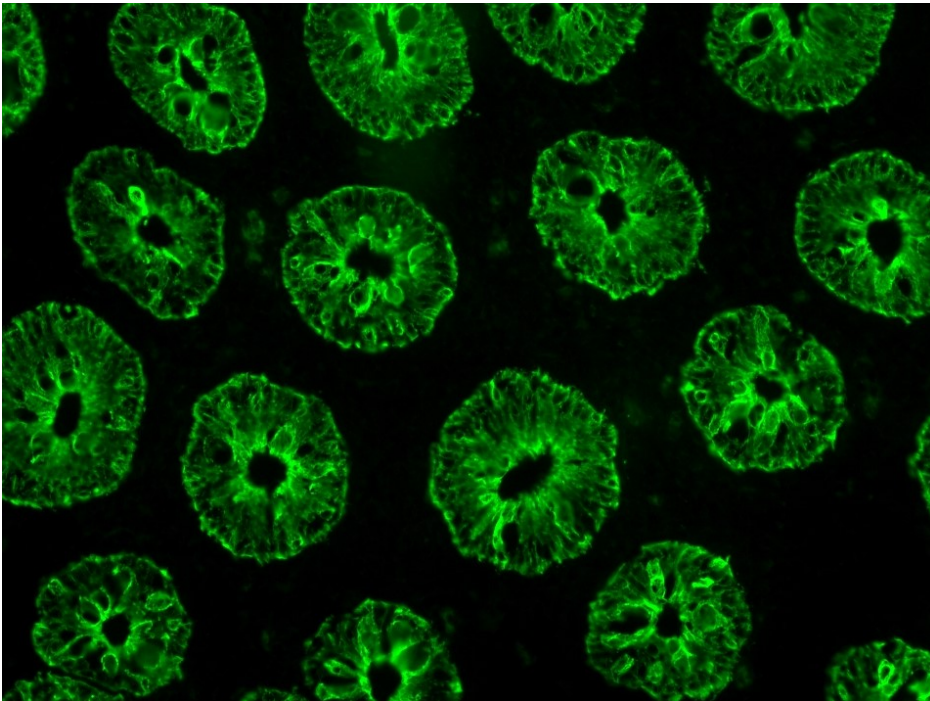




VectaFluor™ Excel Amplified Anti-Rabbit IgG, DyLight™ 488 Antibody Kit

DK-1488

[Product Images](#)



Short Description

VectaFluor Excel Amplified DyLight™ 488 Anti-Rabbit IgG 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 | Rabbit |
| Format | Ready-to-Use |

