



Ultra-sensitive Multiplex Detection

MULTIVIEW® PLUS (mouse-HRP/rabbit-AP) IHC Kit (Brown/Green)

Superior Multiplex Detection

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.

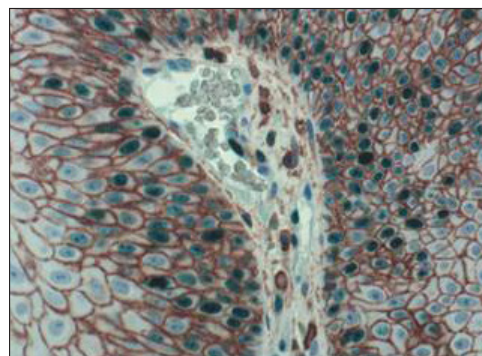
MULTIVIEW PLUS (mouse-HRP/rabbit-AP) IHC Kit (Brown/Green) contains both MULTIVIEW PLUS HRP (Anti-Mouse) Reagent and MULTIVIEW PLUS AP (Anti-Rabbit) Reagent, which have been developed by directly labeling anti-mouse immunoglobulins with enzymes using a proprietary labeling technology. This ensures consistent 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.

Key Features

- 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)

Multiplex IHC Membrane and Nuclear Proteins



Mouse anti-CD44 clone SFF-304 and Rabbit anti-Ki-67 clone MIB-1 was used with MULTIVIEW PLUS (mouse-HRP/rabbit-AP) IHC Kit (Brown/Green) to perform multiplex IHC.

PRODUCT NAME	SIZE	PRODUCT #
MULTIVIEW® PLUS (Mouse-HRP/Rabbit-AP) IHC Kit (Brown/Green)	150 tests	ENZ-KIT181
POLYVIEW® PLUS HRP (anti-Rabbit) Reagent	150 tests	ENZ-ACC103
POLYVIEW® PLUS HRP (anti-Mouse) Reagent	150 tests	ENZ-ACC104
POLYVIEW® PLUS AP (anti-Rabbit) Reagent	150 tests	ENZ-ACC110
HIGHDEF® Blue IHC chromogen (AP)	30 mL	ADI-950-150
HIGHDEF® DAB Chromogen/Substrate Set	200 mL	ENZ-ACC105
HIGHDEF® Hematoxylin	100 mL	ENZ-ACC106
Antigen Retrieval Reagent, pH 6 (10X)	100 mL	ENZ-ACC112

Advance Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com