

HIGHDEF[®] IHC mount

Permanent, high stability mounting medium
for high definition visualization of IHC.

HIGHDEF[®] IHC mount is an aqueous-based permanent mounting medium designed for use with chromogenic labeling. Use for mounting of tissue specimens stained with peroxidase and alkaline phosphatase-based systems and various fluorescent dyes.

AEC and Fast Red are the most commonly used red chromogens for peroxidase and alkaline phosphatase based immunostaining systems. Slides stained with these chromogens cannot be stored permanently because of the lack of availability of an aqueous permanent mounting medium with high refractive index. HIGHDEF[®] IHC mount is designed to fill in these limitations and it is an aqueous mounting medium with a very high refractive index. It allows permanent storage of tissue specimens without fading of the chromogens. Because of its superior refractive index, tissues mounted in this medium look like dehydrated specimens. No coverslipping is required. If coverslipping is required, dried HIGHDEF[®] mount can be post mounted using an organic-based mounting medium. Advantages of HIGHDEF[®] mount are numerous: no coverslip is required, there is no exposure to organic fumes, it allows permanent storage of slides, and it provides a high resolution of tissue specimens. HIGHDEF[®] IHC mount is compatible with chromogens like AEC, DAB, Fast Red, BCIP/NBT, BCIP/INT, and fluorescent dyes like FITC and phycobiliproteins. HIGHDEF[®] IHC mount's high pH ensures increased stability of fluorescence.

Application Note

[Tumor Phenotyping and Immune Cell Regulation using Multiplex IHC](#)

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-950-261-0030

30ml

Manuals, SDS & CofA

[View Online »](#)

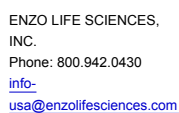
- Aqueous-based permanent mounting medium designed for chromogenic use and ideal for long-term storage of processed tissue
- Flexible formulation optimized for AEC, DAB, Fast Red, BCIP/NBT, and BCIP/INT
- High definition visualization for IHC, IF, ISH, FISH

Long Term Storage Ambient

Regulatory Status RUO - Research Use Only

Formulation	Liquid. Mounting medium containing <0.1% sodium azide.
--------------------	--

Last modified: May 29, 2024



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)