



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

CoverGrip™ Coverslip Sealant

Seals the edges of wet-mounted coverslips for fluorescence microscopy, with no ingredients that can leach into aqueous mounting medium and affect specimen fluorescence.



Product attributes

Storage Conditions
Store at room temperature

Product Description

CoverGrip™ Coverslip Sealant is the first product specifically designed for sealing the edges of wet-mounted coverslips for fluorescence microscopy. Unlike nail polish, CoverGrip™ contains no ingredients that can leach into aqueous mounting medium and affect specimen fluorescence.

- Seals the edges of wet-mounted coverslips for fluorescence microscopy
- Dries hard and clear
- Less damaging to samples than nail polish
- Won't leach into aqueous mounting medium and affect fluorescence
- Contains d-limonene

CoverGrip™ is made with d-limonene, a natural, environmentally-friendly solvent with a pleasant citrus aroma, and dries to form a clear, hard, durable seal along the edges of the coverslip. Available in 15 mL brush-cap applicator bottles or in 100 mL dropper-cap refill bottles. CoverGrip™ is designed to replace nail polish for permanent sealing of wet-mounted coverslips. It can be removed by soaking slides in buffer (see product information sheet for details). However, it is not designed to be a removable temporary seal for hybridization, for example.

This datasheet was generated on January 16, 2026 at 04:46:25 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/covergrip-coverslip-sealant/>