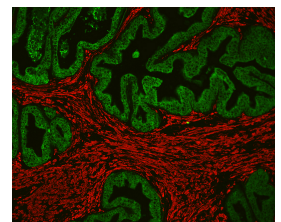
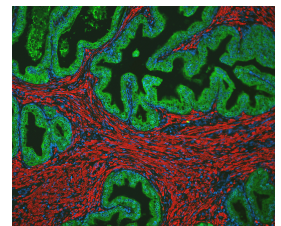
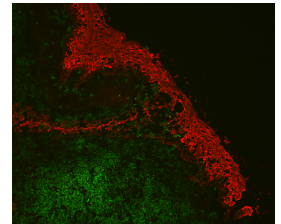
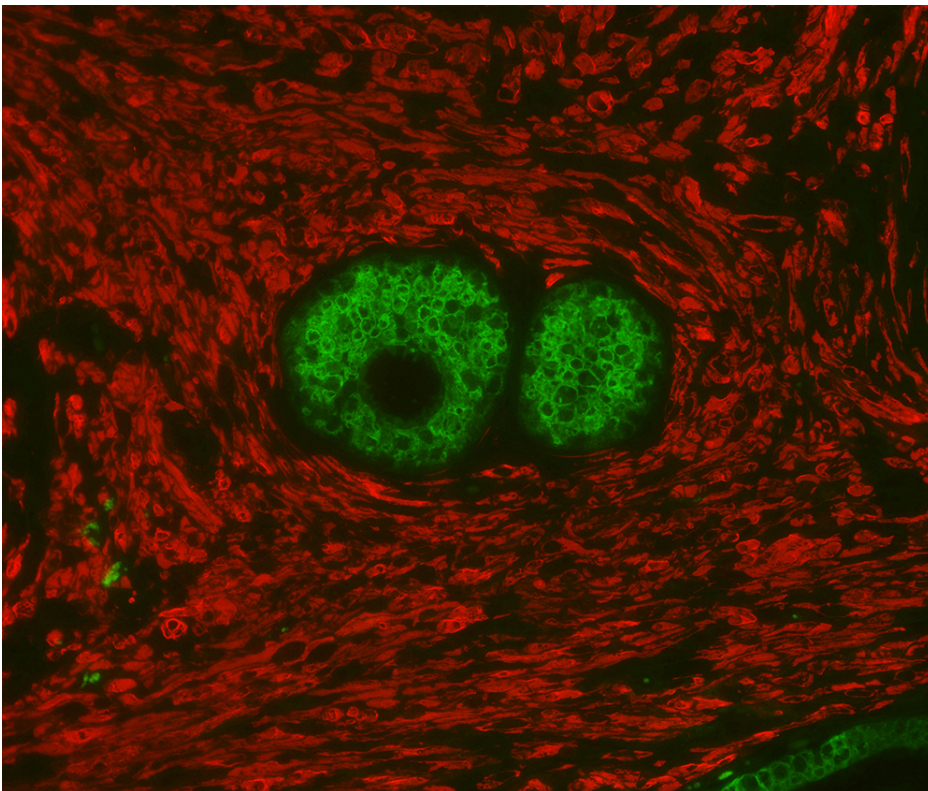




# VectaFluor™ Duet Immunofluorescence Double Labeling Kit, DyLight™ 488 Anti- Rabbit (green), DyLight™ 594 Anti- Mouse (red)

## DK-8818

[Product Images](#)





## Short Description

---

VectaFluor Duet Kit offers maximum convenience to achieve immunofluorescence double labeling in one step. This kit includes a ready-to-use detection reagent that will detect mouse and rabbit primary antibodies with green and red DyLight fluorescent dyes in a single step. Our affinity purified, extensively cross-adsorbed secondary antibodies are conjugated to DyLight dyes in a manner that ensures maximum degree of labeling without compromising antibody affinity or specificity. VectaFluor Duet is compatible with fluorescence staining of cells and tissues.

### Features:

- Robust cocktail formulation of DyLight anti-mouse IgG/anti-rabbit IgG secondary antibodies
- Affinity-purified, ultrapure, high affinity antibodies
- Secondary detection reagent optimally labeled with DyLight dyes to provide the brightest label for fluorescence microscopy

## Additional Information

---

Unit Size	1 kit
Applications	Immunofluorescence
Target Species	Universal (Mouse/Rabbit)
Conjugate	DyLight 488/DyLight 594
Color of Fluorescence	Kit (Green, Red)
Host Species	Horse
Format	Ready-to-Use

