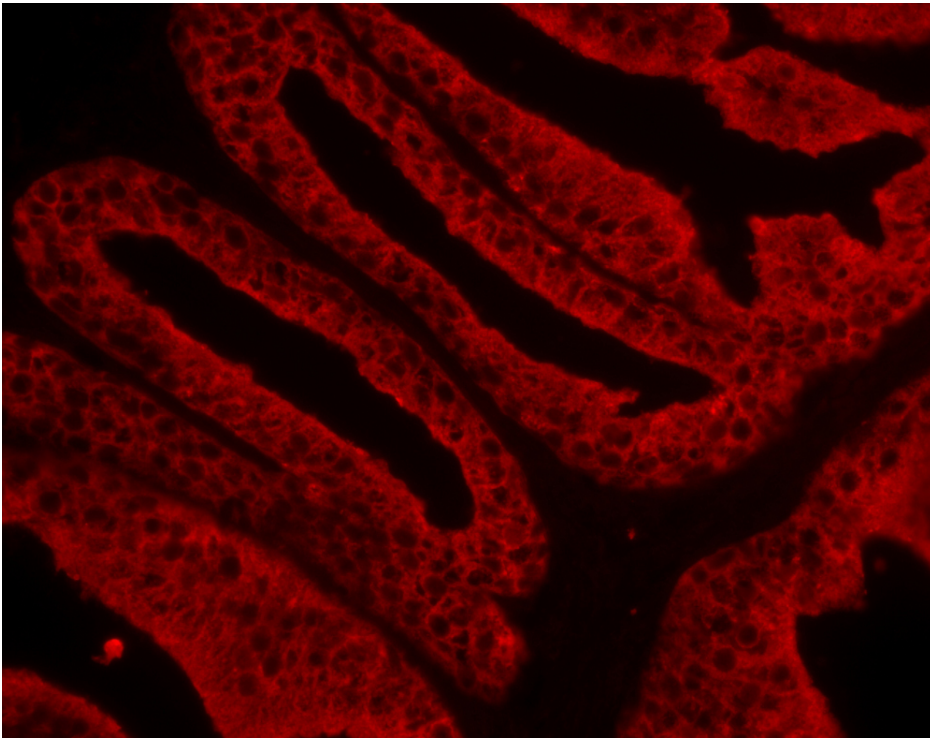




VectaFluor™ Horse Anti-Rabbit IgG, DyLight™ 594 Antibody Kit, R.T.U.

DI-1794

[Product Images](#)



Short Description

VectaFluor R.T.U. DyLight 594 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 594 to provide the brightest label for fluorescence microscopy
- Excitation: 592 nm
- Emission: 617 nm
- Color: Red

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	617 nm
Maximum Excitation	592 nm
Target Species	Rabbit
Conjugate	DyLight 594
Color of Fluorescence	Red
Host Species	Horse
Format	Ready-to-Use

