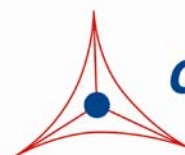

Product Manual

CytoSelect™ 24- Well Cell Migration and Invasion Assay (8 μ m, Colorimetric Format)

Catalog Number

CBA- 100- C	2 x 12 assays (12 migration + 12 invasion)
CBA- 100- C- 5	10 x 12 assays (60 migration + 60 invasion)

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).

The ability of malignant tumor cells to invade normal surrounding tissue contributes in large part to the significant morbidity and mortality of cancers. Invasiveness requires several distinct cellular functions including adhesion, motility, detachment, and extracellular matrix proteolysis. Metastatic cells produce many proteolytic enzymes (e.g. lysosomal hydrolysates, collagenases, plasminogen activators) while the expression of certain cell surface protease receptors is also increased.

Cell Biolabs' CytoSelect™ Cell Migration Assay utilizes polycarbonate membrane inserts (8 µm pore size) to assay the migratory properties of cells. 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. Cell Biolabs' CytoSelect™ Cell Invasion Assay utilizes basement membrane-coated inserts to assay the invasive properties of tumor cells. Each assay contains sufficient reagents for the evaluation of 12 samples.

Related Products

1. CBA-100: CytoSelect™ 24-Well Cell Migration 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-101-C: CytoSelect™ 24-Well Cell Migration and Invasion Assay Combo Kit (8 µm, Fluorometric)
6. CBA-102: CytoSelect™ 24-Well Cell Migration Assay (5µm, Fluorometric)
7. CBA-103: CytoSelect™ 24-Well Cell Migration Assay (3µm, Fluorometric)
8. CBA-106: CytoSelect™ 96-Well Cell Migration Assay (8µm, Fluorometric)
9. CBA-106-C: CytoSelect™ 96-Well Cell Migration and Invasion Assay Combo Kit (8µm, Fluorometric)
10. CBA-110: CytoSelect™ 24-Well Cell Invasion Assay (Basement Membrane, Colorimetric)
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. Invasion Chamber Plate (Part No. 11001): One 24-well plate containing 12 ECM-coated cell culture inserts.
3. Cell Stain Solution (Part No. 11002-C): One 20 mL bottle
4. Extraction Solution (Part No. 11003-C): One 20 mL bottle
5. Cotton Swabs (Part No. 11004): 40 each
6. Forceps: (Part No. 11005) One each

Materials Not Supplied

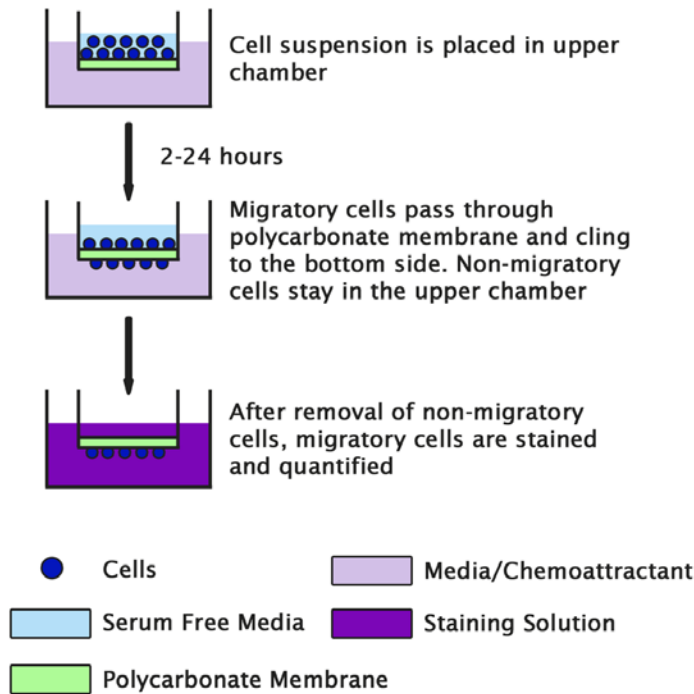
1. Migratory or invasive cell lines
2. Cell culture medium
3. Serum free medium, such as DMEM containing 0.5% BSA, 2 mM CaCl_2 and 2 mM MgCl_2
4. Cell culture incubator (37°C, 5% CO_2 atmosphere)
5. Light microscope
6. 96-well microtiter plate

Storage

Store all components at 4°C.

