

# Cell Invasion Assay Kit

## (Collagen I, 8 µm, 96-well)

### Instruction Manual

<b>Catalog Number</b>	PK-CA577-K916																							
<b>Description</b>	<p>Cell invasion is the ability of cells to migrate from one area to another through an extracellular matrix. Cell invasion is exhibited by both normal cells as well as cancerous cells in response to specific external signals, including chemical &amp; mechanical stimuli. During invasion, extracellular matrix is enzymatically degraded by cellular proteases before cells migrate to the new location. Cell invasion is required for normal processes such as wound repair, vasculature formation and the inflammatory response as well as the abnormal invasion of tissues by tumor cells during metastasis.</p> <p>PromoCell's <i>Cell Invasion Assay Kit</i> (Collagen I, 8 µm, 96-well) utilizes a Boyden chamber coated with Collagen I, where the cells invade the matrix and then migrate through a semipermeable membrane in the Boyden chamber in response to stimulants or inhibitory compounds. The percent cell invasion can be analyzed directly in a plate reader. Our assay is easy to use, sensitive and adaptable to high-throughput systems.</p>																							
<b>Quantity</b>	96 assays																							
<b>Applications</b>	<ul style="list-style-type: none"> <li>• Measure cell invasion in response to stimuli</li> <li>• Screen and characterize compounds that influence cell invasion</li> </ul>																							
<b>Sample Type</b>	<ul style="list-style-type: none"> <li>• Invasive cell lines</li> <li>• Invasion inhibitor or stimuli</li> </ul>																							
<b>Kit Components</b>	<table border="1"> <thead> <tr> <th>Components</th> <th>Quantity</th> <th>Cap code</th> </tr> </thead> <tbody> <tr> <td>Cell Invasion Chamber</td> <td>1 each</td> <td>Plate</td> </tr> <tr> <td>Collagen I</td> <td>5 ml</td> <td>Clear</td> </tr> <tr> <td>Control Invasion Inducer</td> <td>300 µl</td> <td>Red</td> </tr> <tr> <td>Cell Dissociation Solution</td> <td>10 ml</td> <td>NM</td> </tr> <tr> <td>Wash Buffer</td> <td>50 ml</td> <td>NM</td> </tr> <tr> <td>Cell Dye</td> <td>1 vial</td> <td>Green</td> </tr> </tbody> </table>	Components	Quantity	Cap code	Cell Invasion Chamber	1 each	Plate	Collagen I	5 ml	Clear	Control Invasion Inducer	300 µl	Red	Cell Dissociation Solution	10 ml	NM	Wash Buffer	50 ml	NM	Cell Dye	1 vial	Green		
Components	Quantity	Cap code																						
Cell Invasion Chamber	1 each	Plate																						
Collagen I	5 ml	Clear																						
Control Invasion Inducer	300 µl	Red																						
Cell Dissociation Solution	10 ml	NM																						
Wash Buffer	50 ml	NM																						
Cell Dye	1 vial	Green																						
<b>User Supplied Reagents &amp; Equipment</b>	<ul style="list-style-type: none"> <li>• Fluorescence Plate Reader</li> <li>• Cell Culture Media</li> <li>• Cotton Swabs</li> <li>• Centrifuge to spin 96-well plate</li> <li>• 96-well clear bottom white plate</li> </ul> <p>For human primary cells and the corresponding optimized cell culture media visit <a href="http://www.promocell.com">www.promocell.com</a>. For Cytokines/Chemokines and Growth Factors visit <a href="http://www.promokine.info">www.promokine.info</a>.</p>																							
<b>Applications / Assay Protocol</b>	<p><b>A. General Consideration and Reagent Preparation:</b></p> <p>Store kit at -20°C, protected from light. Briefly centrifuge small vials prior to opening. Assay is performed under sterile conditions. Read entire protocol before performing the experiment.</p> <p><b>Cell Invasion Chamber:</b> Open under sterile conditions. Keep at room temperature.</p> <p><b>Collagen I:</b> Aliquot under hood and store at 4°C, if needed.</p> <p><b>Control Invasion Inducer, Cell Dissociation Solution and Wash Buffer:</b> Store at -20°C. Bring to 37°C before use. Stable for six months.</p> <p><b>Cell Dye:</b> Add 100 µl of DMSO (Cat. No. PK-CA707-90082, not provided) to the vial to dissolve the dye. Aliquot and store at -20°C, protected from light.</p> <p><b>B. Cell Invasion Assay Protocol:</b></p> <ol style="list-style-type: none"> <li>1. Add 50 µl Collagen I to coat desired wells of the Top Chamber. Incubate plate at room temperature for 2-3 hours in flow hood or overnight 2-8°C to form a thin film of Collagen I. Check the chamber from the side to make sure the plates are dried. Incubate for a longer duration if needed. Wash the coated plate three times with 100 µl 1 M Tris (not provided). Aspirate.</li> <li>2. Grow cells of interest in desired media and culture conditions. Grow enough cells to perform a Cell</li> </ol>																							

