

CULTREX[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Cultrex[®] Cell Staining Kit

Catalog #: 3437-100-K

Kit contents: 3437-100-01 CS Solution **Size:** 100 ml
3437-100-02 10X CS Wash Solution **Size:** 100 ml

Description: The Cell Staining Kit provides a ready-to-use staining solution, CS Solution (cat# 3437-100-01), and Wash Solution cat# 3437-100-02. The CS Solution contains a mixture of Azur A and Methylene Blue specially formulated to provide optimized staining of cells and structures grown on Basement Matrix Extract (BME). Ideal staining contrast is achieved with minimum background since excess stain can be easily washed away using the wash buffer. The blue coloration allows optimal contrast between cells and the growth matrix.

Storage Conditions: Store at room temperature.

Applications:

Staining of cells grown on Cultrex Basement Membrane Extract (BME) or other extra-cellular membrane protein extracts after application such as:

- Tube Assay: Basement Membrane Extract promotes differentiation of a mouse endothelial cell line derived from axillary lymph node (SVEC4-10), into capillary-like structures.
- Neurite outgrowth assay: Basement Membrane Extract promotes differentiation of neurite from chick dorsal root ganglion.
- Angiogenesis assay such as the Ring Assay: Basement Membrane Extract promotes differentiation of mouse aorta tissue to form capillary-like structures.

Staining Procedure:

Note: This procedure requires methanol and distilled or deionized water to dilute the 10X CS Wash Solution)

- 1- Remove culture media from well.
- 2- Add enough methanol to cover sample and incubate 1 minute at room temperature.
- 3- Remove methanol and wash 3 times with equal volumes of 1X CS Wash Solution.
- 4- Add enough CS Solution to cover samples in culture wells.
- 5- Leave stain on samples from 10 seconds to 1 minute.
- 6- Remove excess dye with pipette.
- 7- Wash samples 3 times with 1X CS Wash Solution or until desired destaining is achieved.
- 8- Discard excess and observe under microscope.
- 9- Samples can be stored for 24 hours at 4°C covered with 1X CS Wash Buffer.

Related product available from Trevigen.

Catalog#	Description	Size
3470-096-K	Cultrex [®] In Vitro Angiogenesis Assay Tube Formation Kit	96 tests

© 2010 Trevigen, Inc. All Rights Reserved. Trevigen and Cultrex are registered trademarks of Trevigen, Inc. E7/2/10v1

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com



Cell Staining Kit

Catalog#: 3437-100-K

TREVIGEN[®]

1-800-873-8443