

Cell Migration Assay, TC

Catalog #: ENZ-KIT113

ENZ-KIT113-0001

1 x 96 wells

96-well, 2-D Assay for Investigating Cell Migration of Adherent Cell Lines

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Introduction	2
Plate Dimensions	3
Materials Provided	4
Materials Required	4
Cell Migration Assay Protocol	5
Data Acquisition	9
Appendix I: Determining Optimal Cell Seeding Concentration	11
Appendix II: Determining Optimal Fluorescence Plate Reader Settings	12
Contact Information.....	14

INTRODUCTION

The Cell Migration Assay is a reproducible, sensitive, and flexible assay that can be used to monitor cell migration. Formatted for a 96-well plate, the assay utilizes Cell Seeding Stoppers made from a medical-grade silicone to restrict cell seeding to the outer annular regions of the wells. Removal of the stoppers reveals a 2mm diameter unseeded region in the center of each well, i.e., the detection zone, into which the seeded cells may then migrate. The Detection Mask is applied to the plate bottom and restricts visualization to the detection zones, allowing only cells that have migrated to be detected (see **Figure 1**). The Cell Migration Assay is designed to be used with any commercially available stain or labeling technique. Readout can be performed by microscopy or use of a microplate reader.

The Cell Migration Assay system has been designed for use with adherent cell cultures. This assay has been successfully used with HT-1080, PC-3, A549, NCI H1650, MDA-MB-231, NMuMG, 3T3-Swiss albino, HCEC, HUVEC, and MCF10A cell lines.

Using the Cell Migration Assay offers the following benefits:

- **Membrane-free Migration:** perform studies without manipulating transmembrane inserts.
- **Reproducible Results:** obtain well-to-well CV's <12% due to the unique design.
- **Preserves Cell Morphology:** monitor changes in cell structure in real-time.
- **Versatile:** analyze data using multiple probes in a single well by using a microscope, digital imager, or fluorescence microplate reader.
- **Flexible:** perform kinetic or endpoint Cell Migration Assays without the use of special instrumentation.



Fig 1. Schematic of Cell Migration Assay

PLATE DIMENSIONS

Diameter of Well	6.5mm
Diameter of Stopper Space (Detection Zone)	2mm
Suggested Media Volume per Well (populated with Stoppers)	100µl
Effective Area of Outer Annular Region (seeding region) per Well	30.03mm ²
Effective Area of Central Detection Zone per Well	3.14mm ²
Plate Height	14.9mm
Plate Height with Lid (with Cell Seeding Stoppers)	17.9mm
Offset of Wells (A-1 location, X)	14.4mm
Offset of Wells (A-1 location, Y)	11.2mm
Distance between Wells	9mm (on center)
Well Depth	12.2mm
Thickness of Well Bottom	0.25mm
Storage Conditions	Refrigerate (4°C)



Important: Read instructions before performing any Assay.

MATERIALS PROVIDED

- 96-well, Tissue Culture Treated (black, clear bottom) Plate with Cell Seeding Stoppers, 1
- Detection Mask, 1
- Stopper Tool, 1

MATERIALS REQUIRED

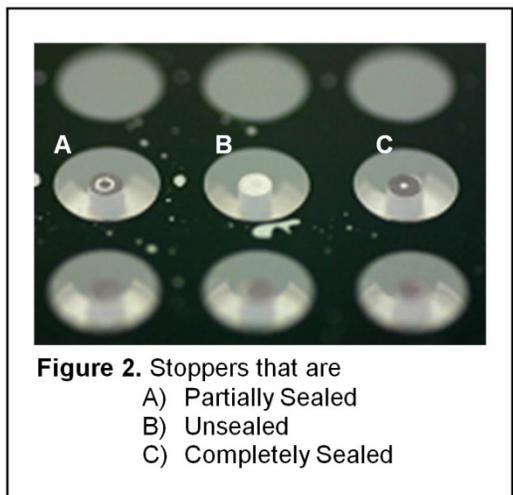
- Biological Cells
- Sterile PBS (containing both Ca⁺⁺ and Mg⁺⁺)
- Complete Cell Culture Growth Medium (containing serum)
- Sterile Pipette Tips/Pipette or Multi-Channel Pipette
- Trypsin or Cell Scraper
- Inverted Microscope (optional)
- Fluorescence Microplate Reader (optional)
- Cell Culture Labeling Medium (phenol red-free/serum-free media)
- Cell Labeling Fluorescent Agent (eg., Cyto-ID[®] Green Cell Tracer Kit, ENZ-51036 or Cyto-ID[®] Red Cell Tracer Kit, ENZ-51037) *(required if performing assay readout via fluorescence analysis)*

CELL MIGRATION ASSAY PROTOCOL

The following steps should be performed in a biological hood using aseptic technique to prevent contamination.

1. Remove the Cell Migration Assay Plate from refrigeration and place on lab bench for ~1 hour to allow it to equilibrate to room temperature.

2. Visually inspect the underside of the populated 96-well plate to ensure that the Cell Seeding Stoppers are firmly sealed against the bottom of the plate. To inspect the stoppers, turn the plate over and examine the stoppers for sealing (see **Figure 2**). If incomplete sealing is observed, return the plate to the upright position and use a sterile instrument to gently push the stopper back into the well until sealing is observed.

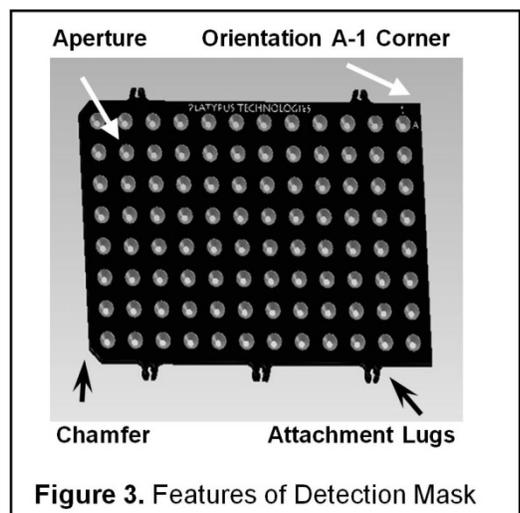


NOTE: the sealing of the stoppers can be most easily observed if the plate is tipped at an angle and viewed under indirect light to reveal the “bullseye” pattern at the bottom of each well.

3. Apply the Detection Mask to the bottom of the 96-well plate if microplate reader data is being collected. The Detection Mask is not necessary if collecting imaging data.

First Time Users: In order to prevent splashing of well contents, familiarize yourself with the attachment and removal of the Detection Mask before any liquids are placed in the wells.

- Orient the chamfered corners of the mask with those of the 96-well plate, ensuring that the A1 corner of the mask is aligned with the A1 well of the plate (see **Figure 3**).



- Align the holes in the attachment lugs with the bosses on the bottom of the 96-well plate.
- Gently press the mask until it is flush with the bottom of the 96-well plate.

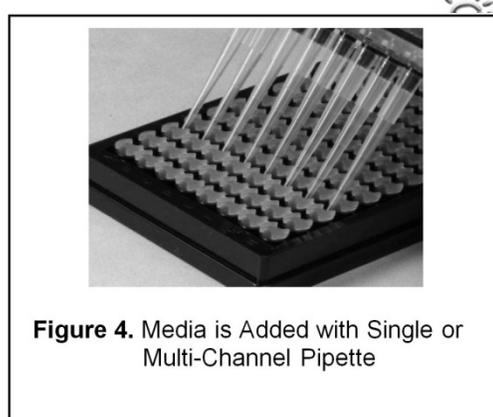


NOTE: It may be necessary to wash the mask with ethanol to remove dust and debris since the mask is not sterile. The mask may be applied at any point during the assay. For kinetic assays, it is often most convenient to apply the mask at the beginning of the assay before any liquids are placed in the well. For endpoint assays, using fixed and stained cells, it is often most convenient to apply the mask just before reading assay results.

4. If performing a kinetic analysis of cell migration, pre-label with a fluorescent stain now.
5. Collect cells and prepare a suspension that is 10-fold greater in density than the optimal seeding concentration.

First Time Users: The optimum seeding density of cells must be determined as an integral part of the design of the Cell Migration Assay. Please refer to **Appendix I** for a discussion of this process.

6. Pipette 100µl of suspended cells into each test well through one of the side ports of the Cell Seeding Stopper.



NOTE: For best results, add or extract media by placing the pipette tip along the wall of the well (see **Figure 4**). Care should be taken not to disturb the Cell Seeding Stopper when introducing the pipette tip into the well. A slender/elongated tip or a gel loading tip may be useful.

7. **IMPORTANT:** Lightly tap the plate on your work surface to evenly distribute well contents (extreme tapping may result in splashing of well contents and lead to contamination).

