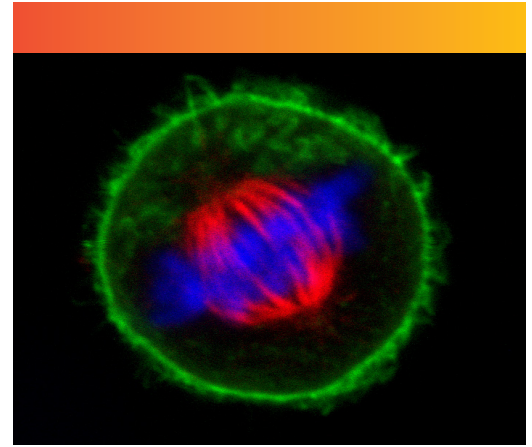




Biotium, Inc.  
 www.biotium.com  
 US orders: 1-800-304-5357



**New!**

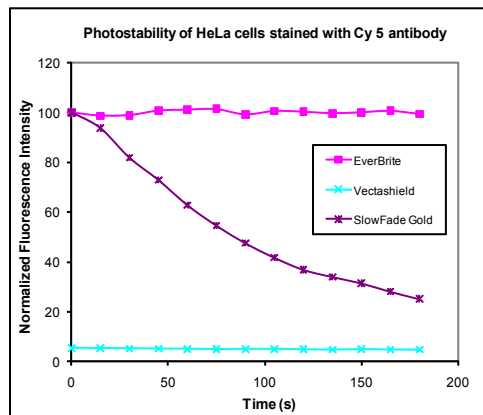
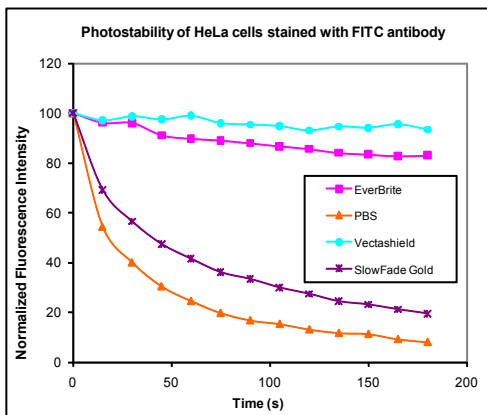
# EverBrite™ Mounting Medium

Biotium's new EverBrite™ mounting medium is an antifade mounting medium uniquely formulated for universally preserving fluorescence of all dyes across the entire visible and near-IR spectra. One major advantage of EverBrite™ mounting media is that they do not decrease the fluorescence of some of the cyanine-based fluorophores (e.g., Cy™2, Cy™5, Alexa Fluor® 647 and DyLight™ 649), such as that observed with Vectashield®. EverBrite™ mounting media may also be purchased with DAPI, which eliminates the need for a separate nuclear counterstaining step. After mounting, slides can be immediately viewed or sealed and stored at 4 °C for at least several months.

\*\*\*Please download EverBrite™ product information sheets from the website for more detailed information.

## FEATURES

- Preserves fluorescence over a wide range of dye colors
- Mounting medium also comes with DAPI to counterstain nuclei
- EverBrite™ mounting media permits immediate imaging of fluorescence and long-term preservation of slides
- Slides can be stored at 4 °C after sealing for at least several months



\*\*\* Vectashield is not shown as normalized data to illustrate the intensity drop of Cy™5 dye.

Cat.#	Product Name	Size
23001	EverBrite™ Mounting Medium	10 mL
23002	EverBrite™ Mounting Medium with DAPI	10 mL