Invasion Assay and a Standard Curve. Adherent cells should be cultured to ~80% confluence.

**3.** Prior to the assay, starve cells for 18-24 hours in a serum-free media (0.5% serum can be used, if needed). After starvation, harvest the cells and centrifuge at 1,000 x g, for 5 minutes to pellet cells. Resuspend cell pellet in Wash Buffer and count the number of cells using hemocytometer or automated cell counter. Resuspend cells at  $1 \times 10^6$  cells/ml in a serum-free media.

**4.** Under sterile conditions, disassemble the Cell Invasion Chamber and carefully remove the plate cover and the top chamber.

**Bottom Chamber:** Add 200  $\mu$ l of medium per well containing the desired chemoattractant to the bottom chamber. In control well(s), we recommend omitting the chemoattractant. For Positive Control, add 20  $\mu$ l of Control Invasion Inducer to 180  $\mu$ l of medium in the bottom chamber. Reassemble the top and bottom chambers while ensuring no air bubbles are trapped between them.

**Top Chamber:** Add 50  $\mu$ l (~50,000 cells) of cell suspension to each well of the top chamber. Add desired stimulator or inhibitor to the top well, and gently mix. Make up the volume to 100  $\mu$ l with media. Carefully replace the plate cover and incubate the Cell Invasion Chamber at 37°C in CO<sub>2</sub> incubator for 2-48 hours.

**Note:** Invasive cells pass through the Collagen I and cling to the outer side of the top chamber. Non-invasive cells stay in the upper chamber.

#### 5. Standard Curve:

**A.** Each cell type requires a separate Standard Curve. Prepare a Standard Curve by adding 50  $\mu$ l cells ( $1 \times 10^6$  cells/ml, ~50,000 cells) in desired well(s) in a 96-well plate (white plate clear bottom). Serially dilute the cells 1:1 in Wash Buffer and generate a Standard Curve with different cell numbers (50,000; 25,000; 12,500; 6,250; 3,125; 1,562; 781 and 390) in 100  $\mu$ l total volume. As blank, use 100  $\mu$ l of Wash Buffer.

**B.** Dilute Cell Dye 1:250 in PBS and add 50  $\mu$ l of diluted Cell Dye to each well. Incubate at 37°C for 30 minutes. Read the fluorescence at Ex/Em = 485/530 nm. Plot the Standard Curve (Number of Cells Vs RFU obtained). Fit the data points using a linear trend line with zero intercept. The equation for the straight line and R-squared value are used for data analysis of samples.

**Note:** The Cell Invasion RFU reading should fall in the linear range of the Standard Curve. We recommend using triplicates for Standard Curve.

#### 6. Separation of Invasive Cells:

**A.** After the desired incubation with cell invasion inducers/inhibitors, carefully remove the plate cover and aspirate media from the top chamber without puncturing the membrane and matrix.

**B.** Remove cells from the top chamber using a cotton swab. Disassemble the Cell Invasion Chamber by removing the top chamber. Invert the top chamber and set it aside. Place the plate cover on top of bottom chamber and centrifuge the plate at 1,000 x g for 5 minutes at room temperature.

**C.** Carefully aspirate the media from the bottom chamber and wash the chamber with 200  $\mu$ l Wash Buffer. Centrifuge the plate at 1,000 x g for 5 minutes at room temperature and aspirate the Wash Buffer from the bottom chamber.

#### 7. Count Invasive Cells:

**A.** Dilute Cell Dye 1:500 in Cell Dissociation Solution. Mix well. Make desired amount of Cell Dye solution depending on the number of wells.

