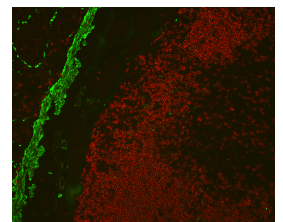
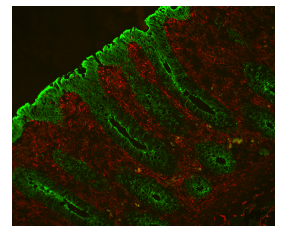
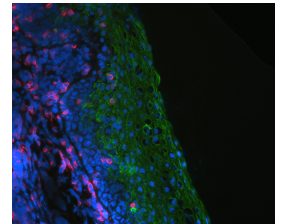
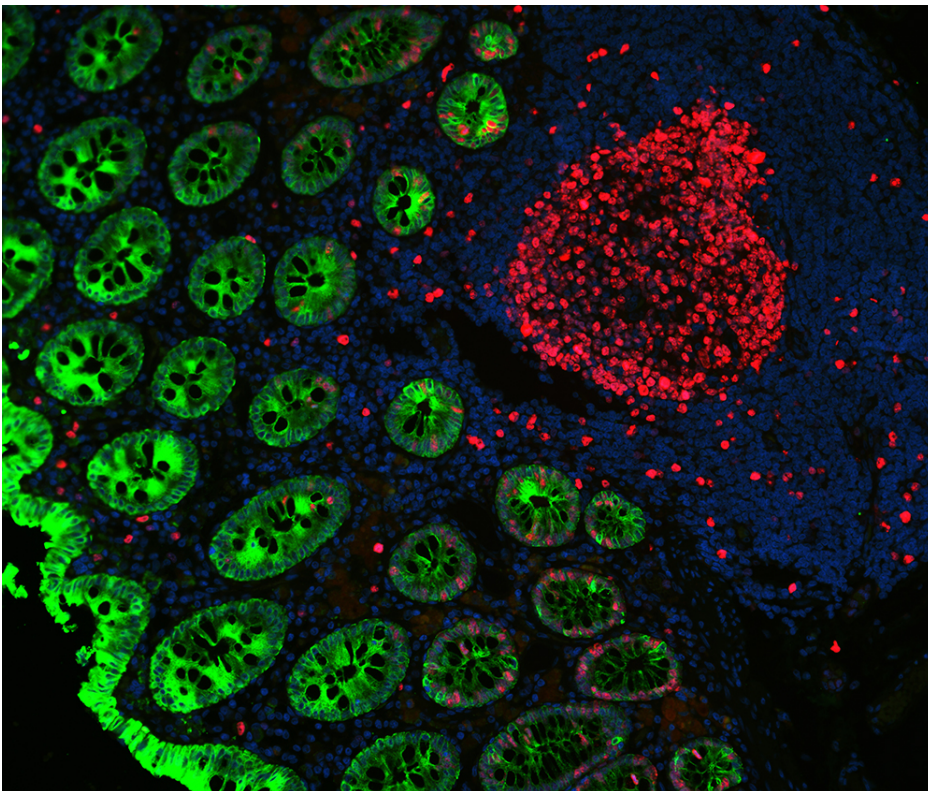


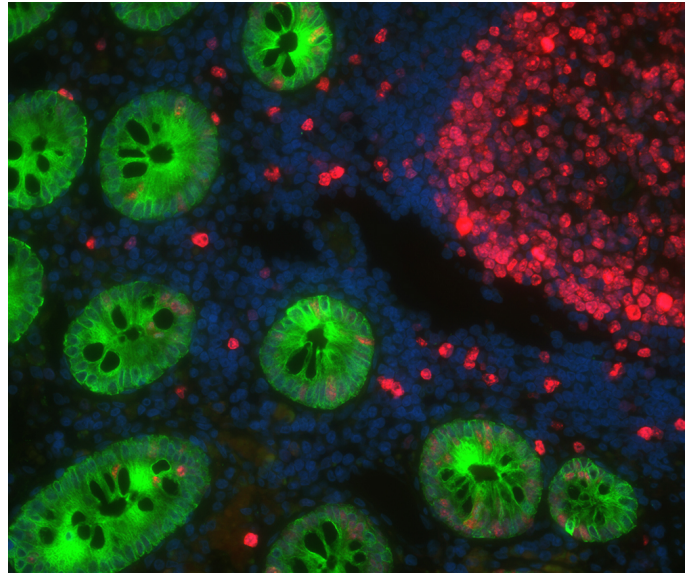
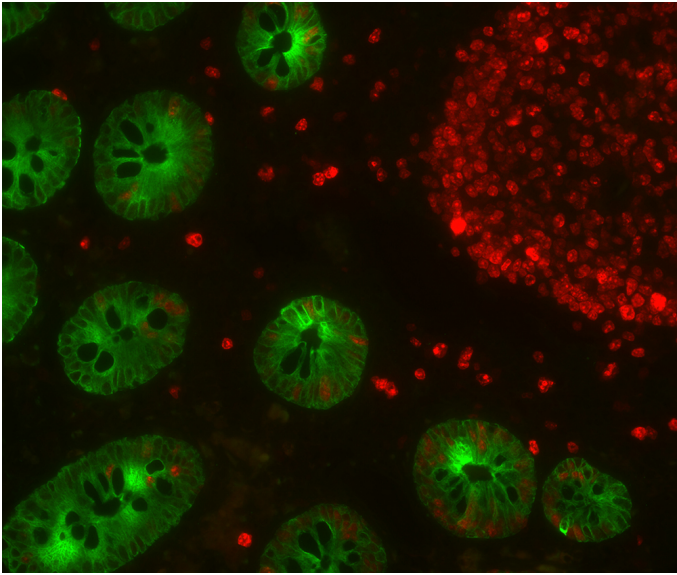


VectaFluor™ Duet Immunofluorescence Double Labeling Kit, DyLight™ 594 Anti- Rabbit (red), DyLight™ 488 Anti- Mouse (green)

DK-8828

[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

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 |

