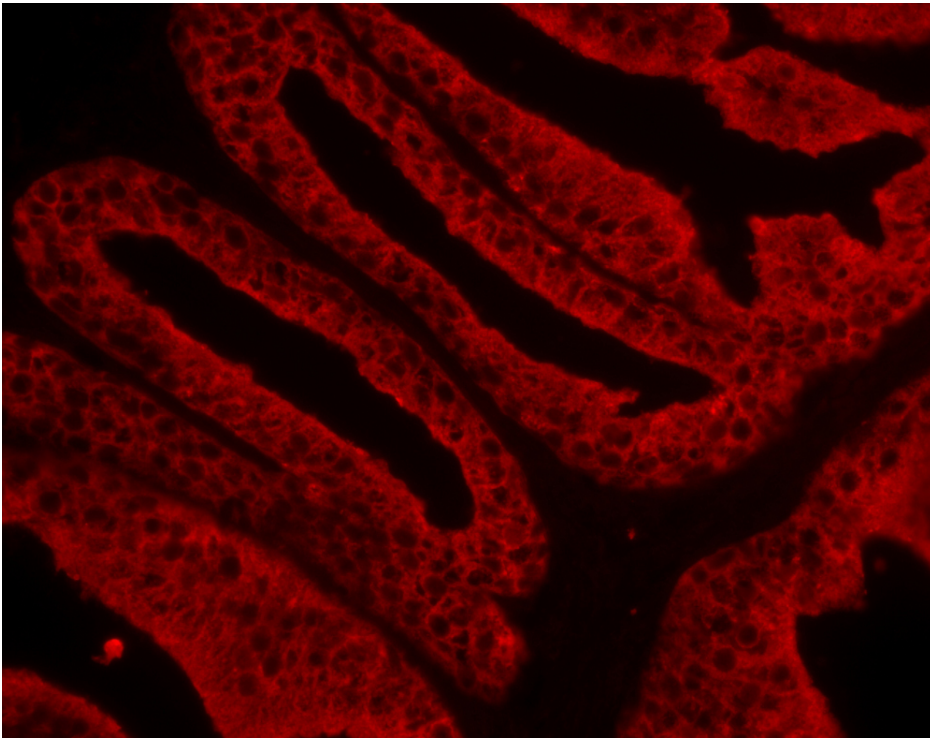




VectaFluor™ Horse Anti-Mouse IgG, DyLight™ 594 Antibody Kit, R.T.U.

DI-2794

[Product Images](#)



Short Description

VectaFluor Ready-to-Use (R.T.U.) DyLight dye-labeled Secondary Antibodies offer maximum convenience for fluorescence staining of cells and tissues. 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. DyLight dyes offer several advantages such as brighter fluorescence, greater photostability, and pH independence.

Features:

- 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	Mouse
Conjugate	DyLight 594
Color of Fluorescence	Red
Host Species	Horse
Format	Ready-to-Use

