
Product Manual

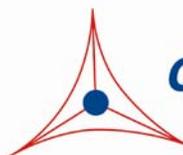
CytoSelect™ 24- Well Cell Migration Assay (8 μm, Fluorometric Format), Trial Size

Catalog Number

CBA- 101- T

4 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

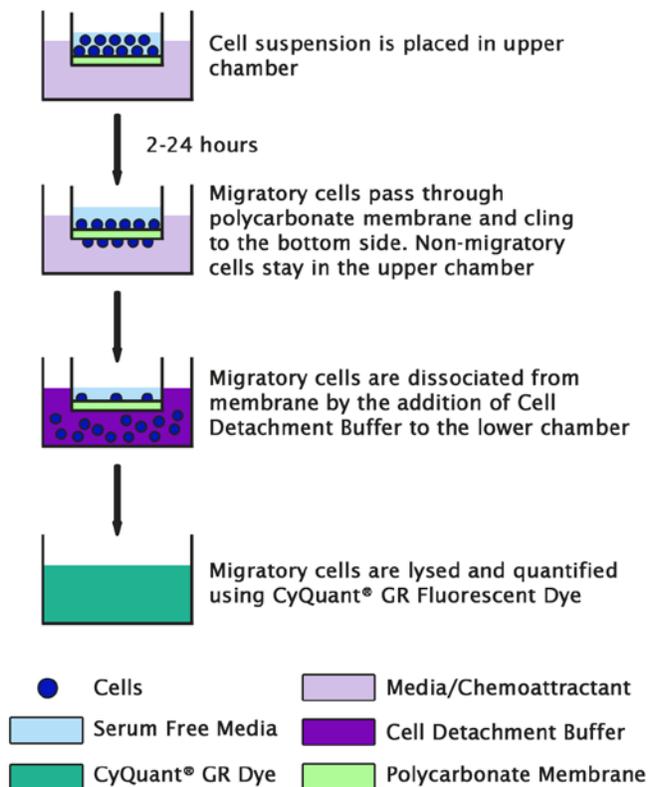
Cell migration is a highly integrated, multistep process that orchestrates embryonic morphogenesis, tissue repair and regeneration. It plays a pivotal role in the disease progression of cancer, mental retardation, atherosclerosis, and arthritis. The initial response of a cell to a migration-promoting 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 does not require you to prelabel the cells with Calcein AM or remove non-migratory cells (i.e. cotton swabbing). Any migratory cells are first dissociated from the membrane, then lysed and detected by the patented CyQuant® GR Dye (Invitrogen).

Cell Biolabs CytoSelect™ Cell Migration Assay Kit provides a robust system for the quantitative determination of cell migration. This Trial Size kit contains sufficient reagents for the evaluation of 4 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.

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. These migratory cells are then dissociated from the membrane and subsequently detected by the patented CyQuant® GR Dye (Invitrogen).

Assay Principle



Related Products

1. CBA-100: CytoSelect™ 24-Well Cell Migration Assay (8µm, Colorimetric)
2. CBA-101-COL: CytoSelect™ 24-Well Cell Haptotaxis Assay (Collagen I, Fluorometric)
3. CBA-101-FN: CytoSelect™ 24-Well Cell Haptotaxis Assay (Fibronectin, Fluorometric)
4. CBA-102: CytoSelect™ 24-Well Cell Migration Assay (5µm, Fluorometric)
5. CBA-103: CytoSelect™ 24-Well Cell Migration Assay (3µm, Fluorometric)
6. CBA-104: CytoSelect™ 96-Well Cell Migration Assay (3 µm, 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-T): One 24-well plate containing 4 cell culture inserts (8 μm pore size)
2. Cell Detachment Solution (Part No. 10101-T): One 2 mL tube
3. 4X Lysis Buffer (Part No. 10102-T): One 2 mL tube
4. CyQuant® GR Dye (Part No. 10103-T): One 10 μL tube
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_2 and 2 mM MgCl_2
4. Cell culture incubator (37°C, 5% CO_2 atmosphere)
5. Light microscope
6. 96-well plate suitable for a fluorescence plate reader
7. Fluorescence 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 $0.5\text{-}1.0 \times 10^6$ 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. Transfer the insert to a clean well containing 225 μL of Cell Detachment Solution. Incubate 30 minutes at 37°C.
7. Completely dislodge the cells from the underside of the membrane by gently tilting the insert several times in the detachment solution. Remove and discard the insert.
8. Prepare sufficient 4X Lysis Buffer/CyQuant® GR dye solution for all samples by diluting the dye 1:75 in 4X Lysis Buffer (for example, add 5 μL dye to 370 μL of 4X Lysis Buffer).