Cell Migration Assay Principle

The Cell Migration portion of this kit uses 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.



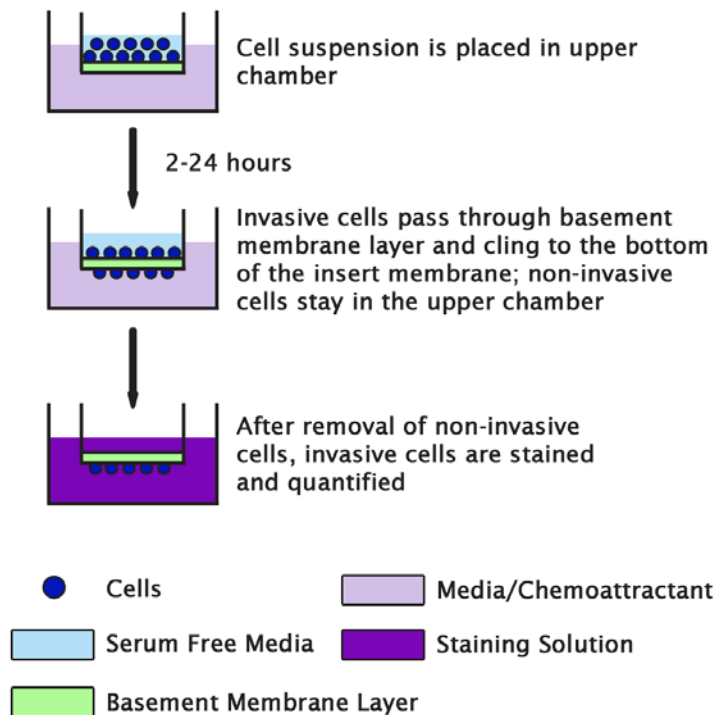
Cell Migration Assay Protocol

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 µL of media containing 10% fetal bovine serum or desired chemoattractant(s) to the lower well of the migration plate.
4. Add 300 µL 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 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 µL 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 μL of Extraction Solution per well, then incubating 10 minutes on an orbital shaker.
11. Transfer 100 μL from each sample to a 96-well microtiter plate and measure the OD 560nm in a plate reader.

Cell Invasion Assay Principle

The Cell Invasion Assay portion of this kit uses a 24-well plate containing polycarbonate membrane inserts (8 μm pore size); the upper surface of the insert membrane is coated with a uniform layer of dried basement membrane matrix solution. This basement membrane layer serves as a barrier to discriminate invasive cells from non-invasive cells. Invasive cells are able to degrade the matrix proteins in the layer, and ultimately pass through the pores of the polycarbonate membrane. Finally, the cells are removed from the top of the membrane and the invaded cells are stained and quantified.



Cell Invasion Assay Protocol

1. Under sterile conditions, allow the invasion chamber plate to warm up at room temperature for 10 minutes.
2. Rehydrate the basement membrane layer of the cell culture inserts by adding 300 μL of warm, serum-free media to the inner compartment. Incubate at room temperature for 1 hour.
3. Prepare a cell suspension containing $0.5\text{-}1.0 \times 10^6$ cells/ml in serum free media. Agents that inhibit or stimulate cell invasion can be added directly to the cell suspension.

Note: Overnight starvation may be performed prior to running the assay

4. Carefully remove the rehydration medium (step 2) from the inserts without disturbing the basement membrane layer.
Note: It will not affect the assay performance if a small amount of rehydration medium is left in the compartment
5. Add 500 μL of media containing 10% fetal bovine serum or desired chemoattractant(s) to the lower well of the migration plate.
6. Add 300 μL of the cell suspension solution to the inside of each insert.
7. Incubate for 12-48 hours in a cell culture incubator.
8. 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-invasive cells. Take care not to puncture the polycarbonate membrane. Be sure to remove cells on the inside perimeter of the insert.
9. Transfer the insert to a clean well containing 400 μL of Cell Stain Solution and incubate for 10 minutes at room temperature.
10. Gently wash the stained inserts several times in a beaker of water. Allow the inserts to air dry.
11. (optional) Count invasive cells with a light microscope under high magnification objective, with at least three individual fields per insert.
12. Transfer each insert to an empty well, adding 200 μL of Extraction Solution per well, then incubating 10 minutes on an orbital shaker.
13. Transfer 100 μL 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 and Invasion Assay Kit. One should use the data below for reference only. This data should not be used to interpret actual results.

