

Drop-n-Stain EverBrite™ Mounting Medium

TOTAL COMMENTS OF THE PROPERTY OF THE PROPERTY

Wet-set antifade mounting medium for preserving fluorescence of our CF® dyes and other fluorochromes, in a dropper bottle for easy dispensing.

Product attributes

Call us: 800-304-5357

Format	Without DAPI, With DAPI
For live or fixed cells	For fixed cells/tissue sections
Probe cellular localization	Nucleus
Storage Conditions	Store at -10 to -35 °C

Product Description

Drop-n-Stain EverBriteTM Mounting Medium is a wet-set antifade mounting medium optimally formulated for preserving fluorescence of our CF® dyes and other fluorochromes. The dropper bottle makes for quick and easy dispensing.

- Convenient dropper bottle for pipette-free mounting
- Protects fluorescent dyes from rapid photobleaching
- Compatible with cyanine-based fluorophores like Cy®3, Cy®5, and Alexa Fluor® 647, unlike VECTASHIELD®
- Available with or without DAPI for nuclear counterstaining
- Works as well as original EverBrite™, but less viscous for rapid DAPI staining

Drop-n-Stain EverBrite™ Mounting Medium is a wet-set medium that remains liquid. After mounting, coverslips can be sealed with nail polish or Biotium's CoverGrip™ Coverslip Sealant (23005). Mounted slides can be stored at 4°C or -20°C, protected from light. Antibody-stained slides can be preserved for a year or longer (note: some labels, like phalloidin, are less stable during storage). Drop-n-Stain EverBrite™ Mounting Medium has a refractive index of 1.42.

We also offer <u>EverBriteTM Hardset Mounting Medium</u>, which cures overnight to form a hard seal, eliminating the need to seal the edges of the coverslips.

EverBrite™ Mounting Media

Product	Counterstain	Cat. No.
EverBrite TM Mounting Medium	None	23001
EverBrite™ Mounting Medium with DAPI	DAPI	23002
<u>Drop-n-Stain EverBrite™</u> <u>Mounting Medium</u>	None	23008
Drop-n-Stain EverBrite™ Mounting Medium with DAPI	DAPI	23009
EverBrite TM Hardset Mounting Medium	None	23003
EverBrite TM 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

Alexa Fluor is a registered trademark of Thermo Fisher Scientific; Cy Dye is a registered trademark of Cytiva; VECTASHIELD is a registered trademark of Vector Laboratories.

This datasheet was generated on November 22, 2025 at 11:08:34 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/drop-n-stain-everbrite-mounting-medium/.