



HIGHDEF[®]

Hematoxylin

Ready-to-use histologic stain with enhancers

HIGHDEF[®] Hematoxylin is intended for use in the histologic demonstration of nuclear staining. This staining technique is sometimes used to make the critical distinction between a normal nucleus and an abnormal one.

Citations: 7

[View Online »](#)

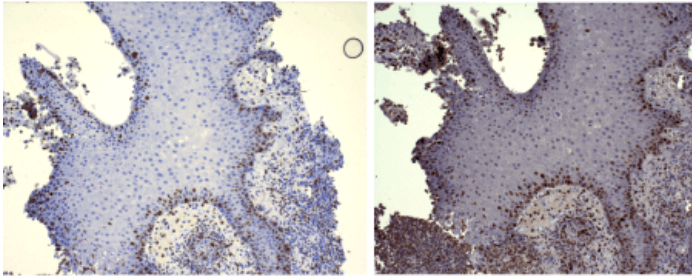
Ordering Information

[Order Online »](#)

ENZ-ACC106-0100	100ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Enzo Hematoxylin

Competitor Hematoxylin

Enzo's hematoxylin allows for clearer staining and less background than a leading competitor. Both slides were stained with Enzo's POLYVIEW® PLUS HRP-DAB Kit, but the competitor's hematoxylin was used on the slide on the right, showing more background and a less clear result.

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Formulation Enhancers added to improve signal.

Technical Info / Product Notes

Hematoxylin is one of the most common staining compounds used in histology.

The protocols for specific applications can vary. These include, but are not limited to: fixation, heat-retrieval methods, incubation times, tissue section thickness, and detection kit used. Due to the superior sensitivity of these unique reagents, the recommended incubation times and titers listed are not applicable to other detection systems, as results may vary. Ultimately, it is the responsibility of the investigator to determine optimal conditions. These products are tools that can be used for interpretation of morphological findings in conjunction with other diagnostic tests and pertinent clinical data by a qualified pathologist.

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)