

---

Product Manual

# CytoSelect™ 24-Well Cell Invasion Assay (Collagen I, Fluorometric Format)

Catalog Number

CBA-111-COL

12 assays

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

---



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

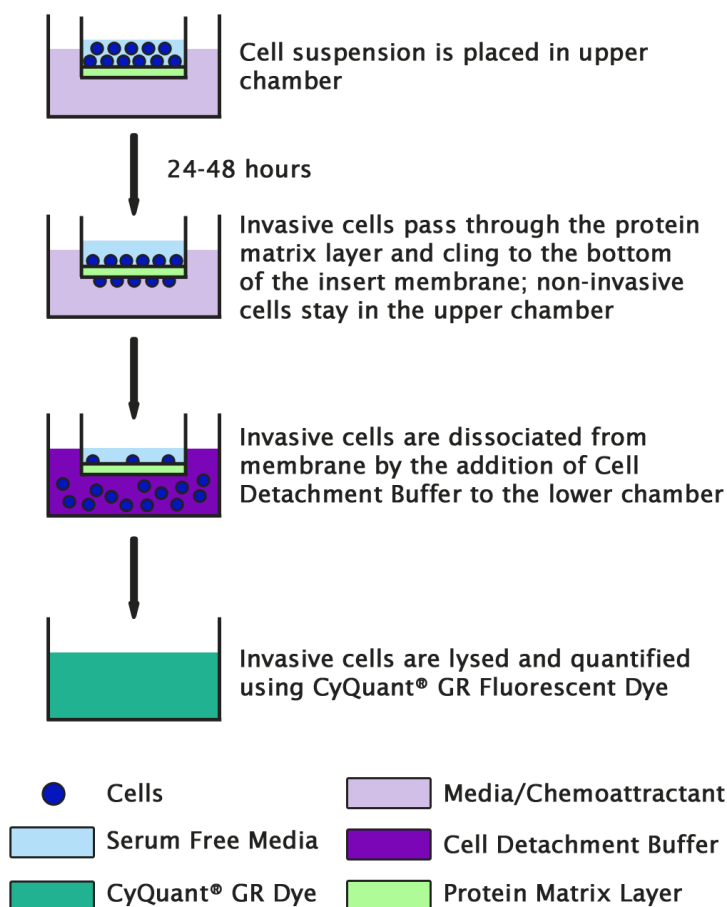
## **Introduction**

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™ Collagen Cell Invasion Assay Kit utilizes Bovine Type I Collagen-coated inserts to assay the invasive properties of tumor cells. It contains sufficient reagents for the evaluation of 12 samples.

## **Assay Principle**

The CytoSelect™ Collagen Cell Invasion Assay Kit contains polycarbonate membrane inserts (8 µm pore size) in a 24-well plate. The upper surface of the insert membrane is coated with a uniform layer of dried Bovine Type I Collagen matrix. This collagen matrix layer serves as a barrier to discriminate invasive cells from non-invasive cells. Invasive cells are able to degrade the collagen matrix layer, and ultimately pass through the pores of the polycarbonate membrane. Finally, these invaded cells are then dissociated from the membrane and subsequently detected by the patented CyQuant® GR Dye (Invitrogen).



## **Related Products**

1. CBA-106: CytoSelect™ 96-Well Cell Migration Assay (8µm, Fluorometric)
2. CBA-106-C: CytoSelect™ 96-Well Cell Migration and Invasion Assay (8µm, Fluorometric)
3. CBA-110: CytoSelect™ 24-Well Cell Invasion Assay (Basement Membrane, Colorimetric)
4. CBA-110-COL: CytoSelect™ 24-Well Cell Invasion Assay (Collagen I, Colorimetric)
5. CBA-110-LN: CytoSelect™ 24-Well Cell Invasion Assay (Laminin, Colorimetric)
6. CBA-111: CytoSelect™ 24-Well Cell Invasion Assay (Basement Membrane, Fluorometric)
7. CBA-111-LN: CytoSelect™ 24-Well Cell Invasion Assay (Laminin, Fluorometric)
8. CBA-112: CytoSelect™ 96-Well Cell Invasion Assay (Basement Membrane, Fluorometric)
9. CBA-112-COL: CytoSelect™ 96-Well Cell Invasion Assay (Collagen I, Fluorometric)
10. CBA-112-LN: CytoSelect™ 96-Well Cell Invasion Assay (Laminin, Fluorometric)
11. CBA-130: CytoSelect™ 96-Well Cell Transformation Assay (Soft Agar Colony Formation)

## **Kit Components**

1. Collagen Invasion Chamber Plate (Part No. 111001-COL): One 24-well plate containing 12 collagen-coated cell culture inserts
2. Cell Detachment Solution (Part No. 10101): One 5 mL bottle
3. 4X Lysis Buffer (Part No. 10102): One 5 mL bottle
4. CyQuant® GR Dye (Part No. 10103): One 25 µL tube
5. Forceps (Part No. 11005): One each

## **Materials Not Supplied**

1. Invasive 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 plate suitable for a fluorescence plate reader
7. Fluorescence plate reader

## **Storage**

Store all components at 4°C.

