WESTERNVIEW® Dual Detection Kit (AntiMouse/Anti-Rabbit)

Chromogenic Western Blot kit with increased signal and sensitivity

The WESTERNVIEW[®] Dual Detection Kit (Anti-Mouse/Anti-Rabbit) can be used to identify primary antibodies on a Western Blot using immunodetection. The kit contains a complete reagent set, with optimized ready-to-use or ready-to-dilute reagents.

Western Blotting is a common lab technique used to detect proteins in a sample of tissue homogenate or extract. Due to its flexibility, it can be used for a wide range of applications within molecular biology and biochemistry, including detection of post-translational modifications and verification of protein cloning. Using the size and color intensity of the protein band, a semi-quantitative estimation of protein can be derived.

Ordering Information

Order Online »

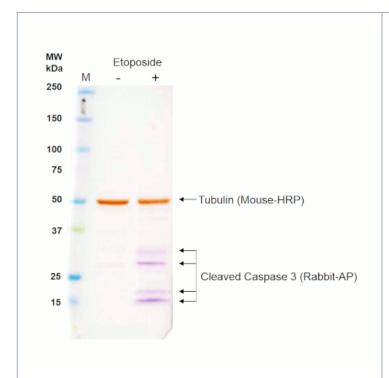
ENZ-KIT163-0001

1Kit

Manuals, SDS & CofA

View Online »

- More sensitive detection able to detect lower concentrations of protein
- Easy to use all reagents are ready-to-use or ready-to-dilute, saving precious time
- Amenable to multiplex analysis determine source of protein bands with differentiated colors
- Easily adaptable works with a wide range of mouse-derived proteins



PANC-1 cells treated with etoposide. The band at 50 kDa is clearly revealed to be tubulin, while the smaller bands are known to be cleaved caspase.

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Assay Time 2 hours

Contents

- Antibody Diluent, 4 x 90 mL
 Diluent solution for dilution of primary and secondary antibodies.
- NBT/BCIP, Ready-to-Use, 100 mL
 Ready-made solution for chromogenic blotting to be used when developing anti-rabbit secondary.
- DAB Chromogen (25X), 4 mL
 Concentrated chromogen for chromogenic blotting to be used when developing anti-mouse secondary.
- DAB Substrate Buffer, 100 mL Buffer used to dilute DAB Chromogen (25X).
- SignaSure[®] Buffer Salts, 2 packets
 Contents used to make SignaSure[®] Wash Buffer.
- Tween[®]-20, 2 mL Detergent for SignaSure[®] Wash Buffer.
- Secondary Antibody (HRP Anti-Mouse), 2 mL
 Anti-Mouse secondary antibody for detection of mouse-based primary antibodies in Western Blots.
 Use DAB Substrate Buffer and Chromogen to develop.
- Secondary Antibody (AP Anti-Rabbit), 2 mL
 Anti-Rabbit secondary antibody for detection of rabbit-based primary antibodies in Western Blots.
 Use NBT/BCIP solution to develop.

Species Reactivity



Last modified: May 29, 2024

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Mouse, Rabbit

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com