

CULTREX[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

100X Phenol Red Solution

Catalog #: 3430-50-01

Size: 50 µl

Description: Trevigen provides 100X Phenol Red Solution as a supplement to our Cultrex[®] Basement Membrane Extract (BME) products.

Specifications:

Concentration: 1.5 mg/ml.

Storage buffer: Dulbecco's Modified Eagle's medium without phenol red, with 10 µg/ml gentamicin sulfate.

Storage: Store frozen, at -20°C or below.

Material Qualification:

Sterility testing:

- No bacterial or fungal growth detected after incubation at 37°C for 14 days following USP sterility testing guidelines.

References:

1. Berthois, Y., Katzenellenbogen, JA., and Katzenellenbogen BS. 1986. Phenol red in tissue culture media is a weak estrogen: Implications concerning the study of estrogen-responsive cells in culture. *Proc. Natl. Acad. Sci. USA* **83**: 2496-500.
2. Ernst, M., Schmid, C., and Froesch, ER. 1989. Phenol red mimics bio-logical actions of estradiol: Enhancement of osteoblast proliferation in vitro and of type I collagen gene expression in bone and uterus of rats in vivo. *J Steroid Biochem.* **33**:907-14.
3. Guy, R., Ridden, C., and Kealey T. 1996. The improved organ maintenance of the human sebaceous gland: Modeling in vitro effects of epidermal growth factor, androgens, estrogens, 13-cis retinoic acid, and phenol red. *J. Invest. Dermatol.* **106**:454-60.

Related Products:

Catalog#	Description	Size
3432-005-01	Cultrex [®] Basement Membrane Extract, PathClear [®]	5 ml
3433-005-01	Cultrex [®] Reduced Growth Factor BME, PathClear [®]	5 ml
3532-005-02	Cultrex [®] Basement Membrane Extract, Type 2, PathClear [®]	5 ml
3533-005-02	Cultrex [®] Reduced Growth Factor BME, Type 2, PathClear [®]	5 ml
3632-005-02	Cultrex [®] Basement Membrane Extract, Type 3, PathClear [®]	5 ml
3445-005-01	Cultrex [®] 3-D Culture Matrix [™] BME, PathClear [®]	5 ml
3446-005-01	Cultrex [®] 3-D Culture Matrix [™] Laminin I	5 ml
3447-020-01	Cultrex [®] 3-D Culture Matrix [™] Collagen I	100 mg
3434-005-02	Cultrex [®] Stem Cell Qualified RGF BME, PathClear [®]	5 ml
3415-001-03	Cultrex [®] Stem Cell Qualified Human BME, PathClear [®]	1 mg
3400-010-03	Cultrex [®] Stem Cell Qualified Laminin I, PathClear [®]	1 mg
3420-001-03	Cultrex [®] Stem Cell Qualified Human Fibronectin, PathClear [®]	1 mg
3420-001-03	Cultrex [®] Stem Cell Qualified Human Vitronectin, PathClear [®]	200 µg

© 2013 Trevigen, Inc. All Rights Reserved. Trevigen, Cultrex, CultreCoat and PathClear are registered trademarks, and 3-D Culture Matrix, DIVAA and CellSpere are trademarks of Trevigen, Inc. E8/15/13v1

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

Related Assays and Kits:

Catalog#	Description	Size
3500-096-K	Cultrex® 3D Spheroid Cell Invasion Assay	96 samples
3510-096-K	Cultrex® 3D Spheroid Fluorometric Proliferation/Viability Assay	96 samples
3511-096-K	Cultrex® 3D Spheroid Colorimetric Proliferation/Viability Assay	96 samples
3470-096-K	Cultrex® In Vitro Angiogenesis Assay, Tube Formation Kit	96 samples
3471-096-K	Cultrex® In Vitro Angiogenesis Assay, Endothelial Cell Invasion Kit	96 samples
3450-048-SK	Cultrex® Directed In Vivo Angiogenesis Assay (DIVAA™) Starter Kit	48 samples
3450-048-K	Cultrex® DIVAAT™ Kit	48 samples
3450-048-IK	Cultrex® DIVAAT™ Inhibition Kit	48 samples
3445-096-K	Cultrex® 3-D Culture BME Cell Proliferation Kit	96 samples
3448-020-K	Cultrex® 3-D Culture Cell Harvesting Kit	20 samples



100X Phenol Red Solution

Cat#: 3430-50-01
 Storage: ≤ -20°C
 (Manual Defrost)
 1-800-873-8443