

Product datasheet

Cell Migration/Chemotaxis Assay Kit (96-well, 8 μm) ab235673

★★★★☆ [1 Abreviews](#) [1 References](#) [2 Images](#)

Overview

Product name	Cell Migration/Chemotaxis Assay Kit (96-well, 8 μm)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Product overview	Cell Migration/Chemotaxis Assay Kit (96-well, 8 μm) (ab235673) measures cell migration in response to stimuli in adherent and suspension cells. It allows you to screen, study, or characterize compounds that influence chemotaxis/cell migration. It utilizes a Boyden chamber, where the cells migrate through a semi-permeable membrane under different stimuli. Cell migration can be analyzed directly by reading fluorescence (Ex/Em = 530/590 nm) in a plate reader. Our assay is easy to use, sensitive and adaptable to high-throughput systems.

8 μm semi-permeable membrane maybe too large to use with tiny cell types, such as the cell types which diameter less than 8 μm .

Notes This product is manufactured by BioVision, an Abcam company and was previously called K906 EZCell™ Cell Migration/Chemotaxis Assay Kit (96-well, 8 μm). K906-100 is the same size as the 100 test size of ab235673.

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 and 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.

Platform Microplate reader

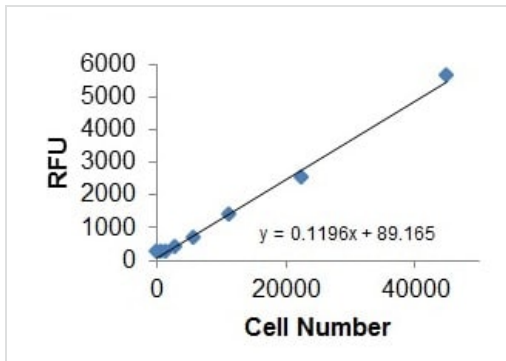
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests	100 tests
Cell Chamber (96 x 8 μm)	1 unit	1 unit

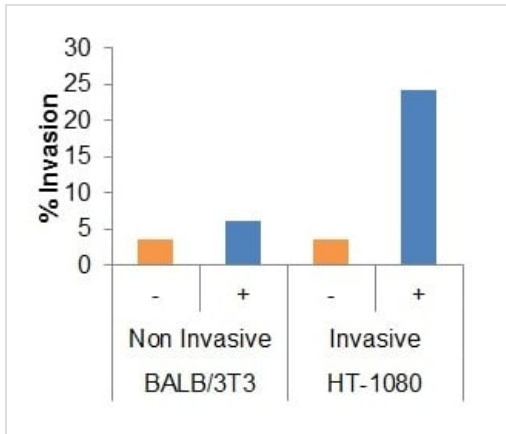
Components	100 tests	100 tests
Cell Dissociation Solution I	1 x 15ml	1 x 10ml
Cell Dye I	1 x 1.5ml	1 x 1.5ml
Control Migration Inducer	1 x 300µl	1 x 300µl
Wash Buffer II	1 x 50ml	1 x 50ml

Images



Standard Curve.

HT-1080 cells were harvested, counted and serially diluted to obtain desired cell number. Cells were incubated according to the protocol.



Cell Invasion.

3T3-NIH 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 compare to 3T3-NIH control cells.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors