

Product Information

EverBrite™ Hardset Mounting Medium and EverBrite™ Hardset Mounting Medium with DAPI

Catalog Numbers:

23003: EverBrite™ Hardset Mounting Medium

23004: EverBrite™ Hardset 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 for at least 6 months from the date it is received when stored as recommended. 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™ Hardset Mounting Medium is an anti-fade mountant for preserving fluorescence during microscopy. EverBrite™ Hardset Mounting Medium prevents rapid photobleaching of fluorescein, rhodamine, and cyanine based dyes and is optimally formulated for use with Biotium's CF™ dye conjugates. Unlike Vectashield, EverBrite™ mounting media is compatible with Cy™ dyes such as Cy™2 and Cy™5. Please note: EverBrite™ Hardset Mounting Medium may not be compatible with all fluorescent proteins or dyes and test sections are recommended to ensure compatibility with uncommon fluorochromes.

EverBrite™ Hardset Mounting Medium cures to form a hard, permanent seal with the coverslip, eliminating the need to seal coverslip edges with nail polish or other sealants. EverBrite™ Hardset Mounting Medium with DAPI allows blue fluorescent nuclear counterstaining without an additional staining step. EverBrite™ Hardset Mounting Medium and EverBrite™ Hardset Mounting Medium with DAPI both have 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.

To mount tissue sections or cells on a slide, dispense approximately 10 - 20 uL of solution depending on the size of the coverslip. Then coverslip and allow EverBrite™ Hardset 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 viewed immediately after mounting, but the coverslip will not be secured in place, so care must be taken to avoid movement of the coverslip. To cure the mounting medium to form a permanent hard seal, incubate slides overnight at room temperature on a flat surface protected from light. Cured slides can be stored at 4°C protected from light for a year or longer.

Coverslips can be removed from cured slides by soaking the slide in PBS or similar buffer in a slide staining jar. Soak the slides for 2.5-3 hours, then gently slide the coverslip off (do not pry the coverslip upwards), dipping the slide frequently in buffer to prevent friction between the coverslip and specimen.

Biotium also offers EverBrite™ wet set mounting media (see Related Products).

Related Products

Catalog Number	Product
40061-T	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO
40043	DAPI in H ₂ O, 10 mg/mL
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
22005	Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses
22006	Super ^{HT} 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 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.