

# CytoSelect™ 96-well Cell Migration Assay (5 µm), Fluorometric

Chemotaxis describes the movement of cells toward or away from a chemical stimulus in their environment. Cell chemotaxis plays a pivotal role in the progression of cancer and other diseases. CytoSelect™ Cell Migration Assays are ideal for determining the chemotactic properties of cells. The 5 µm pore size is ideal for monocytes / macrophages.

## Ordering Information

[Order Online »](#)

CBI-CBA-105	96 assays
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

- Fully quantify chemotaxis with no manual cell counting
- Measure chemotaxis in less than 6 hours with most cell types
- Membrane inserts are uncoated to allow use with any chemoattractant

## Handling & Storage

Long Term Storage      +4°C

Shipping                      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application                      Fluorescent detection

**Technical Info / Product Notes**      For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)