WESTERNVIEW[®] Detection Kit (Anti-Mouse)

Chromogenic Western Blot kit with increased signal and sensitivity

The WESTERNVIEW[®] Detection Kit (Anti-Mouse) 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.

Citations: 1

View Online »

Ordering Information

Order Online »

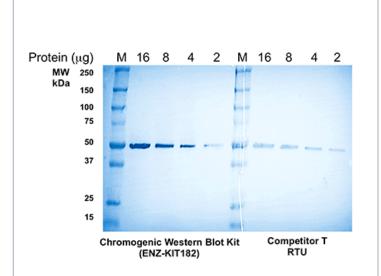
ENZ-KIT182-0001

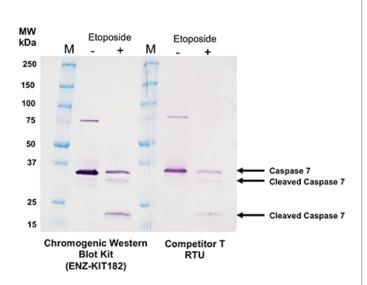
1Kit

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

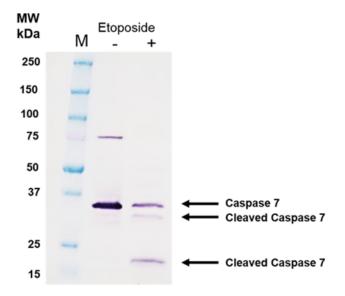




Linear dilution of protein (μg) run in the Chromogenic Western Blot Kit (ENZ-KIT182) and the competitor's kit.

Caspase-7, cleaved and uncleaved, run on the WESTERNVIEW™ Detection Kit (Anti-Mouse) (ENZ-KIT182) and a competitor's kit.

Jurkat cells treated with or without 12.5 µM Etoposide for 18 hours 12 µg total lysate loaded



Western Blot of Jurkat cells using WESTERNVIEW™ Detection Kit

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, WB

Assay Time 2 hours

Contents

- Antibody Diluent, 4 x 90 mL
 Diluent solution for dilution of primary and secondary antibodies.
- NBT/BCIP, 100 mL Ready-to-use solution for chromogenic blotting.
- 3. Wash Buffer (10X), 2 x 100 mL10X Wash Buffer for washing membrane.
- Secondary Antibody (AP Anti-Mouse), 4.5 mL
 Anti-Mouse secondary antibody conjugated to
 alkaline phosphatase for detection of mouse-based primary antibodies in Western Blot.

Species Reactivity Mouse

Last modified: May 29, 2024

