

Cell Invasion Assay (Collagen IV), 96-well, 8 µm ab235888

2 Images

Overview

Product name	Cell Invasion Assay (Collagen IV), 96-well, 8 µm
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Species reactivity	Reacts with: Mouse, Human
Product overview	Cell Invasion Assay (Collagen IV), 96-well, 8 µm (ab235888) utilizes a Boyden chamber coated with Collagen IV, where the cells invade the matrix and then migrate through a semi-permeable 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.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K918 EZCell™ Cell Invasion Assay (Collagen IV), 96-well, 8 µm. K918-100 is the same size as the 100 test size of ab235888.</p> <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 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.</p>
Platform	Microplate reader

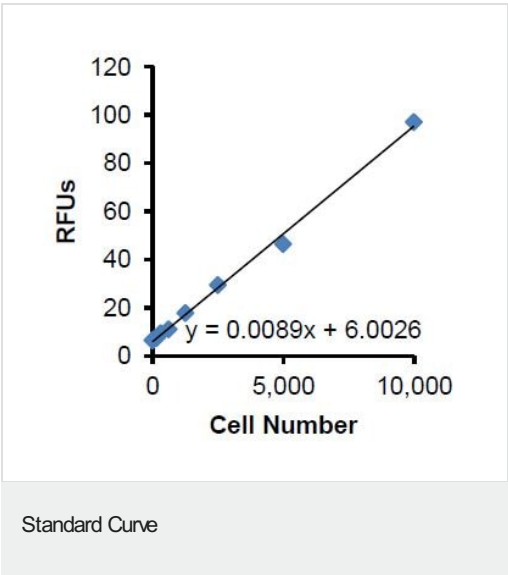
Properties

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

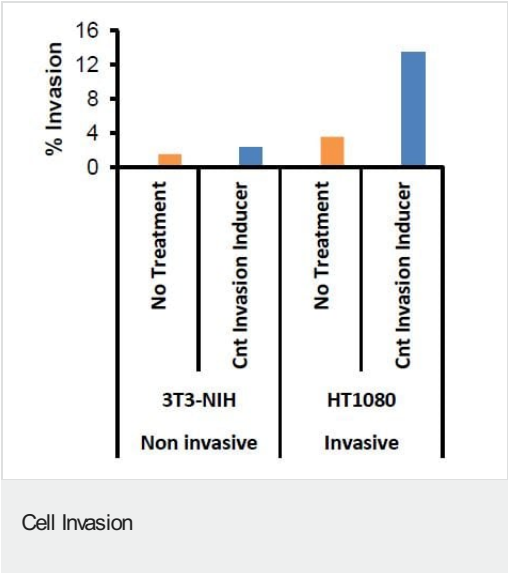
Components	100 tests
Cell Dissociation Solution	1 x 15ml
Cell Dye	1 vial

Components	100 tests
Cell Invasion Chamber	1 unit
Collagen IV	1 x 5ml
Control Invasion Inducer	1 x 300µl
Wash Buffer	1 x 50ml

Images



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



NIH-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 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