8. Incubate the seeded plate containing the Cell Seeding Stoppers in a humidified chamber (37°C, 5% CO₂) for 4 to 18 hours (cell line dependent) to permit cell attachment.
9. Remove plate from incubator.
10. Designate several 'reference' wells in which the stoppers will remain in place until results are read (t=0 pre-migration controls).
11. Using the Stopper Tool, remove all other stoppers (see **Figure 5**).



NOTE: It may be necessary to wash the Stopper Tool with 70% ethanol as the Stopper Tool is not sterile.

- Secure the 96-well plate by holding it firmly against the deck of your work space. Slide the tines of the Stopper Tool under the backbone of the stopper strip, keeping the underside of the Stopper Tool flush with the top surface of the plate.

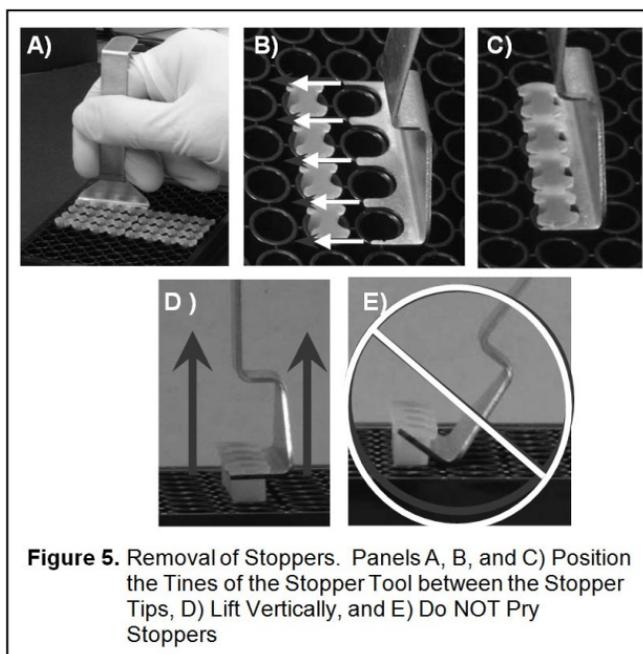


Figure 5. Removal of Stoppers. Panels A, B, and C) Position the Tines of the Stopper Tool between the Stopper Tips, D) Lift Vertically, and E) Do NOT Pry Stoppers

- Lift the Stopper Tool vertically to gently remove the stoppers.



NOTE: DO NOT use the Stopper Tool as a lever to pry the stoppers from the well (see **Figure 5E**), as doing so may cause displacement of seeded cells and may distort the detection zone area.

12. Remove media with a pipette and gently wash wells with 100µl of sterile PBS (or media) to remove any unattached cells. Do not aspirate using an in-house vacuum.
13. Add 100µl of fresh culture media to each well.

14. Incubate plate in a humidified chamber (37°C, 5% CO₂) to permit cell migration. Cells may be examined microscopically throughout the incubation period to monitor progression of migration. Migration time will vary depending upon cell type and experimental design.
15. If performing an endpoint analysis of cell migration, stain cells with a fluorescent stain after sufficient migration has occurred. Refer to next section and **Appendix II** for further information on data acquisition and fluorescence staining technique.



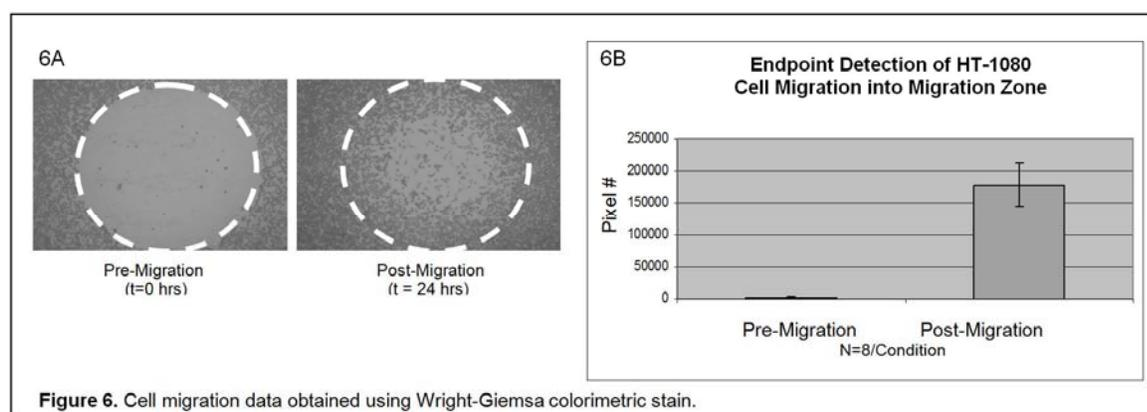
NOTE: Cell Seeding Stoppers are for single use only; Enzo cannot guarantee the integrity of the stopper material after a second sterilization procedure.

DATA ACQUISITION

The readout of the Cell Migration Assay can be conducted at any time, allowing the user to perform a kinetic assay or an endpoint assay. The Cell Migration Assay is designed to be used with any commercially available stain or labeling technique. The readout can be performed by using a microscope, a microplate reader, or a High Content Screening or High Content Imaging Analysis platform.

Microscope Analysis

- Cell counting or image capture / analysis software, such as NIH ImageJ freeware, can be used.
- **Note:** Microscopy observations are possible using phase contrast or bright field microscopy.
- No need to attach the Detection Mask to the plate.
- Sample data using a colorimetric stain is shown in **Figure 6**. Wells populated with Cell Seeding Stoppers were seeded with 50,000 HT-1080 cells (i.e., 100 μ l of 5×10^5 cells/mL) and incubated overnight. The stoppers were removed from test wells, but remained in place in the pre-migration reference wells until the time of the assay readout. The seeded plate was incubated in a humidified chamber for 24 hours to permit cell migration. Stoppers were removed from the reference wells and all cells were fixed and treated with Wright-Giemsa stain. Images were captured using bright field microscopy and then imported to Image J software for analysis using thresholding. The images below, captured without a detection mask in place, illustrate representative data from pre-migration (t=0 hrs) and post-migration (t=24 hrs) wells (**Figure 6A**). The graph depicts the average pixel number in the detection zones for each condition (**Figure 6B**).



Microplate Reader Analysis

- Attach the Detection Mask to the bottom of the plate (refer to **Protocol** section, step 3).
- Optimal settings will vary according to the microplate reader make and model. Consult **Appendix II** and the equipment user manual for your particular instrument.
- The microplate reader **MUST** be set to read from the bottom of the plate.
- Sample data using a fluorescent stain and microplate reader analysis are shown in **Figure 7**. Wells populated with Cell Seeding Stoppers were seeded with 50,000 HT-1080 cells (i.e., 100 μ l of 5x10⁵ cells/mL) and incubated overnight. The stoppers were removed from test wells, but remained in place in the pre-migration reference wells until the time of the assay readout. Cells were fluorescently stained with a green cell tracing dye. The seeded plate was incubated in a humidified chamber for 28 hours and at various time points the fluorescence signals in the detection zones were measured using a plate reader. The images below (**Figure 7A**), captured without a detection mask in place, illustrate representative data from pre-migration (t=0 hrs) and post-migration (t = 21 hrs) wells. The graph depicts a real-time analysis of cell migration that was prepared by transposing the fluorescent signal into cell numbers (**Figure 7B**).

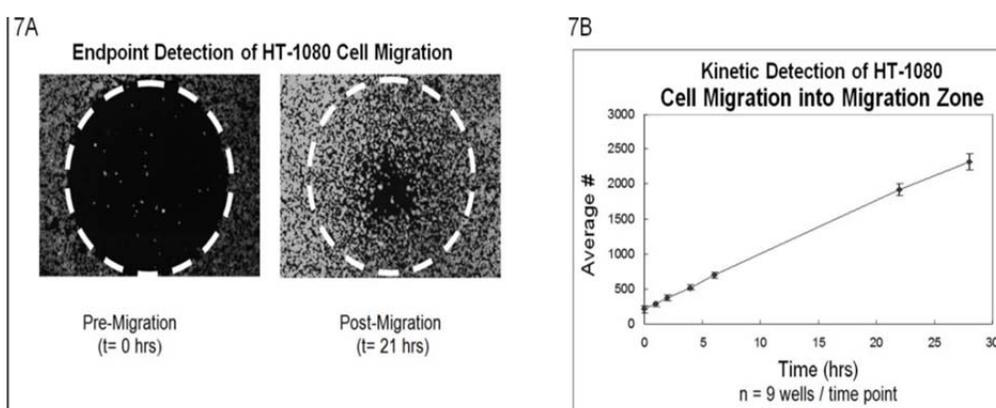


Figure 7. Cell migration data obtained using green fluorescent cell tracing dye.

APPENDIX I: DETERMINING OPTIMAL CELL SEEDING CONCENTRATION

This procedure is intended to assist in determining the cell seeding density needed to achieve confluency of your cell line when using the Cell Migration Assay. The intended goal is to achieve 90-95% confluency of the monolayer surrounding the Cell Seeding Stoppers without overgrowth.

