

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 before use. EverBrite™ is supplied in an amber bottle to protect it from light. Store in an upright position and tightly cap after each use.

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.

To mount tissues or cells on a slide, remove as much excess moisture as possible, and then dispense approximately 10 - 20 µL 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 at least several months.

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.

Related Products

Catalog Number	Product Name	Unit Size
23005	CoverGrip™ Coverslip Sealant	15 mL
23005-1	CoverGrip™ Coverslip Sealant Refill	100 mL

Please visit our website at www.biotium.com to view our full selection of bioconjugates of our exceptionally bright and photostable CF™ dyes, including antibodies, antibody labeling kits, phalloidin, annexin V, nucleotides, and a-bungarotoxin, as well as classic fluorescent nucleic acid dyes.

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