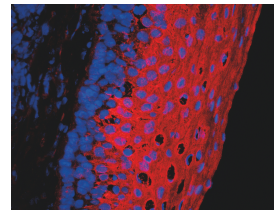




VectaFluor™ Excel Amplified Anti-Mouse IgG, DyLight™ 594 Antibody Kit

DK-2594

[Product Images](#)



Short Description

VectaFluor Excel Amplified Fluorescent Staining System offers a non-biotin amplification method for fluorescence applications. This system uses a ready-to-use (R.T.U.) Amplifier Antibody, followed by a ready-to-use (R.T.U.) VectaFluor DyLight dye-labeled detection system.

Features:

- Ready-to-use
- High sensitivity
- Low background
- VectaFluor DyLight 594 labeled anti-goat IgG antibody for detection
- Excitation: 592 nm
- Emission: 617 nm
- Color: Red

Additional Information

Unit Size	1 kit
Applications	Immunofluorescence
Technology	Amplified Fluorescence Systems
Conjugate	DyLight 594
Color of Fluorescence	Red
Target Species	Mouse
Format	Ready-to-Use

