matrix contraction equally well.

There are several different culture models to study the ability of fibroblasts to reorganize and contract collagen matrices in vitro. In the attached model, a polymerized collagen matrix containing cells remains attached to a culture dish during contraction. Mechanical tension develops during contraction, and cellular stress fibers assemble. The two-step model combines an initial period of attached matrix contraction leading to mechanical loading, followed by release of the matrices, resulting in mechanical unloading and further contraction as mechanical stress dissipates. In the floating matrix contraction model, a freshly polymerized collagen matrix containing cells is released from the culture dish and allowed to float in culture medium, and contraction occurs in the absence of external mechanical load and without appearance of stress fibers in the cells.

Cell Biolabs' CytoSelect™ 24-well Cell Contraction Assay Kit (Floating Matrix Model) provides a simple, in vitro system to assess cell contractivity and screen cell contraction mediators. The kit's proprietary Cell Contraction Plate eliminates the matrix releasing step of the conventional contraction assay, providing a faster, higher-throughput method to assess cell contraction. Each kit contains sufficient quantities to perform up to 24 assays in the provided 24-well plate.

Kit Components

1. 24-Well Cell Contraction Plate (Floating Model) (Part No. 50201B): One 24-well plate (adhesion resistant matrix coated)
2. Collagen Solution (Part No. 20101): One 10 mL sterile bottle of bovine Type I Collagen at 3.0 mg/mL
3. Neutralization Solution (Part No. 20102): One 0.5 mL sterile tube
4. 5X DMEM Medium (Part No. 20103): One 5 mL sterile bottle
5. 5X PBS (Part No. 20104): One 5 mL sterile bottle
6. 100X Cell Contraction Inhibitor (Part No. 20105): One 1 mL sterile tube of 1M 2, 3-Butanedione Monoxime (BDM) in DMSO

Materials Not Supplied

1. Cells such as fibroblasts
2. Cell culture medium
3. 37°C Incubator, 5% CO₂ incubator
4. Light microscope
5. Ruler

Storage

Store all components at 4°C.

Preparation of Collagen Gel Working Solution

This kit is designed for samples in a 24-well plate, and may be modified accordingly to suit other culture plate sizes. Keep all solutions at 4°C during assay preparation. Samples should be assayed at least in duplicate.

1. In a cold sterile tube, add the desired volume of Collagen Solution according to the table below. Next, add corresponding 5X DMEM medium or 5X PBS to the tube, mix well.
2. Add Neutralization solution, IMMEDIATELY mix and keep the Collagen Gel Working Solution on ice. *Note: Try to avoid introducing air bubbles to the mixture.*

Reagents	6 wells	12 wells	24 wells
Collagen Solution	2.385 mL	4.77 mL	9.54 mL
5X Medium or PBS	615 µL	1.23 mL	2.46 mL
Neutralization Solution	85 µL	170 µL	340 µL
Total	3.085 mL	6.17 mL	12.34 mL

Assay Protocol (must be under sterile conditions)

1. Harvest cells and resuspend in desired medium at 2-5 x 10⁶ cells/mL.
2. Prepare the cell contraction matrix by mixing 2 parts of cell suspension and 8 parts of cold Collagen Gel Working Solution.

Note: Try to avoid introducing air bubbles to the mixture. Carefully mix by titrating the solution. Always include negative control wells that contain no cells in the matrix.

3. Add 0.5 mL of the cell contraction matrix to each well of the 24-well Cell Contraction Plate.
4. Incubate the plate at 37°C and 5% CO₂ for 1 hour.
5. After collagen polymerization, carefully add 1.0 mL of culture medium (with/without contraction mediators) atop each collagen gel lattice.

6. Monitor wells for contraction over 2 days at 37°C and 5% CO₂. Media should be changed daily by carefully removing 0.5 mL and replacing with 0.5 mL fresh media (with/without contraction mediators).
7. The collagen gel size change (contraction index) can be measured over time or as a set end point. Results can be quantified with a ruler or with image analysis software, such as NIH Image or Image Pro Plus.

Example of Results

The following figure demonstrates typical contraction results using the Cell Contraction Assay. One should use the data below for reference only. This data should not be used to interpret actual results.

