

---

Product Manual

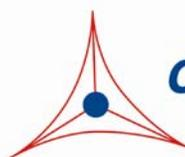
# CytoSelect™ 24- Well Cell Migration Assay (8 $\mu\text{m}$ , Colorimetric Format)

## Catalog Number

CBA- 100	12 assays
CBA- 100- 5	5 x 12 assays

**FOR RESEARCH USE ONLY**  
**Not for use in diagnostic procedures**

---



**CELL BIOLABS, INC.**  
*Creating Solutions for Life Science Research*

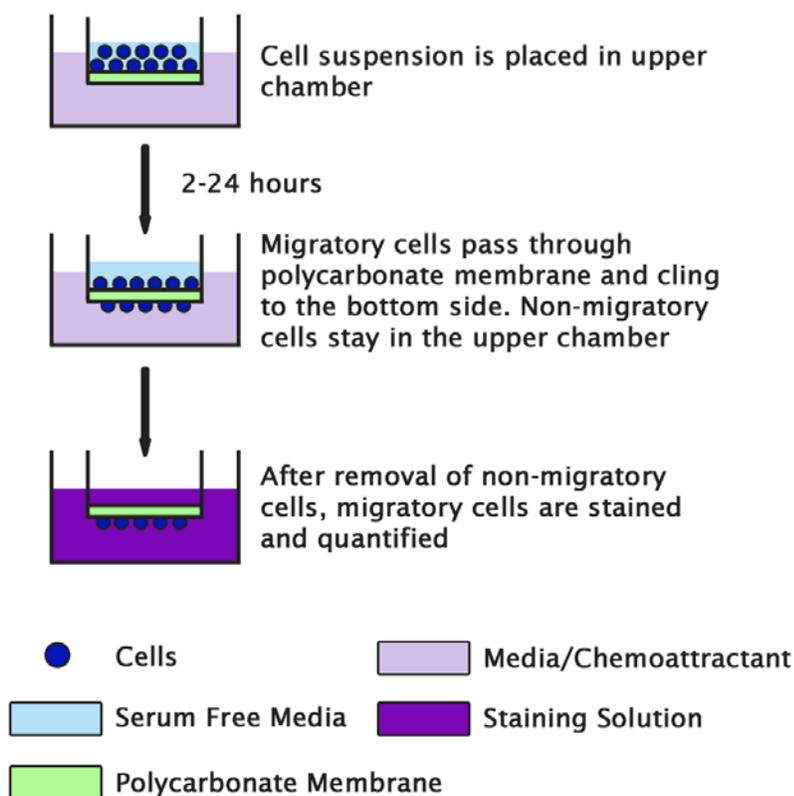
## **Introduction**

Cell migration is a highly integrated, multistep process that orchestrates embryonic morphogenesis, tissue repair and regeneration. It plays a pivotal role in the disease progression of cancer, mental retardation, atherosclerosis, and arthritis. The initial response of a cell to a migration-promoting agent is to polarize and extend protrusions in the direction of the attractant; these protrusions can consist of large, broad lamellipodia or spike-like filopodia. In either case, these protrusions are driven by actin polymerization and can be stabilized by extracellular matrix (ECM) adhesion or cell-cell interactions (via transmembrane receptors).

Cell Biolabs' CytoSelect™ Cell Migration Assay Kit utilizes polycarbonate membrane inserts (8 µm pore size) to assay the migratory properties of cells. The kit contains sufficient reagents for the evaluation of 12 samples. The 8 µm pore size is optimal for epithelial and fibroblast cell migration. However, in the case of leukocyte chemotaxis, a smaller pore size (3 µm) is recommended.

## **Assay Principle**

The CytoSelect™ Cell Migration Assay Kit contains polycarbonate membrane inserts (8 µm pore size) in a 24-well plate. The membrane serves as a barrier to discriminate migratory cells from non-migratory cells. Migratory cells are able to extend protrusions towards chemoattractants (via actin cytoskeleton reorganization) and ultimately pass through the pores of the polycarbonate membrane. Finally, the cells are removed from the top of the membrane and the migratory cells are stained and quantified.



