

# Antigen Retrieval Reagent, pH 6 (10X)

Antigen Retrieval Reagent, pH 6 (10X) is used to recover antigens masked by fixation in cross linking fixatives such as formalin.

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-ACC112-0100	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

**Formulation** Liquid. In citrate buffer.

### Technical Info / Product Notes

In order to perform immunostaining, tissue specimens should be preserved in an appropriate fixative. Fixation stops tissue autolysis, preserves tissue structures, and immobilizes antigens. However, antigens undergo chemical alteration of their primary, secondary and tertiary structures during fixation. Antigenic sites may be masked due to changes induced in the epitope or neighboring proteins. Enzymatic treatment with proteolytic enzymes (i.e. pepsin, trypsin or pronase) has been performed to expose the masked antigens.

Shi et al. (1991) have reported that treating tissue sections with a heavy metal solution in a microwave oven can cover masked antigens significantly. However, heavy metals in the solution increase the risk of toxic exposure to lab personnel. To reduce the risk, we have developed an antigen unmasking solution which is free of heavy metals. The reagent is provided in a convenient 10X concentrate. Use of this antigen recovery buffer avoids unnecessary heavy metal exposure to lab personal and other handling and disposal issues.

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

**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-fr@enzolifesciences.com](mailto:info-fr@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)