

# SEEBRIGHT® Mounting Medium with DAPI

## Antifade mounting medium with DAPI counterstain minimizes photobleaching of

SEEBRIGHT® Mounting Medium with DAPI is an antifade reagent that minimizes photobleaching. Use of the mounting medium significantly extends the lifetime of fluorescent signals. The inclusion of DAPI allows for staining of the nucleus.

All fluorescent dyes 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 with DAPI can be used to prevent photobleaching and prolong fluorescent signals. This formulation contains 4',6-diamidino-2-phenylindole (DAPI), which serves as a nuclear or chromosomal counterstain that produces a blue fluorescence when bound to DNA. SEEBRIGHT® Mounting Medium with DAPI is compatible with many dyes, including Enzo's SEEBRIGHT® Fluorescently-Labeled dUTPs.

## Ordering Information

[Order Online »](#)

ENZ-GEN420-0100	100 tests
-----------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

- Unique antifade formulation inhibits rapid photobleaching of fluorescent proteins and fluorescent dyes
- Ready-to-use non-hardening formulation
- Includes DAPI counterstain at 1.5 µg/ml

## 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 with DAPI contains 200 µL sufficient for 20 tests (ENZ-GEN420-0020) or 1000 µL sufficient for 100 tests (ENZ-GEN420-0100). Apply mounting medium to the sample area as desired.

### Technical Info / Product Notes

SEEBRIGHT<sup>®</sup> Mounting Medium with DAPI 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 stored in the dark at 4°C. SEEBRIGHT<sup>®</sup> Mounting Medium with DAPI 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.

DAPI excites at about 360 nm and emits at about 460 nm. DAPI is a mutagen and appropriate precautions should be taken.

Last modified: September 15, 2025



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)