## **Assay Protocol**

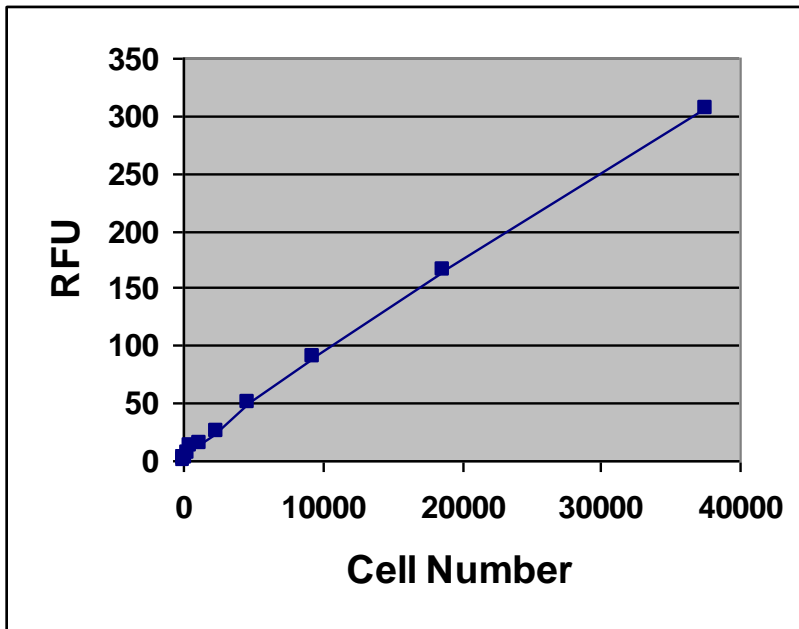
1. Under sterile conditions, allow the collagen invasion chamber plate to warm up at room temperature for 10 minutes.
2. Rehydrate the collagen layer of the cell culture inserts by adding 300  $\mu\text{L}$  of warm, serum-free media to the inner compartment. Incubate at room temperature for 30 minutes.
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*

