

IHC tissue primer

The IHC tissue primer is a universal enzyme blocking reagent to use in horse radish peroxidase (HRP) and alkaline phosphatase (AP) based immunohistochemical staining procedures done on paraffin tissue sections, frozen tissue sections and cell smears to suppress nonspecific staining due to endogenous enzyme activity.

In the evaluation of tissue specimens by IHC using HRP and AP methods, the presence of endogenous enzymes can often obscure the specific staining of the target antigen. Endogenous peroxidase and AP activity can be inhibited by the use of the IHC tissue primer which will suppress endogenous peroxidase, pseudoperoxidase, and alkaline phosphatase in paraffin tissue sections, frozen tissue sections and cell smears.

Ordering Information

[Order Online »](#)

ADI-950-234-0100	100ml
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Store tightly sealed in the dark. Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid. Tissue primer. TRIS-based alkaline solution with 30% H₂O₂.

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)