

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Mounting Medium

Catalog #: 4865-25

Size: 25 ml

Description: Non-aqueous, low viscosity mounting medium for mounting coverslips and long-term storage and preservation of slides.

The formulation of this mounting medium, near the refractive index of fixed proteins, provides low optical interferences and excellent microscopic quality.

Low fading of color (most biological dyes) with age when slides are stored in darkness.

Physical State: Synthetic resin in a slightly viscous toluene based solution.

Color: Clear to very pale yellow.

Storage: Store at room temperature.

Applications: Ready to use with Trevigen's TACS® *in situ* hybridization DAB or Blue Label™ kits and other immunohistology procedures to mount and preserve samples on slides. Not compatible with fluorescent applications.

© 2011 Trevigen, Inc. All Rights Reserved. Trevigen and TACS are registered trademarks and Blue Label is a trademark of Trevigen, Inc. E3/16/11v1

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

Mounting Medium

Size: 25 ml

Catalog #: 4865-25

Storage: Room Temp.

TREVIGEN®

1-800-873-8443