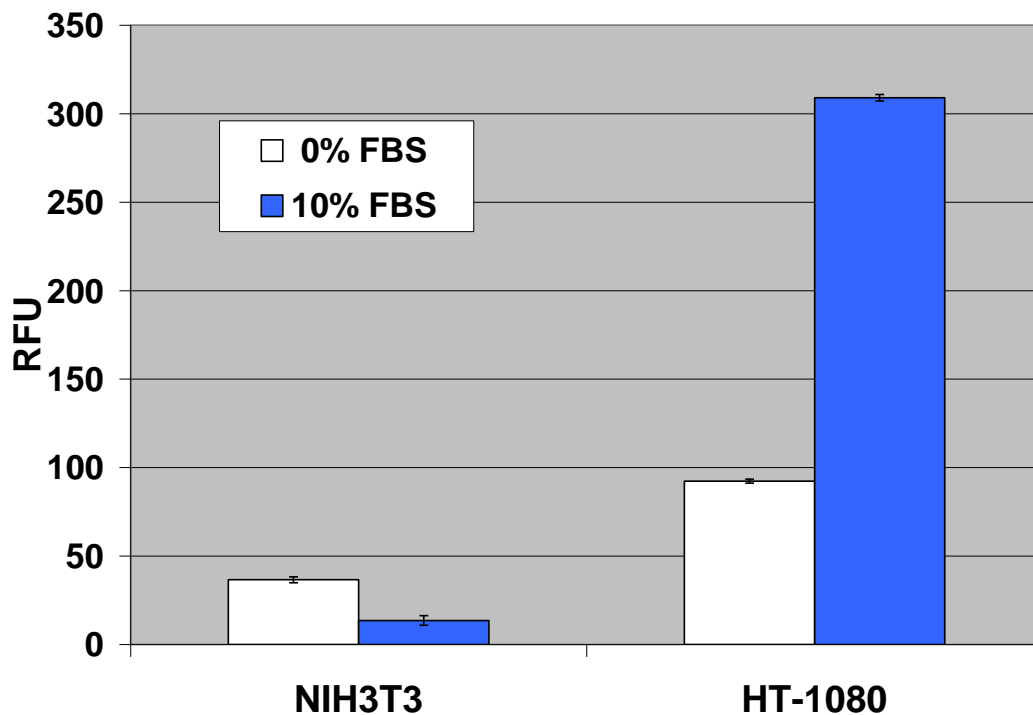
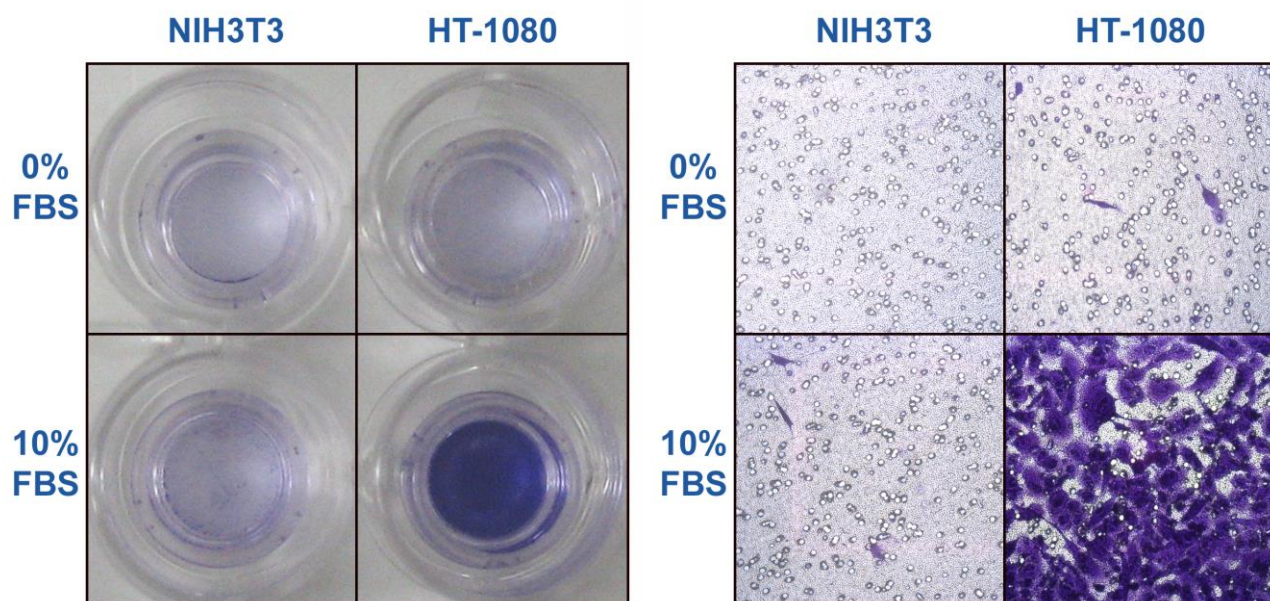
4. Carefully remove 250  $\mu\text{L}$  of rehydration medium (step 2) from the inserts without disturbing the collagen layer (leaving 50  $\mu\text{L}$  inside).
5. Add 250  $\mu\text{L}$  of the cell suspension solution to the inside of each insert.
6. Add 500  $\mu\text{L}$  of media containing 10% fetal bovine serum or desired chemoattractant(s) to the lower well of the invasion plate.
7. Incubate for 12-24 hours in a cell culture incubator.
8. Carefully aspirate the media from the inside of the insert. Transfer the insert to a clean well containing 225  $\mu\text{L}$  of Cell Detachment Solution. Incubate 30 minutes at 37°C.
9. Completely dislodge the invaded cells from the underside of the membrane by gently tilting the insert several times in the detachment solution. Remove and discard the insert.
10. 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  $\mu\text{L}$  to 370  $\mu\text{L}$  of 4X Lysis Buffer).
11. Add 75  $\mu\text{L}$  of 4X Lysis Buffer/CyQuant® GR dye solution to each well containing cells and 225  $\mu\text{L}$  of Cell Detachment Solution. Incubate 20 minutes at room temperature.
12. Transfer 200  $\mu\text{L}$  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™ Collagen Cell Invasion 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 cell suspension was titrated 1:2 in Cell Detachment Buffer. 150  $\mu$ L of each dilution was transferred to a 96-well, fluorometer plate, followed by addition of 50  $\mu$ L 4X Lysis Buffer/CyQuant® GR dye solution (1:75).



**Figure 2. Human Fibrosarcoma HT-1080 Collagen Cell Invasion.** HT-1080 and NIH3T3 (negative control) were seeded at 200,000 cells/well and allowed to invade toward FBS for 24 hrs. Invasive cells on the bottom of the invasion membrane were stained (top pictures) and quantified by CyQuant® GR dye as described in the Assay Protocol (bottom figure).

## **References**

1. Erkell, L. J., Schirmacher, V. (1988) *Cancer Res* **48**, 6933-6937.
2. Montgomery, A. M. P., De Clerck, Y. A., Langley, K. E., Reisfeld, R. A., Mueller, B. M. (1993) *Cancer Res* **53**,693-700.
3. 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. Zhen, Q. et al. (2020). EPAS1 promotes peritoneal carcinomatosis of non-small-cell lung cancer by enhancing mesothelial-mesenchymal transition. *Strahlenther Onkol*. doi: 10.1007/s00066-020-01665-1.
2. Negro, G. et al. (2020). Molecular heterogeneity in breast carcinoma cells with increased invasive capacities. *Radiol Oncol*. **54**(1):103-118. doi: 10.2478/raon-2020-0007.

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

©2006-2020: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.