

Antifade mounting medium with DAPI (hardset)

Antifade reagent for use in multiple applications

Antifade mounting medium with DAPI (hardset) is an antifade reagent that causes little or no quenching. Use of the mounting medium significantly extends the photobleaching half-life of the fluorescent stain. 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 dye, the intensity and the time of illumination. Many products have been developed to prevent fading of fluorescent dyes. These products often work well for one or more dyes, but often cause problems, such as degradation of other dyes. Antifade mounting medium with DAPI (hardset) works well with many stains including fluorescein and rhodamine.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-53003-M010	10ml
----------------	------

Manuals, SDS & CofA

[View Online »](#)

- Inhibits photobleaching of fluorescent dyes and fluorescent proteins
- Hardening formulation
- Includes DAPI counterstain in formulation

Handling & Storage

Handling	Protect from light. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Technical Info / Product Notes

The Antifade mounting medium with DAPI (hardset) is recommended for use on fixed cells and tissues that have been stained with fluorescent stains. Cells can be viewed immediately, or if long term storage is desired, the reagent can be allowed to harden overnight.

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)