9. Add 75 μL of 4X Lysis Buffer/CyQuant® GR dye solution to each well containing cells and 225 μL of Cell Detachment Solution. Incubate 20 minutes at room temperature.
10. Transfer 200 μL of the mixture to a 96-well plate suitable for fluorescence measurement. Read fluorescence with a fluorescence plate reader at 480 nm/520 nm.

Example of Results

The following figures demonstrate typical with the CytoSelect™ Cell Migration Assay Kit. Fluorescence measurement was performed on SpectraMax Gemini XS Fluorometer (Molecular Devices) with a 485/538 nm filter set and 530 nm cutoff. One should use the data below for reference only. This data should not be used to interpret actual results.

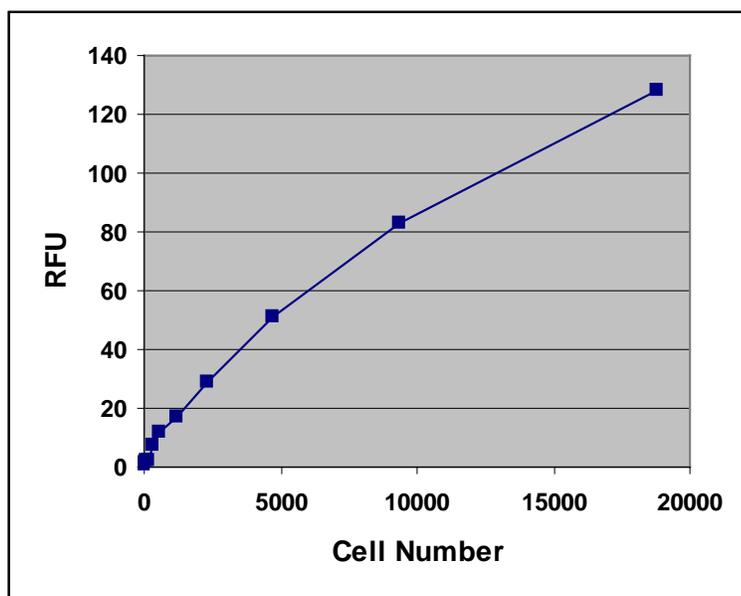


Figure 1: Quantitation of Human HT-1080. HT-1080 cells were titrated in Cell Detachment Buffer, then subsequently lysed and detected with 4X Lysis Buffer/Cyquant® GR Dye (150 μL cell suspension was mixed with 50 μL of 4X Lysis Buffer/dye).