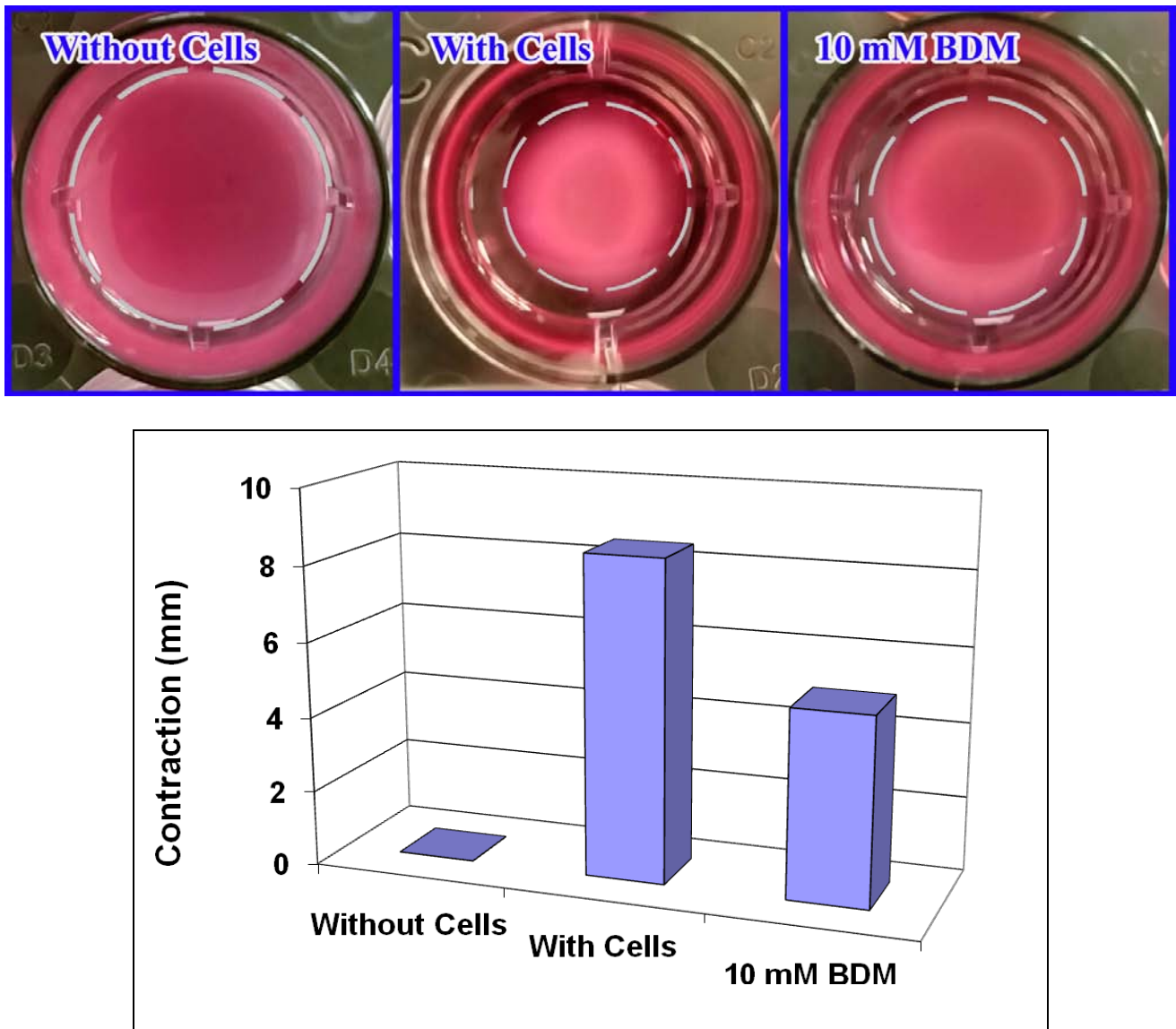


Figure 1. Contraction inhibition by BDM. 0.5×10^6 COS-7 cells in 0.5 mL collagen gel lattice were cultured for two days according to the Assay Protocol (top image). Dashed lines designate the gel edges. The change in matrix diameter size (in millimeters) was determined with a ruler.

References

1. Martin, P. (1997) *Science* **276**, 75-81
2. Bell, E., Ivarsson, B., and Merrill, C. (1979) *Proc. Natl. Acad. Sci. U. S. A.* **76**, 1274-1278
3. Stupak, D., and Harris, A. K. (1982) *Dev. Biol.* **90**, 383-398
4. Mochitate, K., Pawelek, P., and Grinnell, F. (1991) *Exp. Cell Res.* **193**, 198-207
5. Tian, B., Lessan, K., Kahm, J., Kleidon, J., and Henke, C. (2002) *J. Biol. Chem.* **277**, 24667-24675

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2016: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.