

Trevigen Cell Assays, (TCA) a division of Trevigen, Inc, was established in 2008 to conduct contract research work for the pharmaceutical, biotechnology, government and academic segments of the research market. TCA specializes in designing and conducting assays for lead compound and genotoxic screening based on DNA damage and repair and cancer cell behavior.

TREVIGEN®**CELL INVASION/MIGRATION ASSAY**

Cultrex® Cell Invasion and Migration Assays

Cultrex® Cell Invasion Assays

The Cultrex® Cell Invasion Assays were created in an effort to accelerate the screening process for compounds that influence cellular digestion and migration across extracellular matrices, which is a fundamental component of cellular processes such as angiogenesis, embryonic development, immune responses, and tumor cell metastasis. These assays offer a flexible, standardized, high-throughput format for quantitating the degree to which invasive cells penetrate a barrier consisting of basement membrane components in vitro in response to chemoattractants and/or inhibiting compounds.

These assays employ a simplified Boyden chamber design with a polyethylene terephthalate (PET) membrane containing 8 micron pores, which allow access from the input chamber (top) to the assay chamber (bottom) without dismantling the device. The assay chamber may be directly analyzed in a 96-well plate reader, eliminating transfer steps that introduce additional variability to the assay.

Cell invasion is quantified using calcein-acetoxymethyl ester (Calcein-AM), which is internalized by the cells, and cleaved by intracellular esterases to generate fluorescence signals. In conjunction with a standard curve, these signals may be used to quantitate the number of cells that have migrated or invaded – thereby eliminating the need for direct cell counting. Since different cell lines and different treatments can result in a wide range of invasive potentials, the permissiveness of each matrix may also be optimized to fit each experiment by adjusting the coating concentration. Cultrex® Cell Invasion Assays are provided in multiple formats for evaluation against different extracellular matrices and matrix components: Laminin I, Collagen I, Collagen IV, and Basement Membrane Extract (BME).

Cultrex® Cell Migration Assays

The Cultrex® Cell Migration Assays utilize uncoated Boyden chambers to assess cell migration in the absence of extracellular proteins.

Assay	Coating	Catalog #
Invasion	BME	3455
Invasion	Laminin I	3456
Invasion	Collagen I	3457
Invasion	Collagen IV	3458
Migration	None	3465

Trevigen, Inc.
8405 Helgerman Court
Gaithersburg, Maryland 20877
info@trevigen.com
1-800-873-8443 • www.trevigen.com

Getting started is easy. Fax the completed form to 301-560-4973; or email us at TCA@trevigen.com; or complete the form online at www.trevigencellassays.com with the information that we need in order to set up your screening service.

Quotation Request Form - Cell Invasion/Migration Assay

First and Last Name

Email Address

Company

Street Address

City State/Province

Zip/Postal Code Country

Telephone Number Fax Number

What type of assay is required? **Cell Invasion – Cat. #** **Migration – Cat. #**

Please check the appropriate assay and write in the catalog #. If unsure of the proper kit to use, please contact us and we can aid you in identifying the proper kit for your needs.

How many samples do you have to screen?

How many replicates are required?

How many compounds and concentrations do you have to screen?

What type of cells are required?

Will you be providing the cells?

What is the desired reporting format?

How soon is the data required?

Are there any other screening parameters or special conditions that you require?

Compound Handling Instructions

What compound(s) are you screening?

Will you be providing the compound(s)?

If you are not providing the compound(s), where can they be purchased?

Is the compound toxic? If yes, are MSDS available?

What storage conditions are required?

What diluent is required?

Upon receipt, a TCA senior scientist will contact you to go over the desired work and discuss options as appropriate. A proposal and cost will then be prepared. The proposal will include the turn around time and the agreed upon reporting format.

CELL INVASION/MIGRATION ASSAY