## **Related Products**

1. CBA-100-C: CytoSelect™ 24-Well Cell Migration and Invasion Assay (8µm, Colorimetric)
2. CBA-100-COL: CytoSelect™ 24-Well Cell Haptotaxis Assay (Collagen I, Colorimetric)
3. CBA-100-FN: CytoSelect™ 24-Well Cell Haptotaxis Assay (Fibronectin, Colorimetric)
4. CBA-101: CytoSelect™ 24-Well Cell Migration Assay (8µm, Fluorometric)
5. CBA-102: CytoSelect™ 24-Well Cell Migration Assay (5µm, Fluorometric)
6. CBA-103: CytoSelect™ 24-Well Cell Migration Assay (3µm, Fluorometric)
7. CBA-106: CytoSelect™ 96-Well Cell Migration Assay (8µm, Fluorometric)
8. CBA-110: CytoSelect™ 24-Well Cell Invasion Assay (Basement Membrane, Colorimetric)
9. CBA-125: Radius™ 24-Well Cell Migration Assay (Microscopy)
10. CBA-126: Radius™ 96-Well Cell Migration Assay (Microscopy)
11. CBA-130: CytoSelect™ 96-Well Cell Transformation Assay (Soft Agar Colony Formation)

## **Kit Components**

1. 24-well Migration Plate (Part No. 10001): One 24-well plate containing 12 cell culture inserts (8 µm pore size)
2. Cell Stain Solution (Part No. 11002): One 10 mL bottle
3. Extraction Solution (Part No. 11003): One 10 mL bottle
4. Cotton Swabs (Part No. 11004): 40 each.
5. Forceps (Part No. 11005): One each.

## **Materials Not Supplied**

1. Migratory cell lines
2. Cell culture medium
3. Serum free medium, such as DMEM containing 0.5% BSA, 2 mM CaCl<sub>2</sub> and 2 mM MgCl<sub>2</sub>
4. Cell culture incubator (37°C, 5% CO<sub>2</sub> atmosphere)
5. Light microscope
6. 96-well microtiter plate
7. Microtiter plate reader

## **Storage**

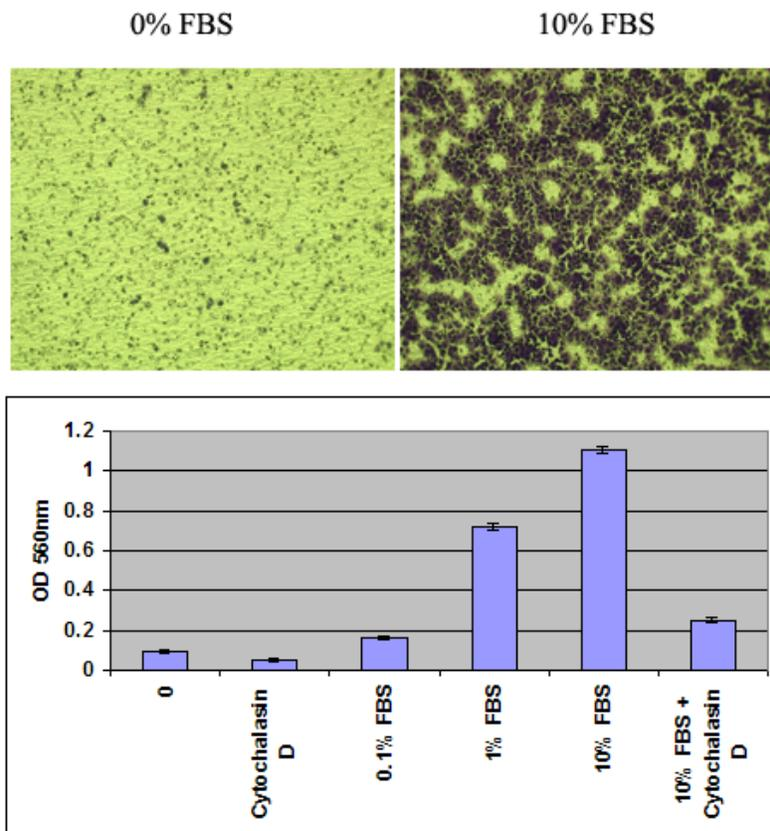
Store all components at 4°C.

