

MULTIVIEW® (mouse-HRP/rabbit-AP) IHC kit

Superior multiplex IHC kit for efficient detection of multiple antigens in a single tissue sample.

The MULTIVIEW® (mouse-HRP/rabbit-AP) IHC kit is a non-biotin one-step detection system suitable for demonstrating antigens in formalin-fixed paraffin-embedded tissues and cryostat sections. The kit may also be used with blood smears, cytosmears, and cell preparations.

Citations: 19

[View Online »](#)

Ordering Information

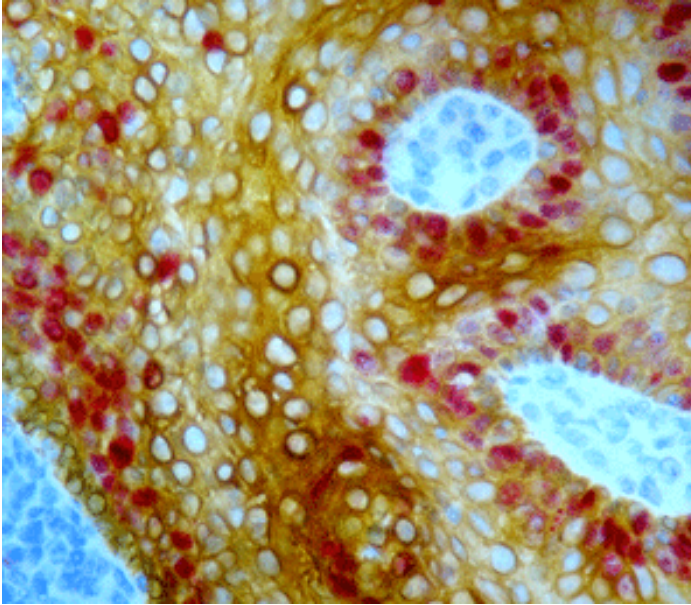
[Order Online »](#)

ADI-950-100-0001	100 tests
------------------	-----------

Manuals, SDS & CofA

[View Online »](#)

- Biotin-free nanopolymer-based system for superior multiplex antigen detection
- Complete kit includes HIGHDEF® Chromogens for sensitive, crisp 2-color development
- Preserves precious tissue while reducing reagent and labor costs
- Expandable with additional chromogens (sold separately)



Immunohistochemical analysis: Formalin fixed paraffin embedded human tonsil double stained using MULTIVIEW[®] (mouse-HRP/rabbit-AP) IHC kit (Prod. no. ADI-950-100). Primary rabbit anti-Ki67 antibody was visualized with HIGHDEF[®] red IHC chromogen (AP) (Prod. no. ADI-950-140), while primary mouse anti-HMW CK antibody was visualized with HIGHDEF[®] yellow IHC chromogen (HRP) (Prod. no. ADI-950-170, sold separately), producing a distinct golden yellow color contrasting with the bright red.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC

Assay Time ~2 hours

Contents Peroxidase Block, POLYVIEW[®] IHC reagent (mouse-HRP), POLYVIEW[®] IHC reagent (rabbit-AP), HIGHDEF[®] IHC chromogen substrate (DAB, stable) buffer & chromogen, HIGHDEF[®] red IHC chromogen (AP) buffer & chromogen, Empty mixing bottles

Technical Info / Product Notes

The MULTIVIEW[®] (mouse-HRP/rabbit-AP) IHC kit has been developed by directly labeling anti-rabbit and anti-mouse 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 readily detected. The MULTIVIEW[®] (mouse-HRP/rabbit-AP) IHC kit enables faster staining procedures than traditional two-step methods using biotin and avidin/streptavidin conjugates by being able to directly label mouse antibodies with HRP and rabbit antibodies with AP at the same time, all with significantly lower background staining. The kit is suitable for use with mouse and rabbit antibodies, both monoclonal and polyclonal. The reagents can be used for manual staining or with automated staining instruments and is specifically designed for multiplex immunohistochemical staining assays.

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-
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)