

EverBrite™ Hardset Mounting Medium

A revolutionary antifade mounting medium optimally formulated for preserving fluorescence of our CF® dyes and other fluorochromes. EverBrite™ Hardset mounting medium hardens at room temperature completely in 24 hours.



Product attributes

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

Product Description

Biotium's EverBrite™ line of mounting media are a revolutionary antifade mounting media optimally formulated for preserving fluorescence of our CF® dyes and other fluorochromes. EverBrite™ Hardset mounting medium hardens at room temperature completely in 24 hours with excellent protection against photobleaching of fluorescent dyes. Unlike VectaShield®, EverBrite™ is compatible with cyanine-based fluorophores such as Cy®3 and Cy®5, Alexa Fluor® 555, Alexa Fluor® 647, and DyLight® 649.

- Antifade mounting medium optimally formulated to prevent photobleaching of fluorophores
- Hardens completely at room temperature in 24 hours
- Refractive index is well-matched to that of coverslip glass and immersion oil
- Cures to form a hard, permanent seal with the coverslip
- No need to seal coverslip edges with nail polish or other sealants
- Available with or without DAPI

EverBrite™ Hardset mounting medium cures to form a hard, permanent seal with the coverslip at room temperature in 24 hours. This facilitates handling of the slide and eliminates the need to seal coverslip edges with nail polish or other sealants. It is optimally formulated for use with Biotium's CF® dye conjugates but also provides a high level of protection from rapid photobleaching of fluorescein, rhodamine, and cyanine based dyes. Mounted slides can be stored at 4 °C for a year or longer. EverBrite™ Hardset Mounting Medium has a refractive index of 1.38 initially, which increases to 1.42 after 24 hours of curing, and 1.46 after four days of curing, after which it remains constant. EverBrite™ Hardset Mounting Medium is available with or without DAPI for nuclear counterstaining. Mounting media with DAPI gives a blue fluorescent nuclear counterstain without an additional staining step.

We also offer [EverBrite™ Mounting Media](#) in wet-set formulation, 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 VectorLabs; Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific; Cy and CyDye are registered trademarks of Cytiva.