

# IHC/ISH Peroxidase Block

Use IHC/ISH Peroxidase Block to block nonspecific background staining observed during immunostaining.

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-ACC107-0100	100ml
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Technical Info / Product Notes** The IHC/ISH Peroxidase Block is a unique serum-free blocking solution used to reduce nonspecific background staining. It is a universal blocking agent which eliminates the need to match species with the link antibody. This blocking reagent has been shown to be superior to serum-based blocking solutions. In addition to immunohistochemistry, the IHC background blocker can also be used for ISH.

Interpretation of the results will be the sole responsibility of the user.



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)