

POLYVIEW[®] IHC reagent (mouse/rabbit-HRP)

High sensitivity, low background nanopolymer detection reagent for use with

HIGHDEF[®] chromogens.

The POLYVIEW[®] (mouse/rabbit-HRP) IHC detection reagent is a non-biotin one-step detection reagent suitable for demonstrating antigens in formalin-fixed paraffin-embedded tissues and frozen sections. The HRP detection reagent may also be used with blood smears, cytospins, and cell preparations.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-950-113-0100	100 tests
------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Biotin-free nanopolymer detection circumvents endogenous biotin background
- High intensity color development delivers sharp, crisp staining
- Customize with HIGHDEF[®] chromogens of your choice (sold separately)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC

Contents Anti-Mouse/Rabbit HRP detection reagent

Technical Info / Product Notes POLYVIEW[®] IHC detection reagent (mouse/rabbit-HRP) has been developed by directly labeling anti-mouse and anti-rabbit immunoglobulins with enzymes using a proprietary tandem hyperlabelling technology. This ensures consistent and reproducible immunodetection of mouse and rabbit antibodies with a single reagent. Nuclear, cytoplasmic and membrane antigens in different types of tissues can be detected readily. The POLYVIEW[®] IHC detection reagent (mouse/rabbit-HRP) enables faster staining procedures than traditional two-step methods using biotin and avidin/streptavidin conjugates, with significantly lower background. The kit is suitable for use with all mouse and rabbit antibodies, both monoclonal and polyclonal. The reagents can be used for manual staining or with automated staining platforms.



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)