



Revised: May 24, 2016

Product Information

Related Products

CoverGrip™ Coverslip Sealant

Catalog Number: 23005, 23005-1

Unit Size: 23005: 15 mL brush cap applicator bottle 23005-1: 100 mL dropper cap refill bottle

Color and Form: Yellow solution

Storage and Handling

Store at room temperature. Avoid contact with skin and eyes. Flammable; keep away from open flame. For laboratory use only.

Product Description

CoverGrip™ Coverslip Sealant is an organic adhesive medium formulated specifically for sealing the edges of coverslips mounted using wet set antifade mounting medium for immunofluorescence. Traditionally, nail polish has been used to seal the edges of coverslips prior to fluorescence microscopy, to secure the coverslip in place, prevent seepage of mounting medium during imaging, and to keep slides from drying out during storage. However, components in nail polish such as isopropyl alcohol can leach into wet set mounting medium and interfere with fluorescence imaging. In particular, GFP fluorescence is reported to be shown to be sensitive to nail polish used to seal coverslips (1). In addition, because nail polish is not designed to be an adhesive, different brands of nail polish may not provide consistent sealing. Alternative methods of sealing coverslips with paraffin or rubber cement are messy and do not generate a permanent hard seal. Cover-Grip™ Coverslip Sealant is easy to apply, is not miscible with aqueous mounting media, and dries to form clear, hard seal along the edges of the coverslip. Sealing with CoverGrip™ Coverslip Sealant allows coverslips to be cleaned before imaging, prevents coverslip slippage, protects microscope objectives from exposure to mounting medium during imaging, and provides a long lasting hard seal for storing wet-mounted specimens at 4°C or -20°C.

CoverGrip™ Coverslip Sealant is made with limonene, a pleasant-smelling natural hydrocarbon made from citrus peels that is used as a more environmentally friendly and less toxic alternative to organic solvents.

CoverGrip™ Coverslip sealant is supplied in a 15 mL brush cap bottle for ease of application, and also is available in 100 mL dropper-top refill bottles.

Instructions for use

Use CoverGrip™ Coverslip sealant the same way you would use nail polish to seal coverslips. Swirl bottle to mix. Mount coverslips using antifade mounting medium. Carefully aspirate or blot all excess mounting medium from the coverslip edges. When the coverslip edges are completely dry, use the brush attached to the cap to apply CoverGrip™ Coverslip Sealant along all edges of the coverslip, completely overlapping the coverslip edges. Allow CoverGrip™ Coverslip Sealant to dry completely at room temperature, about 1 hour. Drying time can be shortened by drying slides under a fan. When CoverGrip™ Coverslip Sealant is completely dry, gently clean the surface of the coverslip using a low-lint wiper and window cleaner or 70% alcohol. Store slides at 4°C or -20°C.

Like nail polish, CoverGrip™ can break down when exposed to excessive moisture or condensation. Store slides with dessicant in humid storage environments (e.g. deli refrigerators and cold rooms). CoverGrip™ Coverslip Sealant can be removed from slides by soaking slides in PBS or similar buffer for 1-2 hours. The CoverGrip™ Coverslip Sealant will not dissolve, but will loosen to a degree that it can be peeled or gently scraped off of the slide using coverslip forceps.

References

1. Chalfie, M. et al. (1994). Science 263, 802-805.

Catalog Number	Product
40061	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
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
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.

Tween is a registered trademark of Uniqema Americas LLC.