## **Assay Protocol**

1. Under sterile conditions, allow the 24-well migration plate to warm up at room temperature for 10 minutes.
2. Prepare a cell suspension containing  $0.5-1.0 \times 10^6$  cells/ml in serum free media. Agents that inhibit or stimulate cell migration can be added directly to the cell suspension.  
*Note: Overnight starvation may be performed prior to running the assay*
3. Add 500  $\mu$ L of media containing 10% fetal bovine serum or desired chemoattractant(s) to the lower well of the migration plate.
4. Add 300  $\mu$ L of the cell suspension solution to the inside of each insert.
5. Incubate for 2-24 hours in a cell culture incubator.
6. Carefully aspirate the media from the inside of the insert. Wet the ends of 2-3 cotton-tipped swabs with water, flatten the ends of the swabs by pressing them against a clean hard surface, and gently swab the interior of the inserts to remove non-migratory cells. Take care not to puncture the polycarbonate membrane. Be sure to remove cells on the inside perimeter of the insert.
7. Transfer the insert to a clean well containing 400  $\mu$ L of Cell Stain Solution and incubate for 10 minutes at room temperature.
8. Gently wash the stained inserts several times in a beaker of water. Allow the inserts to air dry.
9. (optional) Count migratory cells with a light microscope under high magnification objective, with at least three individual fields per insert.
10. Transfer each insert to an empty well, adding 200  $\mu$ L of Extraction Solution per well, then incubating 10 minutes on an orbital shaker.
11. Transfer 100  $\mu$ L from each sample to a 96-well microtiter plate and measure the OD 560nm in a plate reader.

## Example of Results

The following figures demonstrate typical with the CytoSelect™ Cell Migration Assay Kit. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1. Human Fibrosarcoma HT-1080 Cell Migration.** HT-1080 cells were seeded at 150,000 cells/well and allowed to migrate toward FBS for 4 hrs in the presence or absence of 2  $\mu$ M Cytochalasin D. Migratory cells on the bottom of the polycarbonate membrane were stained (top panel picture) and quantified at OD 560nm after extraction (bottom panel figure).

## References

1. Ridley AJ, Schwartz MA, Burridge K, Firtel RA, Ginsberg MH, Borisy G, Parsons JT, Horwitz AR. (2003) *Science* **302**, 1704-9.
2. Horwitz R, Webb D. (2003) *Curr Biol.* **13**, R756-9.
3. Lauffenburger DA, Horwitz AF. (1996) *Cell* **84**, 359-369.

