

MULTIVIEW® PLUS

Antigen Retrieval Reagent (10X) (For Protease-Free

Workflow)

MULTIVIEW® Antigen Retrieval Reagent (10X) is a high-pH, heat-based epitope and nucleic acid target retrieval solution formulated for **protease-free ISH and IHC workflows** on formalin-fixed paraffin-embedded (FFPE) tissue sections. Unlike conventional ISH protocols that require Proteinase K (protease) digestion to unmask nucleic acid targets, MULTIVIEW® achieves effective target retrieval by preserving protein epitopes for **simultaneous IHC detection**. This makes it the ideal pre-treatment reagent for **ISH/IHC multiplex assays**, enabling visualization of both protein (antibody-based IHC) and nucleic acid (RNA or DNA ISH) targets in the same FFPE tissue section without protease treatment between staining steps. Supplied as a 10X concentrate; dilute 1:10 with deionized water before use. Compatible with AMPIVIEW® RNA ISH (LoopRNA ISH™ technology) and Enzo IHC detection workflows.

Formalin fixation forms protein–protein and protein–nucleic acid cross-links that mask both antibody epitopes and hybridization targets. Traditional ISH protocols use Proteinase K digestion to restore accessibility of nucleic acid targets; however, protease treatment degrades protein epitopes, making subsequent IHC on the same tissue section unreliable or impossible. Heat-induced antigen/target retrieval using a high-pH buffered solution reverses cross-linking through a protease-independent mechanism, unmasking both epitopes and hybridization targets simultaneously. This protease-free approach preserves protein antigenicity, enabling true ISH/IHC multiplexing — the simultaneous or sequential detection of RNA/DNA targets by ISH and protein markers by IHC on the same tissue section without the trade-off imposed by protease digestion.

Ordering Information

[Order Online »](#)

ENZ-ACC185-0100

100ml

Manuals, SDS & CofA

[View Online »](#)

- **Protease-free** — no Proteinase K required; protein epitopes preserved for IHC
- Enables **ISH/IHC multiplexing**: visualize protein and nucleic acid targets simultaneously on the same FFPE section
- Optimal for workflows requiring both RNA/DNA detection (ISH) and protein detection (IHC) without protease steps between staining layers
- High-pH heat-induced retrieval unmask epitopes and hybridization targets in a single step
- Concentrated 10X format; dilute 1:10 for economical, consistent use
- Compatible with AMPIVIEW® ISH (LoopRNA ISH™) and Enzo POLYVIEW® IHC detection workflows

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label.
Handling	Dilute 1:10 with deionized water before use; discard working solution after each use.
Short Term Storage	2–8°C
Long Term Storage	2–8°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Contents MULTIVIEW® PLUS Antigen Retrieval Reagent (10X) (ENZ-ACC185)

Technical Info / Product Notes Dilute 1:10 with deionized water to prepare 1X working solution. Without protease digestion between steps, enabling simultaneous visualization of protein and nucleic acid targets. Optimal retrieval conditions should be verified per tissue type and target combination.



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)