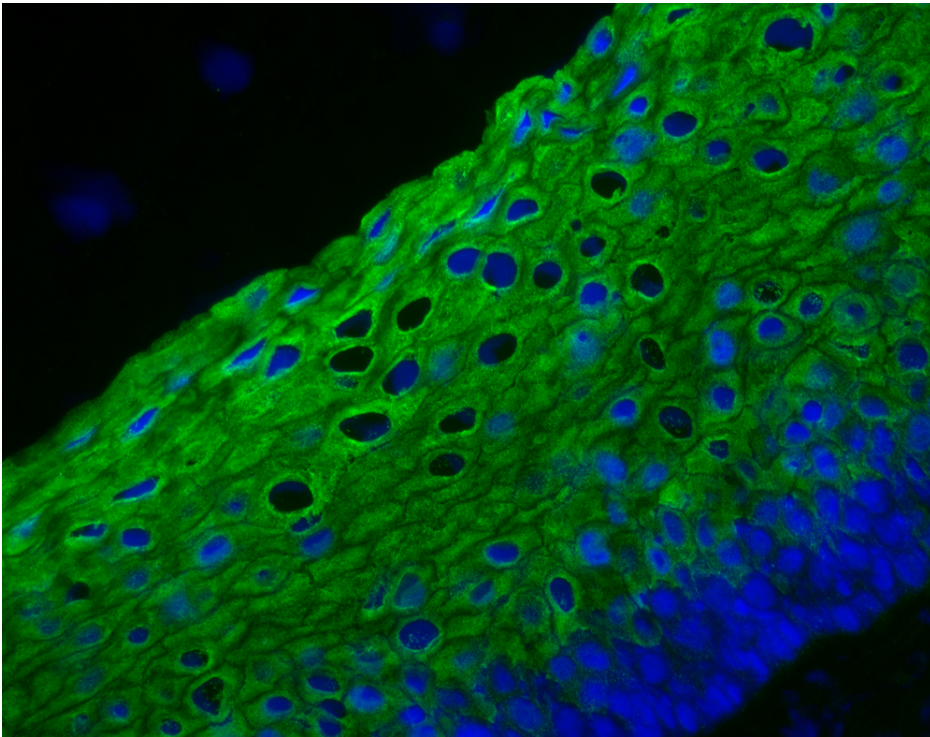




# VectaFluor™ Excel Amplified Anti-Mouse IgG, DyLight™ 488 Antibody Kit

## DK-2488

[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:

- High sensitivity
- Low background
- VectaFluor DyLight 488 labeled anti-goat IgG antibody for detection
- Excitation: 493 nm
- Emission: 518 nm
- Color: Green

## Additional Information

---

Unit Size	1 kit
Applications	Immunofluorescence
Technology	Amplified Fluorescence Systems
Conjugate	DyLight 488
Color of Fluorescence	Green
Target Species	Mouse
Format	Ready-to-Use

