

IHC enzyme antigen retrieval reagent

IHC enzyme antigen retrieval reagent is designed to unmask immunoreactive sites altered by fixation and/or the embedding process.

The most commonly used tissue fixatives are aldehydes, such as formalin, which work by protein cross-linking. Proper fixation of a specimen prevents antigen from being washed out and is a critical aspect of successful IHC staining. Cross-linking can mask antigens, preventing primary antibodies from binding, thus yielding weak staining or false negative results. Immunoglobulins are especially susceptible to formalin fixation. Treatment with a proteolytic enzyme is required to digest excess aldehyde linkages and expose the antigen. The IHC enzyme antigen retrieval reagent is a proteolytic enzyme solution that results in superior staining over traditionally used enzyme pretreatments, such as pepsin and trypsin. Moreover, the IHC enzyme antigen retrieval reagent requires only a brief incubation at room temperature. Performing antigen recovery with the IHC enzyme antigen retrieval reagent will allow you to dilute the primary antibody 5 to 10 fold further and achieve optimal staining.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-950-280-0015	15ml
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Clear slightly viscous solution.

Formulation Liquid. Enzyme antigen retrieval reagent.

Technical Info / Product Notes 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)