

CULTREX[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Phenol Red

Catalog #: 3430-50-06

Size: 6 x 50 µl

Contents: Aqueous Phenol Red, Sterile (concentrated)

Description: Because phenol red is a weak estrogen mimetic, investigators working with estrogen receptor positive cells may want to control for phenol red-associated effects. Trevigen provides phenol red in separate vials for addition to our basement membrane extract (BME) products.

Instructions: Add 50 µl to 5 ml tube of BME (including PathClear and Growth Factor Reduced) for a 1X final working solution.

Performance: No microbial growth detected.

Storage Conditions: Store at 15 to 30°C protected from light. For long term storage, freeze in working aliquots at -80°C.

References: 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.

Ernst, M., Schmid, C., and Froesch, ER. 1989. Phenol red mimics biological 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.

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
3415-001-02	Cultrex [®] Human BME, PathClear [®]	1 ml
3445-048-01	Cultrex [®] 3-D Culture Matrix [™] BME	15 ml
3430-005-01	Cultrex [®] BME with Phenol Red	5 ml
3430-005-02	Cultrex [®] BME with Phenol Red, PathClear [®]	5 ml
3431-005-01	Cultrex [®] BME with Phenol Red; Reduced Growth Factor	5 ml
3431-005-02	Cultrex [®] BME with Phenol Red, Reduced Growth Factor PathClear [®]	5 ml
3438-100-01	Cultrex [®] Poly-L-Lysine	100 ml
3439-100-01	Cultrex [®] Poly-D-Lysine	100 ml
3437-100-K	Cultrex [®] Cell Staining Kit	100 ml
3450-048-05	CellSpense [™]	15 ml

© 2010 Trevigen, Inc. All Rights Reserved. Trevigen, Cultrex and PathClear are registered trademarks, and CellSpense is a trademark of Trevigen, Inc. P4/29/10v1

8405 Helgerman Court, Gaithersburg, MD 20877 USA

TREVIGEN[®]

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

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



Phenol Red

Catalog #: 3430-50-06

Storage: 15 to 30 °C

Protect from light

1-800-873-8443