



# SEEBRIGHT®

## Mounting Medium

**Antifade mounting medium minimizes photobleaching of most fluorophores, dyes and fluorescent proteins.**

SEEBRIGHT® Mounting Medium is an antifade reagent that minimizes photobleaching. Use of this mounting medium significantly extends the lifetime of fluorescent signals.

All fluorescent dyes signal fade (photobleach) when exposed to excitation light. The amount of photobleaching depends on the composition of the molecule and the time of light exposure. SEEBRIGHT® Mounting Medium can be used to prevent photobleaching and prolong fluorescent signals. SEEBRIGHT® Mounting Medium is compatible with many dyes, including Enzo's SEEBRIGHT® Fluorescently-Labeled dUTPs.

### Ordering Information

[Order Online »](#)

ENZ-GEN422-1000	10ml
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)

- Antifade formulation inhibits rapid photobleaching of fluorescent proteins and fluorescent dyes
- No added counterstain allows for custom titration of DAPI as desired
- Ready-to-use non-hardening formulation
- Refractive index: 1.45

## Handling & Storage

Handling	Protect from light. Do not freeze.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

### Quantity

Each vial of SEEBRIGHT<sup>®</sup> Mounting Medium contains 10 mL sufficient for 1000 tests (ENZ-GEN-420-1000). Apply mounting medium to the sample area as desired.

### Technical Info / Product Notes

SEEBRIGHT<sup>®</sup> Mounting Medium is recommended for use on fixed cells and tissues that have been stained with fluorescent dyes or hybridized with fluorescent probes. Cells can be viewed immediately or mounted slides can be stored in the dark at 4°C. SEEBRIGHT<sup>®</sup> Mounting Medium is a non-hardening formula. Hence, slides with coverslips can be stored unsealed or coverslips can be sealed around their perimeter using nail polish or a plastic sealant.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)