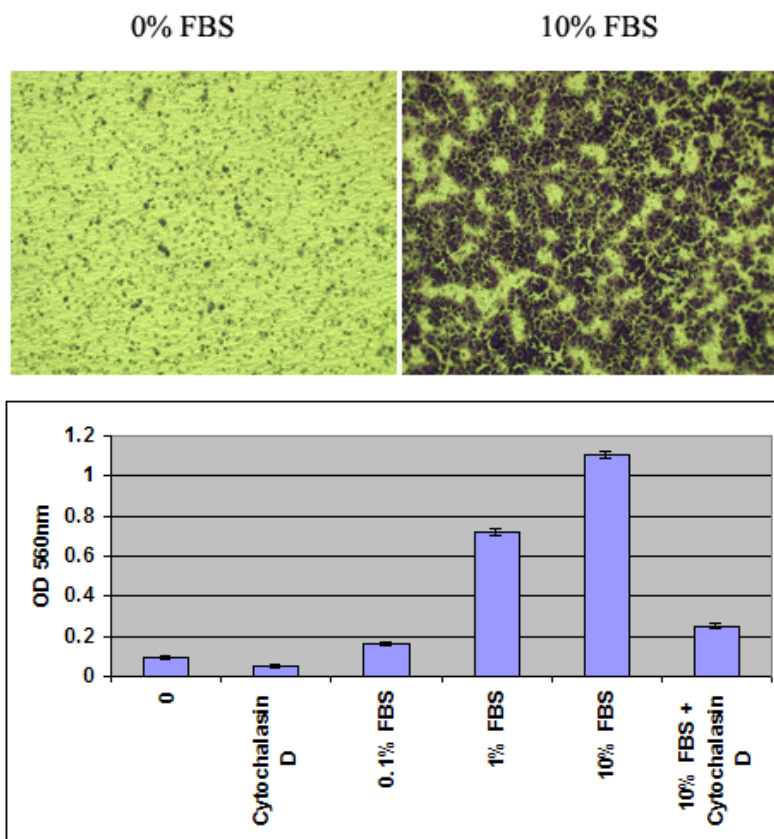


Figure 1. Human Fibrosarcoma HT-1080 Cell Migration. HT-1080 cells were seeded at 150,000 cells/well and allowed to migrate toward FBS for 4 hrs in the presence or absence of 2 μ M Cytochalasin D. Migratory cells on the bottom of the polycarbonate membrane were stained (top panel picture) and quantified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