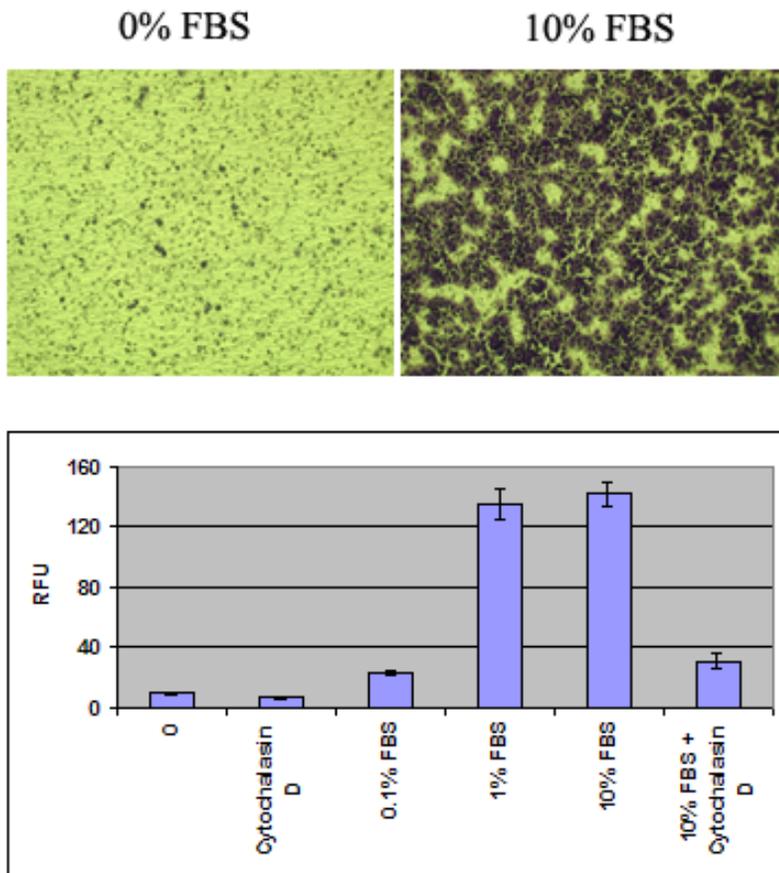


Figure 2. Human Fibrosarcoma HT-1080 Cell Migration. HT-1080 was seeded at 30,000 cells/well and allowed to migrate toward FBS for 4 hrs in the presence or absence of 2 μ M Cytochalasin D. Migratory cells on the bottom of the polycarbonate membrane were stained (top panel picture) and quantified by CyQuant® GR Dye as described in the Assay Protocol.

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. Ibrahim, S.A. et al. (2016). Cancer derived peptide of vacuolar ATPase 'a2' isoform promotes neutrophil migration by autocrine secretion of IL-8. *Sci. Rep.* **6**:36865.
2. Słoniecka, M. et al. (2015). Substance P enhances keratocyte migration and neutrophil recruitment through interleukin-8. *Mol Pharmacol.* doi:10.1124/mol.115.101014.
3. Banerjee, D. et al. (2015). Notch suppresses angiogenesis and progression of hepatic metastases. *Cancer Res.* **75**:1592-1602.

- Izhak, L. et al. (2010). Predominant expression of CCL2 at the tumor site of prostate cancer patients directs a selective loss of immunological tolerance to CCL2 that could be amplified in a beneficial manner. *J. Immunol.* **184**:1092-1101.
- Wang, W. et al. (2009). Netrin-1 increases proliferation and migration of renal proximal tubular epithelial cells via the UNC5B receptor. *Am. J. Physiol. Renal Physiol.* **296**:F723-729.
- Izhak, L. et al. (2009). A novel recombinant fusion protein encoding a 20-amino acid residue of the third extracellular (E3) domain of CCR2 neutralizes the biological activity of CCL2. *J. Immunol.* **183**:732-739.
- Zhao, L. et al. (2008). CD44 expressed on both bone marrow-derived and non-bone marrow-derived cells promotes atherosclerosis in ApoE-deficient mice. *Arterioscler. Thromb. Vasc. Biol.* **28**:1283-1289.
- Fava, G. et al. (2008). Leptin enhances cholangiocarcinoma cell growth. *Cancer Res.* **68**:6752-6761.
- Uddin, M. et al. (2008). Marinobufagenin inhibits proliferation and migration of cytotrophoblast and CHO cells. *Placenta* **29(3)**:266-273.
- Reing, J. et al. (2008). Degradation products of extracellular matrix affect cell migration and proliferation. *Tissue Eng. Part A* 10.1089/ten.tea.2007.0425.

License Information

This product is provided under an intellectual property license from Life Technologies Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased product and components of the product only in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The sale of this product is expressly conditioned on the buyer not using the product or its components, or any materials made using the product or its components, in any activity to generate revenue, which may include, but is not limited to use of the product or its components: (i) in manufacturing; (ii) to provide a service, information, or data in return for payment; (iii) for therapeutic, diagnostic or prophylactic purposes; or (iv) for resale, regardless of whether they are resold for use in research. For information on purchasing a license to this product for purposes other than as described above, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad CA 92008 USA or outlicensing@lifetech.com.

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

©2013-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.