



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

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 December 29, 2025 at 04:39:52 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/everbrite-mounting-medium/>