

# WESTERNVIEW<sup>®</sup> Dual Detection Kit (Anti- Mouse/Anti-Rabbit)

## Chromogenic Western Blot kit with increased signal and sensitivity

The WESTERNVIEW<sup>®</sup> 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

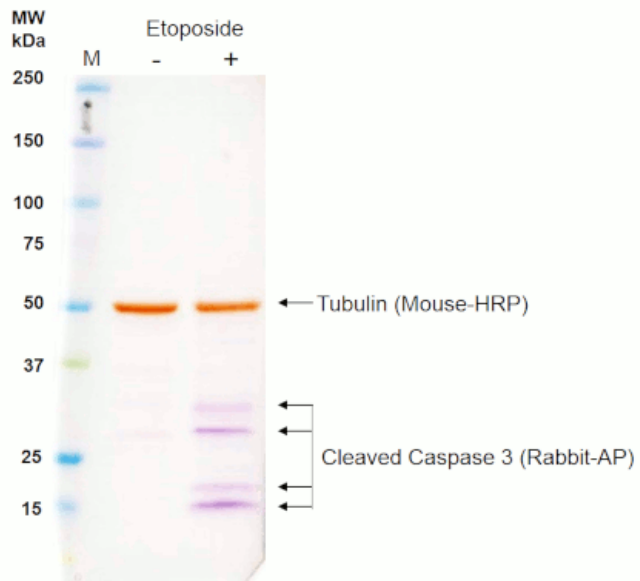
[Order Online »](#)

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

1. Antibody Diluent, 4 x 90 mL  
Diluent solution for dilution of primary and secondary antibodies.
2. NBT/BCIP, Ready-to-Use, 100 mL  
Ready-made solution for chromogenic blotting to be used when developing anti-rabbit secondary.
3. DAB Chromogen (25X), 4 mL  
Concentrated chromogen for chromogenic blotting to be used when developing anti-mouse secondary.
4. DAB Substrate Buffer, 100 mL  
Buffer used to dilute DAB Chromogen (25X).
5. SignaSure<sup>®</sup> Buffer Salts, 2 packets  
Contents used to make SignaSure<sup>®</sup> Wash Buffer.
6. Tween<sup>®</sup>-20, 2 mL  
Detergent for SignaSure<sup>®</sup> Wash Buffer.
7. 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.
8. 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 Mouse, Rabbit



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-  
fr@enzolifesciences.com](mailto:info-fr@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)