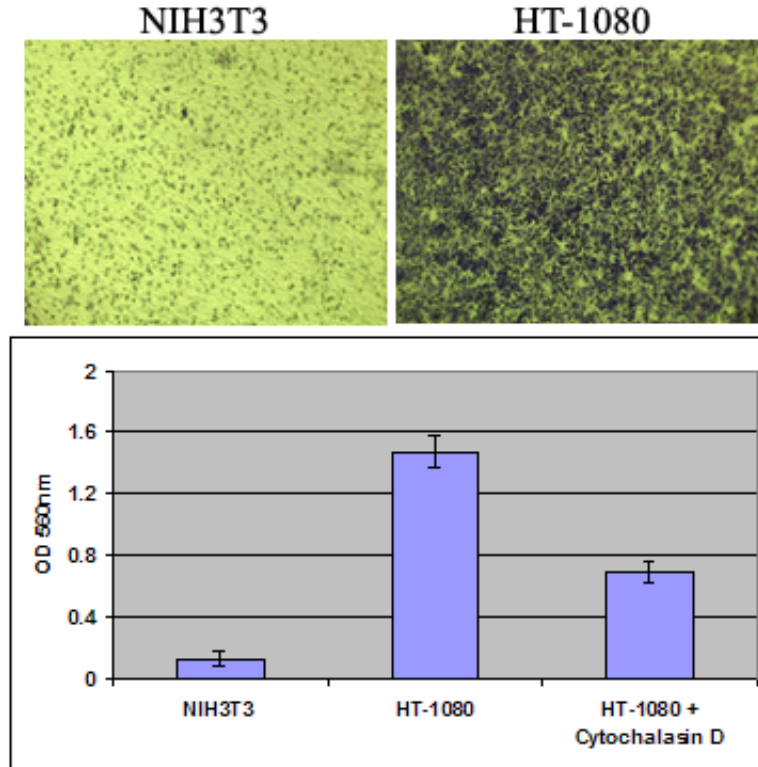


Figure 2. Human Fibrosarcoma HT-1080 Cell Invasion. HT-1080 and NIH3T3 (negative control) were seeded at 300,000 cells/well and allowed to invade toward FBS for 24 hrs in the presence or absence of 2 μ M Cytochalasin D. Invasive cells on the bottom of the invasion membrane were stained (top panel picture) 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.
4. Erkell, L. J., Schirmacher, V. (1988) *Cancer Res* **48**, 6933-6937.
5. Montgomery, A. M. P., De Clerck, Y. A., Langley, K. E., Reisfeld, R. A., Mueller, B. M. (1993) *Cancer Res* **53**,693-700.
6. Monsky, W. L., Lin, C. Y., Aoyama, A., Kelly, T., Akiyama, S. K., Mueller, S. C., Chen, W. T. (1994) *Cancer Res* **54**,5702-5710.

Recent Product Citations

1. Demirci, S. et al. (2017). Cytoglobin inhibits migration through PI3K/AKT/mTOR pathway in fibroblast cells. *Mol. Cell Biochem.* doi:10.1007/s11010-017-3101-2.
2. Pak, J.H. et al. (2017). Clonorchis sinensis excretory-secretory products promote the migration and invasion of cholangiocarcinoma cells by activating the integrin β 4-FAK/Src signaling pathway. *Mol Biochem Parasitol.* **214**:1-9. doi: 10.1016/j.molbiopara.2017.03.002.
3. Coban, E.A. et al. (2017). Characterization of stem-like cells in a new astroblastoma cell line. *Exp Cell Res.* **352**(2):393-402. doi: 10.1016/j.yexcr.2017.02.030.
4. Gong, J. et al. (2017). Inhibition of FASN suppresses migration, invasion and growth in hepatoma carcinoma cells by deregulating the HIF-1 α /IGFBP1 pathway. *Int J Oncol.* **50**(3):883-892. doi: 10.3892/ijo.2017.3867
5. Quinn, S. et al. (2016). Adenylyl cyclase 3/adenylyl cyclase-associated protein 1 (CAP1) complex mediates the anti-migratory effect of forskolin in pancreatic cancer cells. *Mol. Carcinog.* doi:10.1002/mc.22598.
6. Komina, A. et al. (2016). Antiproliferative and pro-apoptotic effects of miR-4286 inhibition in melanoma cells. *PLoS One* **11**:e0168229.
7. Nahar, S. et al. (2016). Systematic evaluation of biophysical and functional characteristics of selenomethylene locked nucleic acid mediated inhibition of miR-21. *Biochemistry* doi:10.1021/acs.biochem.6b00895.
8. Airoidi, I. et al. (2016). Interleukin-30 promotes breast cancer growth and progression. *Cancer Res.* **76**:6218-6229.
9. Li, X. S. & He, X. L. (2016). Kallikrein 12 downregulation reduces AGS gastric cancer cell proliferation and migration. *Genet Mol Res.* doi:10.4238/gmr.15038452.
10. Li, Y. et al. (2016). p27 is a candidate prognostic biomarker and metastatic promoter in osteosarcoma. *Cancer Res.* doi:10.1158/0008-5472.CAN-15-3189.
11. Du, L. et al. (2016). Transcriptome profiling reveals novel gene expression signatures and regulating transcription factors of TGF β -induced epithelial-to-mesenchymal transition. *Cancer Med.* doi:10.1002/cam4.719.
12. Costa, H. et al. (2016). Human cytomegalovirus may promote tumour progression by upregulating arginase-2. *Oncotarget.* doi:10.18632/oncotarget.9722.
13. Yoo, B. et al. (2015). Combining miR-10b-targeted nanotherapy with low-dose doxorubicin elicits durable regressions of metastatic breast cancer. *Cancer Res.* **75**:4407-4415.
14. Bose, D. et al. (2015). Selective inhibition of miR-21 by phage display screened peptide. *Nucleic Acids Res.* doi: 10.1093/nar/gkv185.
15. Cai, X. Z. et al. (2015). iTRAQ-based quantitative proteomic analysis of nasopharyngeal carcinoma. *J Cell Biochem.* doi: 10.1002/jcb.25105.
16. Zheng, X. et al. (2015). Targeting LunX inhibits non-small-cell lung cancer growth and metastasis. *Cancer Res.* doi: 10.1158/0008-5472.
17. Kim, K. S. et al. (2015). Increased expression of endocan in arthritic synovial tissues: Effects of adiponectin on the expression of endocan in fibroblast-like synoviocytes. *Mol Med Rep.* **11**:2695-2702.
18. Liu, Y. et al. (2014). The role of von Willebrand factor as a biomarker of tumor development in hepatitis B virus-associated human hepatocellular carcinoma: A quantitative proteomic based study. *J Proteomics.* **106**:99-112.
19. Chavali, P. L. et al. (2014). TLX activates MMP-2, promotes self-renewal of tumor spheres in neuroblastoma and correlates with poor patient survival. *Cell Death Dis.* **5**:e1502.

