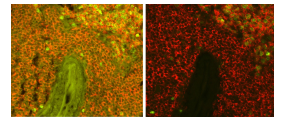
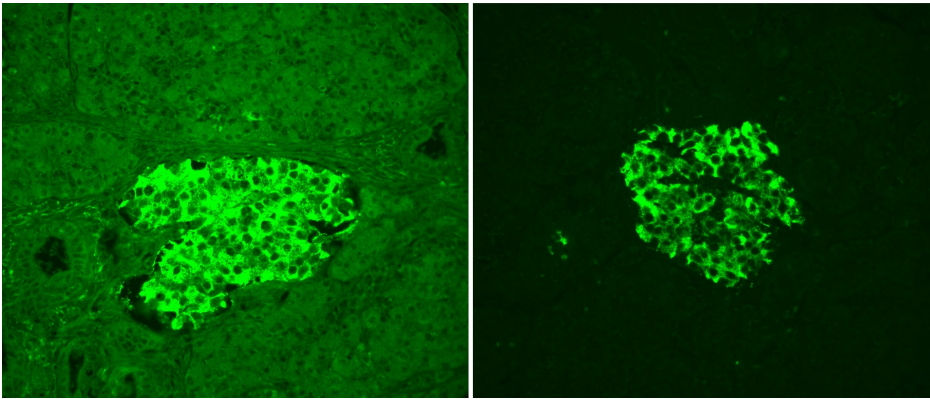




# Vector® TrueVIEW® Autofluorescence Quenching Kit

## SP-8400-15

[Product Images](#)



## Short Description

---

The TrueVIEW Autofluorescence Quenching Kit provides a novel way to diminish unwanted autofluorescence from non-lipofuscin sources and dramatically improve signal-to-noise ratio. TrueVIEW removes unwanted autofluorescence in tissue sections due to aldehyde fixation, the presence of red blood cells, and structural elements such as collagen and elastin. The quenching action of the kit reagents provides a clear, unambiguous, “true view” localization of the target antigen.

### This kit offers you:

- Effectivity with problematic tissues such as kidney and spleen
- An easy-to-use, one-step method
- Compatibility with a wide selection of fluorophores
- Compatibility with standard epifluorescence and confocal laser microscopes
- Availability with DAPI counterstain and without.

One kit is sufficient to treat approximately 100 to 150 tissue sections.

## Additional Information

---

Unit Size	15 ml
Applications	Immunofluorescence, In situ hybridization
Blocking Action	Autofluorescence
Format	Concentrate