1. A suggested starting point is to evaluate three serial dilutions at the cell densities shown below. The cell seeding area of the well with the stopper in place is ~ 0.3 cm². Based on the typical seeding density of your particular cell line, you can infer a different cell number for your first serial dilution and adjust the numbers below accordingly.
2. Prepare a log-phase culture of the cell line to be tested. Collect cells and determine the total number of cells present.
3. Pellet cells by centrifugation. Prepare three serial dilutions at final concentrations of 1.0 x 10⁶, 0.5 x 10⁶ and 0.25 x 10⁶ cells/mL.
4. Dispense 100µl of cell suspension per well into the 96-well plate to result in the following plate layout:

Column	1	2	3
Cells / well	100,000	50,000	25,000
Number of wells	8	8	8

5. Incubate the plate in a humidified chamber (37°C, 5% CO₂) for 4-18 hours (cell line dependent) with cell seeding stoppers in place to allow the cells to firmly attach to the well surface.
6. Following cell attachment, remove the Cell Seeding Stoppers from each well (see **Figure 5**) and gently wash the wells with PBS to remove non-attached cells.
 - Secure the 96-well plate by holding it firmly against the deck of your work space. Slide the tines of the Stopper Tool under the backbone of the stopper strip, keeping the underside of the tool flush with the top surface of the plate.

- Lift the Stopper Tool vertically to gently remove the stopper. Do not use the Stopper Tool as a lever to pry the stoppers from the well as doing so may cause displacement of the seeded cells.
7. Without a Detection Mask in place, use a microscope to visually inspect each well to determine the minimum cell seeding concentration that yields a confluent monolayer at the perimeter of the detection zone.

At this point, if you plan to obtain the results of the Cell Migration Assay via colorimetric or microscopy analysis, you have successfully determined the optimal cell seeding concentration to be used in Step 5 of the **Cell Migration Assay Protocol**.

APPENDIX II: DETERMINING OPTIMAL FLUORESCENCE MICROPLATE READER SETTINGS

This procedure is intended to assist in optimizing your instrument settings when using a fluorescence microplate reader to capture data from the Cell Migration Stopper Assay.

1. Using the optimal cell seeding concentration determined in **Appendix I**, perform a Cell Migration Assay per the **Cell Migration Assay Protocol** using culture conditions expected to result in robust cell migration. Be sure to include equal numbers of pre-migration reference wells (stoppers left in place until staining) and post-migration test wells (stoppers removed after cell attachment period). A minimum of 8 wells per condition are recommended.
2. Perform the desired fluorescent staining technique.

The Cell Migration Assay has been designed to work with all types of fluorescent stains and staining techniques. The precise method for staining cells with fluorescent stains varies according to the nature of the individual stain. It is important to stain cells using a fluorescent reagent that uniformly stains cells. Probes affected by experimental conditions will increase variability of results and reduce correlation between fluorescence signal and cell migration. Please consult the manufacturer of your fluorescent stain for specific considerations.

The following is an example Fluorescent Staining Protocol for using Calcein AM:

- a. To stain one fully-seeded 96-well plate, combine 5 μ l of Calcein AM (1mg/mL in dry DMSO) with 10ml of phenol red-free and serum-free media or 1x PBS (containing both Ca⁺⁺ and Mg⁺⁺). Protect diluted Calcein AM solution from light until ready to use in step d.
 - b. Carefully remove culture medium from wells.
 - c. Wash wells with 100 μ l of PBS (containing both Ca⁺⁺ and Mg⁺⁺).
 - d. Add 100 μ l of diluted Calcein AM solution to each well.
 - e. Incubate plate at 37°C for 30-60 minutes.
 - f. Attach mask and read promptly with microplate reader using appropriate filter set and sensitivity/gain settings (for a BioTek Synergy™ HT microplate reader, use 485/528nm excitation/emission filters, sensitivity 55nm).
3. If not already in place, apply the Detection Mask to the plate. Using the bottom probe of a fluorescence microplate reader, obtain the fluorescence reading from each well. To achieve the optimal dynamic range, adjust the instrument settings (e.g., gain) to result in the greatest difference in fluorescence signal between pre-migration and post-migration wells. Refer to the instrument manual for your microplate reader for further guidance on instrument settings.

You have now successfully determined the optimal cell seeding concentration (to be used in Step 5 of the **Cell Migration Assay Protocol**) and microplate reader settings for analysis of cell migration using a fluorescence microplate reader.



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com