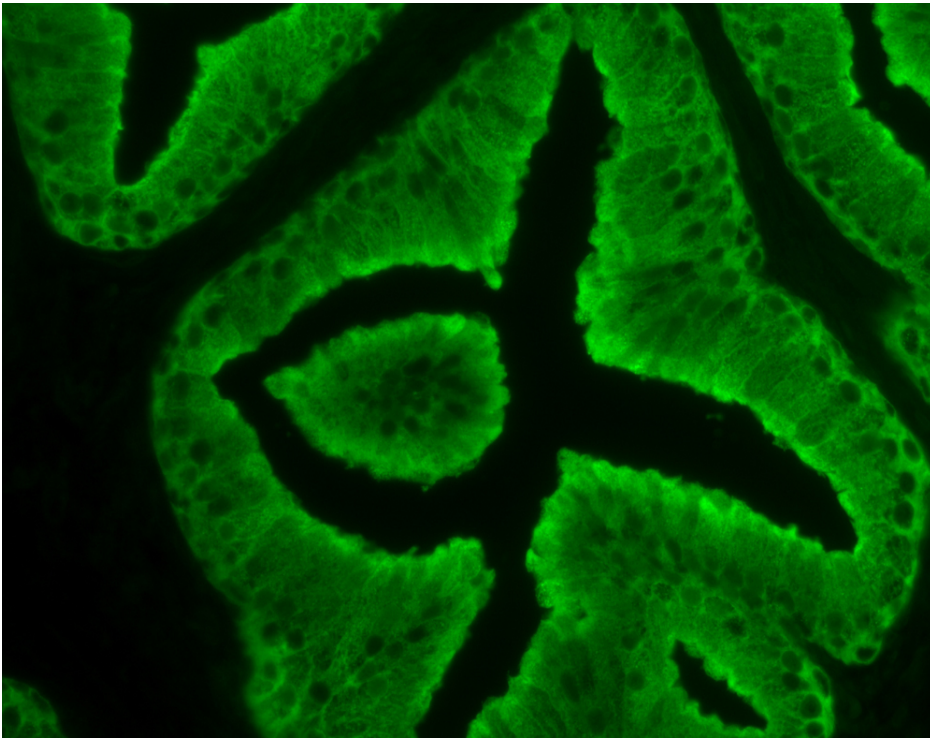




# VectaFluor™ Horse Anti-Rabbit IgG, DyLight™ 488 Antibody Kit, R.T.U.

## DI-1788

[Product Images](#)



## Short Description

---

VectaFluor R.T.U. DyLight 488 Anti-Rabbit IgG can be used for immunofluorescence and other applications. Optimal F/P ratios have been established for each conjugate to ensure maximum fluorescence with minimal background staining.

### Features:

- Ready-to-use format offers maximum convenience for fluorescence staining of cells and tissues
- Recognizes both heavy and light chains (H+L)
- Optimally labeled with DyLight 488 to provide the brightest label for fluorescence microscopy
- Excitation: 493 nm
- Emission: 518 nm
- Color: Green

### Kit Contents:

- 15 ml pre-diluted, stabilized secondary antibody solution
- 15 ml 2.5% normal horse serum for blocking

## Additional Information

---

Unit Size	1 kit
Applications	Immunofluorescence, In situ hybridization, Flow Cytometry/Cell Separation
Maximum Emission	518 nm
Maximum Excitation	493 nm
Target Species	Rabbit
Conjugate	DyLight 488
Color of Fluorescence	Green
Host Species	Horse
Format	Ready-to-Use