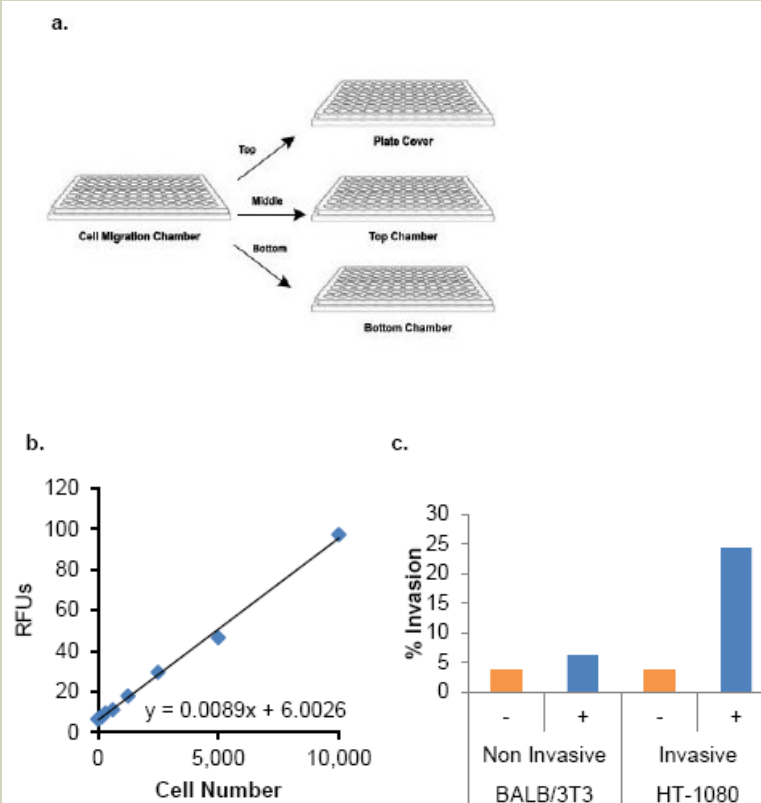
**B.** Add 100  $\mu$ l of the mix to each well of the bottom chamber. Reassemble the Cell Invasion Chamber by placing the top chamber into the bottom chamber. Incubate at 37°C in CO<sub>2</sub> incubator for 30 minutes\*. After incubation, disassemble the Cell Invasion Chamber and remove the top chamber.

**C.** Read the plate at Ex/Em = 485/530 nm. Calculate the number of cells invaded using the equation of the straight line obtained from Standard Curve. Percentage Invasion can be calculated as follows:

#### Calculation:

$$\% \text{ Invasion} = \left( \frac{\text{Number of Cells in Lower Chamber}}{\text{Total Number of Cells added to Top Chamber}} \right) \times 100$$

**\*Note:** During incubation with Cell Dissociation Solution/Cell Invasion Dye, gently tap the plate on the side to ensure optimal dissociation of the invasive cells that cling to the outer side of the top chamber.



**Figure 1:** (a) Cell Invasion plate: The cells are added to the Top Chamber and the Control Invasion Inducer or chemoattractant are added to the Bottom Chamber. (b) Standard Curve: HT-1080 cells were harvested, counted and serially diluted to obtain desired cell number. Cells were incubated according to the protocol. (c) Cell Invasion: BALB/3T3 and HT-1080 cells were starved overnight and treated with Control (Cnt) Invasion Inducer or remain untreated (No Treatment). Treatment with Control Invasion Inducer demonstrated a significant increase in invasion of HT 1080 cells as compared to BALB/3T3 control cells.

#### Intended Use

For in vitro research use only. Not for diagnostic or therapeutic procedures.

#### Storage & Stability

Store kit at -20°C, protected from light. Briefly centrifuge small vials prior to opening.

#### PromoCell GmbH

Sickingenstr. 63/65  
69126 Heidelberg  
Germany

Email: [info@promokine.info](mailto:info@promokine.info)  
[www.promokine.info](http://www.promokine.info)

#### USA/Canada

Phone: 1 – 866 – 251 – 2860 (toll free)  
Fax: 1 – 866 – 827 – 9219 (toll free)

#### Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)  
Fax: 0800 – 100 83 06 (gebührenfrei)

#### France

Téléphone: 0800 90 93 32 (ligne verte)  
Téléfax: 0800 90 27 36 (ligne verte)

#### United Kingdom

Phone: 0800 – 96 03 33 (toll free)  
Fax: 0800 – 169 85 54 (toll free)

#### Other Countries

Phone: +49 6221 – 649 34 0  
Fax: +49 6221 – 649 34 40