

# HIGHDEF<sup>®</sup> IHC fluoromount

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

HIGHDEF<sup>®</sup> IHC fluoromount is for mounting tissue specimens stained with fluorescent dyes. Designed for use in fluorescence, but can be used with other applications.

Citations: 42

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-950-260-0025	25ml
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

- Aqueous-based permanent mounting medium designed for fluorogenic labeling and ideal for long-term storage of processed tissue
- High definition visualization for IHC, IF, ISH, FISH

## Handling & Storage

Long Term Storage                      Ambient

Shipping                                      Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

### Formulation

Liquid. Mounting medium containing <0.1% sodium azide.

### Technical Info / Product Notes

HIGHDEF<sup>®</sup> IHC fluoromount is an aqueous-based mounting medium for applying coverslips to tissues stained with fluorescent dyes. This product is compatible with FITC, phycoerythrin, phycocyanin, and allophycocyanin. It can also be used with other fluorescent markers such as Texas Red and rhodamine.

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)