

Vector® TrueVIEW™ Autofluorescence Quenching Kit

Cat. No.: SP-8400

Storage: 2-8 °C

DESCRIPTION

Vector® TrueVIEW™ Autofluorescence Quenching Kit provides a novel way to remove unwanted fluorescence in tissue sections due to aldehyde fixation, red-blood cells, and structural elements such as collagen and elastin. This unique formulation binds and effectively quenches the autofluorescent elements in even the most problematic tissues, such as kidney, spleen and pancreas. The use of Vector® TrueVIEW™ leads to significant enhancement in overall signal-to-noise in most immunofluorescence assays.

KIT COMPONENTS

Product Name	Volume
Vector® TrueVIEW™ Reagent A	5 ml
Vector® TrueVIEW™ Reagent B	5 ml
Vector® TrueVIEW™ Reagent C	5 ml
VECTASHIELD® HardSet™ Antifade Mounting Medium	2 ml

STORAGE:

- Store reagents in original bottles at 2-8 °C
- Avoid storing reagents or working solution in strong direct light

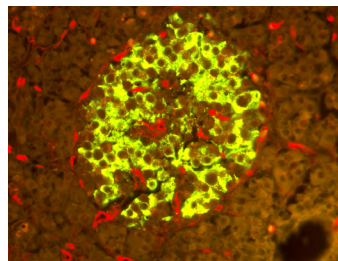
SIGNAL TO NOISE OPTIMIZATION:

The primary antibody and detection reagents should be optimized (titered) in conjunction with Vector® TrueVIEW™ Quenching Kit to achieve maximum signal to noise. With some sections, it may be useful to reduce the incubation time of the TrueVIEW™ Reagent to optimize signal to noise.

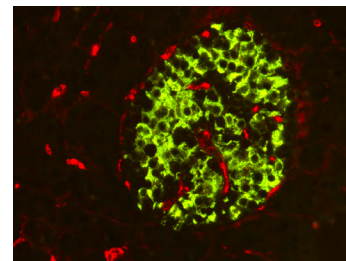
Related Reagents

Product Name	Unit Size	Cat. No.
VECTASHIELD® HardSet™ Antifade Mounting Medium	10 ml	H-1400
VECTASHIELD® HardSet™ Antifade Mounting Medium with DAPI	10 ml	H-1500

Without Treatment



With Vector® TrueVIEW™ Treatment



Human Pancreas (FFPE): Stained for CD34 (using anti-mouse DyLight® 594, red) and Insulin (using anti-guinea pig, fluorescein, green). Coverslipped with VECTASHIELD® HardSet™ Antifade Mounting Medium. Note significant reduction of autofluorescence in treated section (right) with the retention of specific staining.

INSTRUCTIONS FOR USE:

A) Reagent Preparation

For each standard tissue section in your assay, you will need approximately 150 µL of Vector® TrueVIEW™ Reagent (i.e., 50 µL A + 50 µL B + 50 µL C).

To prepare Vector® TrueVIEW™ Reagent, a ratio of 1:1:1 of Reagents A, B and C is required. The order of mixing is important.

- 1) Add equal volumes of Reagent A and Reagent B in a clean test tube. Mix for 10 seconds.
- 2) Add Reagent C to the mixture (ensuring a 1:1:1 volume ratio) and mix again for 10 seconds.

Vector® TrueVIEW™ Reagent is now ready to use. Once prepared, Vector® TrueVIEW™ Reagent is stable for at least 2 hours at room temperature.

B) Tissue Treatment Procedure

Following completion of the immunofluorescent staining:

- 1) Drain excess buffer from tissue section.
- 2) Add Vector® TrueVIEW™ Reagent to cover tissue section completely (~150 µl); and incubate for 2 - 5 minutes.
- 3) Wash in PBS buffer for 5 minutes.
- 4) Drain excess buffer from section.
- 5) Dispense VECTASHIELD® HardSet™ Antifade Mounting Medium onto the tissue section. Coverslip and allow VECTASHIELD® HardSet™ Antifade Mounting Medium to disperse over the entire section.
- 6) Slides can be visualized immediately after mounting, but the coverslip will not be immobilized until mounting media is cured at room temperature for 1-2 hours.
- 7) For optimal results, slides should be evaluated within 48 hours of mounting.

Patent Pending Formulation

For Laboratory Use Only