20. Barui, S. et al. (2014). Simultaneous delivery of doxorubicin and curcumin encapsulated in liposomes of pegylated RGDK-lipo-peptide to tumor vasculature. *Biomaterials*. **35**:1643-1656.
21. Peng, Y. et al. (2014). microRNA-503 inhibits gastric cancer cell growth and epithelial-to-mesenchymal transition. *Oncol Lett*. **7**:1233-1238.
22. Fatemi, M. et al. (2014). Epigenetic silencing of CHD5, a novel tumor-suppressor gene, occurs in early colorectal cancer stages. *Cancer*. **120**:172-180.
23. Li, M. et al. (2013). Physiological role of the interaction between CARMIL1 and capping protein. *Hum. Reprod*. **28**:2822-2831.
24. Majid, S. et al. (2013). miRNA-34b inhibits prostate cancer through demethylation, active chromatin modifications, and AKT pathways. *Clin. Cancer Res*. **19**: 73-84.
25. Majid, S. et al. (2012). miR-23b represses proto-oncogene Src kinase and functions as methylation-silenced tumor suppressor with diagnostic and prognostic significance in prostate cancer. *Cancer Res*. **72**:6435-6446.
26. Shin, S.Y. et al. (2012). Transcriptional regulation of the Interleukin-11 gene by oncogenic Ras. *Carcinogenesis* 10.1093/carcin/bgs297.
27. Saini, S. et al. (2012). miRNA-708 control of CD44+ prostate cancer-initiating cells. *Cancer Res*. **72**: 3618-3630.
28. Dennis, M. et al. (2012). Snail controls the mesenchymal phenotype and drives erlotinib resistance in oral epithelial and head and neck squamous cell carcinoma cells. *Otolaryngology--Head and Neck Surgery*. **147**: 726-732.
29. Dalezis, P. et al. (2012). Dexamethasone plus octreotide regimen increases anticancer effects of docetaxel on TRAMP-C1 prostate cancer model. *In Vivo* **26**:75-86.
30. Kahlert, C. et al. (2011). Overexpression of ZEB2 at the invasion front of colorectal cancer is an independent prognostic marker and regulates tumor invasion *in vitro*. *Clin. Cancer Res*. **17**:7654-7663.

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

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