

MULTIVIEW[®] PLUS

(mouse-HRP/rabbit-AP) IHC Kit

(Brown/Green)

The MULTIVIEW[®] PLUS (mouse-HRP/rabbit-AP) IHC Kit is a non-biotin one-step detection system suitable for detecting and localizing antigens in formalin-fixed paraffin-embedded tissues.

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

Application Note

[Three-color IHC Staining On Free-floating Sections From Human Brain Tumor Phenotyping and Immune Cell Regulation using Multiplex IHC](#)

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

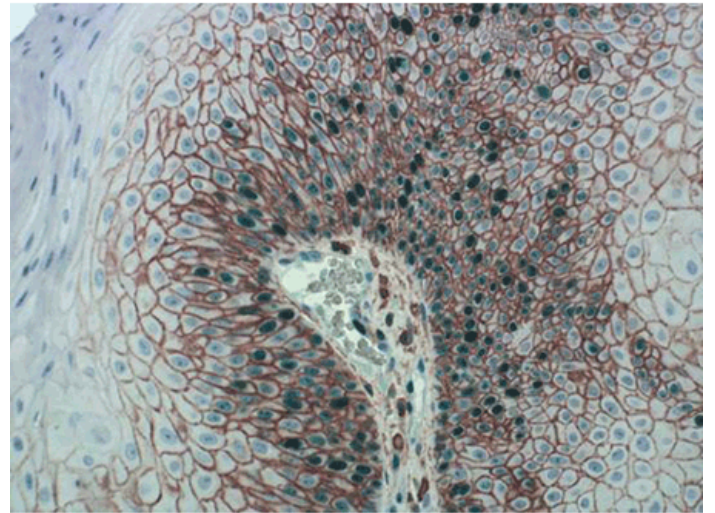
ENZ-KIT181-0150

150 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)



Mouse anti-CD44 clone SFF-304 and mouse anti-Ki-67 clone MIB-1 with Rabbit anti-Mouse Linker (Cat. No. ENZ-ACC116) was used with MULTIVIEW[®] PLUS (mouse-HRP/rabbit-AP) Kit (Prod. No. ENZ-KIT181) to perform multiplex immunohistochemistry staining with hematoxylin used as a counterstain. Tissue used is human buttock.

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 Antigen Retrieval Reagent, pH 9 (10X), ISH/IHC Peroxidase Block, Antibody Blocker/Diluent, MULTIVIEW[®] PLUS HRP (Anti-Mouse) Reagent, MULTIVIEW[®] PLUS AP (Anti-Rabbit) Reagent, HIGHDEF[®] DAB Chromogen, HIGHDEF[®] DAB Substrate, HIGHDEF[®] Green AP Chromogen, HIGHDEF[®] Green AP Substrate, Hematoxylin, Empty Mixing Bottle

Technical Info / Product Notes

The MULTIVIEW[®] PLUS (mouse-HRP/rabbit-AP) IHC Kit (Brown/Green) provides reagents and materials for the ultrasensitive multiplex detection of mouse and rabbit antibodies used in immunohistochemistry (IHC) staining. The kit is a non-biotin detection system suitable for identifying antigens in formalin-fixed paraffin-embedded tissues and frozen sections. This kit may also be used with blood smears, cytosmeas, and cell preparations.

MULTIVIEW[®] PLUS (mouse-HRP/rabbit-AP) IHC Kit (Brown/Green) contains both MULTIVIEW[®] PLUS HRP (Anti-Mouse) Reagent and MULTIVIEW[®] PLUS AP (Anti-Rabbit) Reagent, which have been developed by directly labeling anti-mouse immunoglobulins with enzymes using a proprietary tandem hyper-labelling technology in a ready-to-use format. This ensures consistent and reproducible immunodetection of mouse and rabbit antibodies against nuclear, cytoplasmic, and membrane antigens in different types of tissues and cells. The MULTIVIEW[®] PLUS HRP (Anti-Mouse) Reagent and MULTIVIEW[®] PLUS AP (Anti-Rabbit) Reagent enable faster staining procedures compared to traditional two-step methods using biotin and avidin/streptavidin conjugates, with significantly lower background.



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)