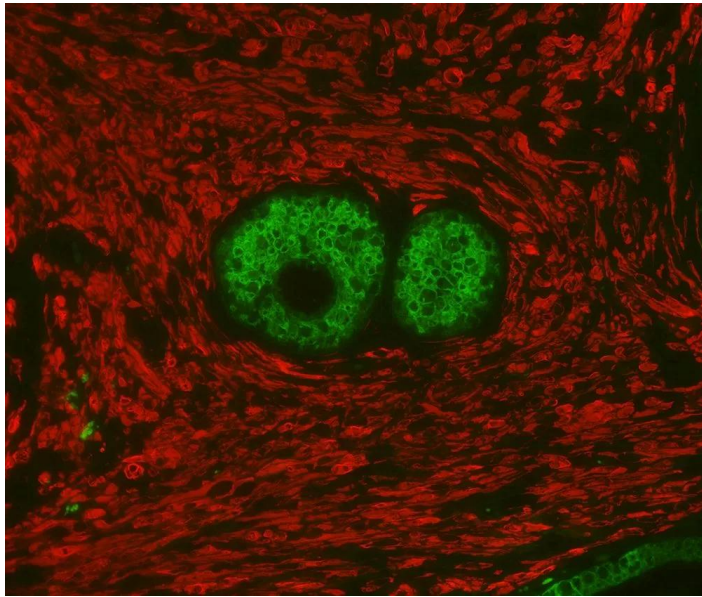




VECTAFLUOR DUET IMMUNOFLOURESCENCE DOUBLE LABELING KIT, DYLIGHT 488 ANTI-RABBIT (GREEN), DYLIGHT 594 ANTI-MOUSE (RED)

SKU: DK-8818



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

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



- Affinity-purified, ultrapure, high affinity antibodies
- Secondary detection reagent optimally labeled with DyLight dyes to provide the brightest label for fluorescence microscopy

SPECIFICATIONS

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

KIT COMPONENTS

- 15 ml 2.5% Normal Horse Serum (ready-to-use) for blocking
- 15 ml VectaFluor Duet Reagent [DyLight 488 Anti-Rabbit IgG and DyLight 594 Anti-Mouse IgG cocktail (rabbit-green, mouse-red)] made in horse, ready-to-use in 5% NHS.

DyLight® 488:

- Excitation maximum: 493 nm
- Emission maximum: 518 nm
- Color: Green

DyLight® 594:

- Excitation maximum: 592 nm
- Emission maximum: 617 nm
- Color: Red

TECHNICAL INFORMATION

Kit Contents:

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



- 15 ml 2.5% Normal Horse Serum (ready-to-use) for blocking
- 15 ml VectaFluor Duet Reagent [DyLight 488 Anti-Rabbit IgG and DyLight 594 Anti-Mouse IgG cocktail (rabbit-green, mouse-red)] made in horse, ready-to-use in 5% NHS.

DyLight 488:

- Excitation maximum: 493 nm
- Emission maximum: 518 nm
- Color: Green

DyLight 594:

- Excitation maximum: 592 nm
- Emission maximum: 617 nm
- Color: Red

DyLight fluorescent dyes are direct alternatives to traditional fluorophores such as fluorescein (FITC) and rhodamine. The excitation and emission spectra parallel that of other commercially available fluorescent reagents allowing for easy substitution into an existing protocol without requiring any further instrumentation or filter sets.

DyLight dyes offer a number of potential advantages including greater photostability and brighter fluorescence. DyLight dyes are completely stable over a pH range of 4-9 making them compatible with many aqueous-based buffers and diluents. The DyLight dyes can be applied as single labels or in combination with other DyLight dyes and fluorophores as part of a multiple immunofluorescent antigen staining methodology. The DyLight dyes currently offered are DyLight 488 (green), DyLight 549 (orange), DyLight 594 (red), and DyLight 649 (far red).

DOCUMENTS

- [Safety Data Sheet](#)
- [User Guide](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES

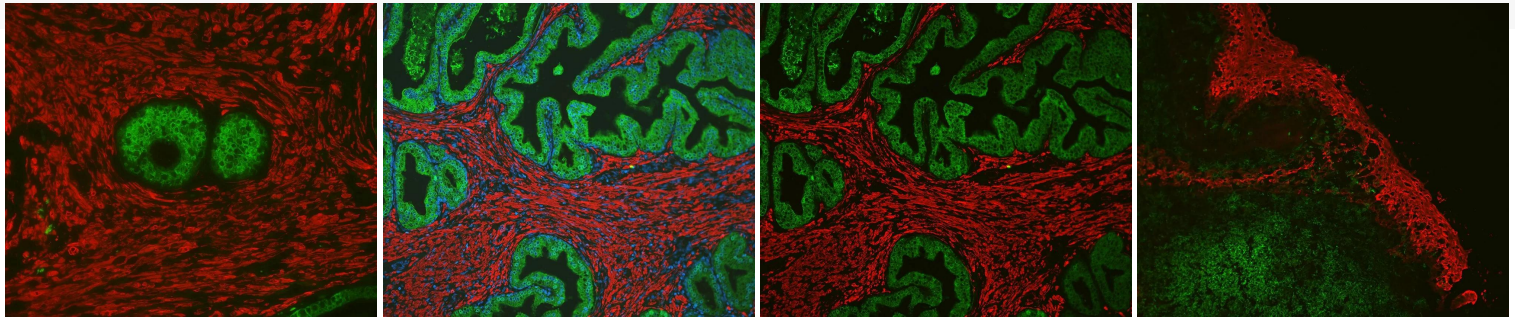
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



www.vectorlabs.com

Email: customerservice@vectorlabs.com

Telephone: [\(650\) 697-3600](tel:(650)697-3600)



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

[VectaFluor Duet Immunofluorescence Double Labeling Kit, DyLight 488 Anti-Rabbit \(green\), DyLight 594 Anti-Mouse \(red\)](https://vectorlabs.com/products/vectafluor-duet-kit-rabbit-gr-mouse-red/)

<https://vectorlabs.com/products/vectafluor-duet-kit-rabbit-gr-mouse-red/>