

EverBrite™ Mounting Medium

A revolutionary antifade mounting medium optimally formulated for preserving fluorescence of our CF® dyes and other fluorochromes. Available with or without DAPI.



Product attributes

Format	Without DAPI, With DAPI
For live or fixed cells	For fixed cells/tissue sections
Probe cellular localization	Nucleus
Storage Conditions	Store at 2 to 8 °C, Protect from light

Product Description

Biotium's EverBrite™ mounting medium is a revolutionary antifade mounting medium optimally formulated for preserving fluorescence of our CF® dyes and other fluorochromes for your immunofluorescence (IF) imaging applications.

- Antifade mounting medium optimally formulated to prevent photobleaching of fluorophores
- Refractive index is well-matched to that of coverslip glass and immersion oil
- Available with or without DAPI

EverBrite™ provides a high level of protection from photobleaching to prolong fluorescence over a wide range of wavelengths. EverBrite™ Mounting Medium has a refractive index of 1.46, which is well-matched to the refractive indices of coverslip glass and immersion oil (1.5). The mounting medium remains liquid and permits immediate imaging of slides and tissues after sealing the cover slip edges with nail polish or Biotium's [CoverGrip™ Coverslip Sealant](#). After sealing, the slides can be stored at 4 °C for a year or longer. EverBrite is available with or without DAPI for nuclear counterstaining. Unlike VectaShield®, EverBrite™ is compatible with cyanine-based fluorophores like Cy® dyes.

We also offer [EverBrite™ Hardset Mounting Media](#) with or without DAPI. Also see our [NucSpot® Nuclear Stains](#) and our other [accessory products](#) for cell biology research.

EverBrite™ Mounting Media

Product	Counterstain	Cat. No.
EverBrite™ Mounting Medium	None	23001
EverBrite™ Mounting Medium with DAPI	DAPI	23002
Drop-n-Stain EverBrite™ Mounting Medium	None	23008
Drop-n-Stain EverBrite™ Mounting Medium with DAPI	DAPI	23009
EverBrite™ Hardset Mounting Medium	None	23003
EverBrite™ Hardset Mounting Medium with DAPI	DAPI	23004
EverBrite™ Hardset with NucSpot® 640	NucSpot® 640	23016
EverBrite TrueBlack® Hardset Mounting Medium	None	23017
EverBrite TrueBlack® Hardset with DAPI	DAPI	23018
EverBrite TrueBlack® Hardset with NucSpot® 640	NucSpot® 640	23019

EverBrite is a trademark of Biotium, Inc.; CF is a registered trademark of Biotium, Inc.; VectaShield is a registered trademark of Vector Laboratories. Cy Dye is a registered trademark of Cytiva.

This datasheet was generated on June 4, 2026 at 01:07:37 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/everbrite-mounting-medium/>