## **Recent Product Citations**

1. Yamauchi, K., et al. (2017). Structure-Activity Relationships of Methylquercetin on Anti-migration and Anti-proliferation Activity in B16 Melanoma Cells. *Anticancer Res.* **37**(4):1575-1579.
2. Colden, M. et al. (2017). MicroRNA-466 inhibits tumor growth and bone metastasis in prostate cancer by direct regulation of osteogenic transcription factor RUNX2. *Cell Death Dis.* **8**(1):e2572. doi: 10.1038/cddis.2017.15.
3. Afghani, N. et al. (2017). Microtubule actin cross-linking factor 1, a novel target in glioblastoma. *Int. J. Oncol.* **50**:310-316.
4. Yu, X. M. et al. (2016). Notch1 signaling regulates the aggressiveness of differentiated thyroid cancer and inhibits SERPINE1 expression. *Clin Cancer Res.* doi:10.1158/1078-0432.CCR-15-1749.
5. Kestens, C. et al. (2016). BMP4 Signaling is able to induce an epithelial-mesenchymal transition-like phenotype in Barrett's esophagus and esophageal adenocarcinoma through induction of SNAIL2. *PLoS One.* **11**:e0155754.
6. Kim, J. et al. (2016). Vimentin filaments regulate integrin–ligand interactions by binding to the cytoplasmic tail of integrin  $\beta$ 3. *J Cell Sci.* **129**:2030-2042.
7. Slusser-Nore, A. et al. (2016). SPARC expression is selectively suppressed in tumor initiating urospheres isolated from As+ 3-and Cd+ 2-transformed human urothelial cells (UROtsa) stably transfected with SPARC. *PLoS One.* **11**:e0147362.
8. Paris, J. L. et al. (2016). Decidua-derived mesenchymal stem cells as carriers of mesoporous silica nanoparticles. In vitro and in vivo evaluation on mammary tumors. *Acta Biomater.* doi:10.1016/j.actbio.2016.01.017.
9. Barui, S. et al. (2015). Systemic co-delivery of a homo-serine derived ceramide analog and curcumin to tumor vasculature inhibits mouse tumor growth. *Mol Pharm.* doi:10.1021/acs.molpharmaceut.5b00644.
10. Wang, K. P. et al. (2015). Morphine protects SH-SY5Y human neuroblastoma cells against Dickkopf1-induced apoptosis. *Mol Med Rep.* **11**:1174-1180.
11. Song, G. Q. & Zhao, Y. (2015). Kisspeptin-10 inhibits the migration of breast cancer cells by regulating epithelial-mesenchymal transition. *Oncol Rep.* **33**:669-674.
12. Ran, X. et al. (2015). A quantitative proteomics study on olfactomedin 4 in the development of gastric cancer. *Int J Oncol.* **47**:1932-1944.
13. Wang, C. J. et al. (2015). RhoBTB2 (DBC2) functions as a multifunctional tumor suppressor in thyroid cancer cells via mitochondrial apoptotic pathway. *Int J Clin Exp Med.* **8**:5954-5958.
14. König, J. et al. (2015). Placental mesenchymal stromal cells derived from blood vessels or avascular tissues: what is the better choice to support endothelial cell function? *Stem Cells Dev.* **24**:115-131.
15. Bliley, J. M. et al. (2015). Ethylene oxide sterilization preserves bioactivity and attenuates burst release of encapsulated glial cell line derived neurotrophic factor from tissue engineering nerve guides for long gap peripheral nerve repair. *ACS Biomater Sci Eng.* doi:10.1021/ab5001518.
16. Choi, S. I. et al. (2014). TM4SF4 overexpression in radiation-resistant lung carcinoma cells activates IGF1R via elevation of IGF1. *Oncotarget.* **5**:9823.
17. Kwon, S. M. et al. (2014). Cross talk with hematopoietic cells regulates the endothelial progenitor cell differentiation of CD34 positive cells. *PLoS One.* **9**:e106310.
18. Kanan, Y. et al. (2014). Fibulin 2, a tyrosine O-sulfated protein, is up-regulated following retinal detachment. *J Biol Chem.* **289**:13419-13433.

19. Al-Wadei, M.H. et al. (2013). Gamma-amino Butyric Acid (GABA) prevents the induction of nicotinic receptor-regulated signaling by chronic ethanol in pancreatic cancer cells and normal duct epithelia. *Cancer Prevention Research*. **6**:139-148.
20. Al-Wadei, M.H. et al. (2012). Effects of chronic nicotine on the autocrine regulation of pancreatic cancer cells and pancreatic duct epithelial cells by stimulatory and inhibitory neurotransmitters. *Carcinogenesis*. **33**:1745-7153.
21. Chen, Z. et al. (2012). The iron chelators Dp44mT and DFO inhibit TGF- $\beta$ -induced epithelial-mesenchymal transition via up-regulation of N-Myc downstream-regulated gene 1 (NDRG1). *J. Biol. Chem.* **287**:17016-17028.
22. Isakova, I. et al. (2007). Age and dose-related effects on MSC engraftment levels and anatomical distribution in the CNS of non-human primates: Identification of novel MSC subpopulations that respond to guidance cues in brain. *Stem Cells* **25**: 3261-3270.
23. Schuller, H. et al. (2007). GABA B receptor is a novel drug target for pancreatic cancer. *Cancer* **112(4)**:767-768.
24. Eckstein, N. et al. (2008). EGFR-pathway analysis identifies amphiregulin as a key factor for cisplatin resistance of human breast cancer cells. *J. Biol. Chem.* **283**:739-750.
25. Liu, S. et al. (2008). Matrix metalloproteinase-activated anthrax lethal toxin demonstrates high potency in targeting tumor vasculature. *J. Biol. Chem.* **283**:529-540.
26. Awasthi, N. et al. (2008). Downregulation of MMP-2 and -9 by proteasome inhibition: A possible mechanism to decrease LEC migration and prevent posterior capsular opacification. *Invest. Ophthalmol. Vis. Sci.* **49**:1998-2003.

Please see the complete list of product citations: <http://www.cellbiolabs.com/24-well-chemotaxis-assays-8-um>.

## **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](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2004-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.