

# Product Information

## EverBrite™ Mounting Medium EverBrite™ Mounting Medium with DAPI

### Catalog Numbers:

23001 EverBrite™ Mounting Medium

23002 EverBrite™ Mounting Medium with DAPI

**Unit Size:** 10 mL**Color and Form:** Light yellow to clear solution.

### Storage and Handling

Store at 4°C. Product is stable at least for 6 months from the date it is received. Warm to room temperature and mix well by gently swirling or inverting bottle before use; avoid vortexing or rapid shaking, which may introduce air bubbles. Store in an upright position and tightly cap after each use. Protect from light.

### Product Description

EverBrite™ Mounting Medium is a unique, stable formula for preserving fluorescence. EverBrite™ Mounting Medium prevents rapid photobleaching of fluorescein, rhodamine, and cyanine based dyes and is optimally formulated for use with our CF™ dye conjugates. EverBrite™ Mounting Medium is available with or without the blue fluorescent DNA dye DAPI for nuclear counterstaining. EverBrite™ Mounting Medium and EverBrite™ Mounting Medium with DAPI both have a refractive index of 1.46, which is well-matched to the refractive indices of coverslip glass and immersion oil (1.5).

To mount tissues or cells on a slide, remove as much excess moisture as possible, and then dispense approximately 10 - 20  $\mu$ L of the mounting medium depending on the size of the coverslip. Coverslip and allow EverBrite™ Mounting Medium to spread over the entire section. Applying an excess of mounting medium is not recommended. Excess mounting medium can be carefully aspirated or blotted away. Slides can be view immediately after mounting.

EverBrite™ Mounting Medium remains liquid. After mounting, coverslips can be sealed around the perimeter with nail polish or Biotium's CoverGrip™ Coverslip Sealant. Mounted slides should be stored at 4°C protected from light. Slides can be preserved for a year or longer.

EverBrite™ Mounting Medium is superior to Vectashield® and SlowFade® Gold. For example, EverBrite™ does not dissipate or weaken some of the cyanine-based fluorophores, such as that observed with Vectashield®. Please note: EverBrite™ Mounting Medium may not be compatible with all fluorescent proteins or dyes or with certain tissue sections. We recommend using EverBrite™ on a test sample to ensure compatibility with tissue sections or with uncommon fluorochromes.

Biotium also offers EverBrite™ Hardset Mounting Medium, which cures overnight to form a hard seal, eliminating the need to seal the edges of the coverslip (see Related Products).

### Related Products

Catalog Number	Product
40061	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO
40043	DAPI in H <sub>2</sub> O, 10 mg/mL
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher, 20X in DMF
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer
22010	10% Fish Gelatin Blocking Buffer
22011	Fish Gelatin Powder
22014	30% Bovine Serum Albumin Solution
22002	Tween®-20

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including fluorescent CF™ dye secondary antibodies and other conjugates, antibody labeling kits, apoptosis reagents, fluorescent probes, and kits and accessories for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

EverBrite is a trademark of Biotium, Inc.; Vectashield is a trademark of Vector Labs; SlowFade is a trademark of Invitrogen, Corp., Tween is a registered trademark of Uniqema Americas LLC.