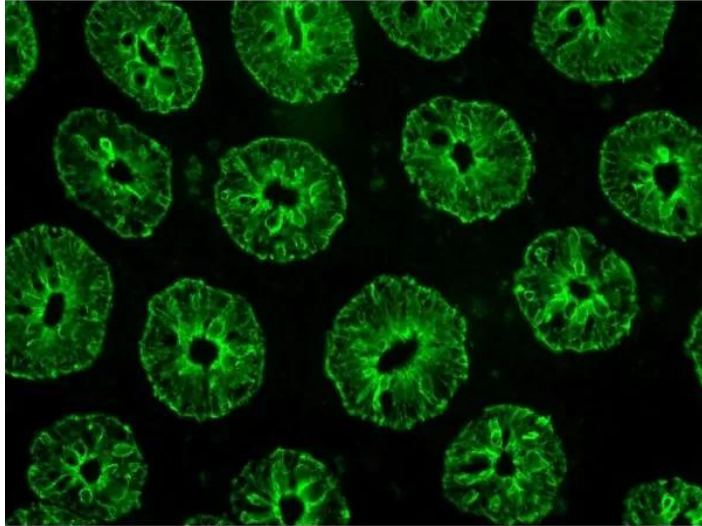




VECTAFLUOR EXCEL AMPLIFIED ANTI-RABBIT IGG, DYLIGHT 488 ANTIBODY KIT

SKU: DK-1488



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

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



SPECIFICATIONS

Color of Fluorescence	Green
Format	Ready-to-Use
Isolation Technology	Amplified Fluorescence Systems
Unit Size	1 Kit
Applications	Immunofluorescence
Target Species	Rabbit
Conjugate	DyLight 488
Platform	Amplified Fluorescence Systems

KIT COMPONENTS

- 15 ml 2.5% Normal Horse Serum for blocking
- 15 ml Amplifier Antibody (Goat anti-rabbit IgG)
- 15 ml VectaFluor DyLight labeled Horse Anti-Goat IgG

TECHNICAL INFORMATION

All the reagents in the VectaFluor Excel Kit are affinity purified and extensively cross-adsorbed to ensure high sensitivity and low background. In addition, the VectaFluor™ Reagents are conjugated to the DyLight dyes using technology that ensures maximum labeling without compromising antibody affinity or specificity. DyLight dyes offer several advantages such as brighter fluorescence, greater photostability, and pH independence.

All reagents are supplied in stabilized, R.T.U. solutions for maximum convenience.

Kit Contents:

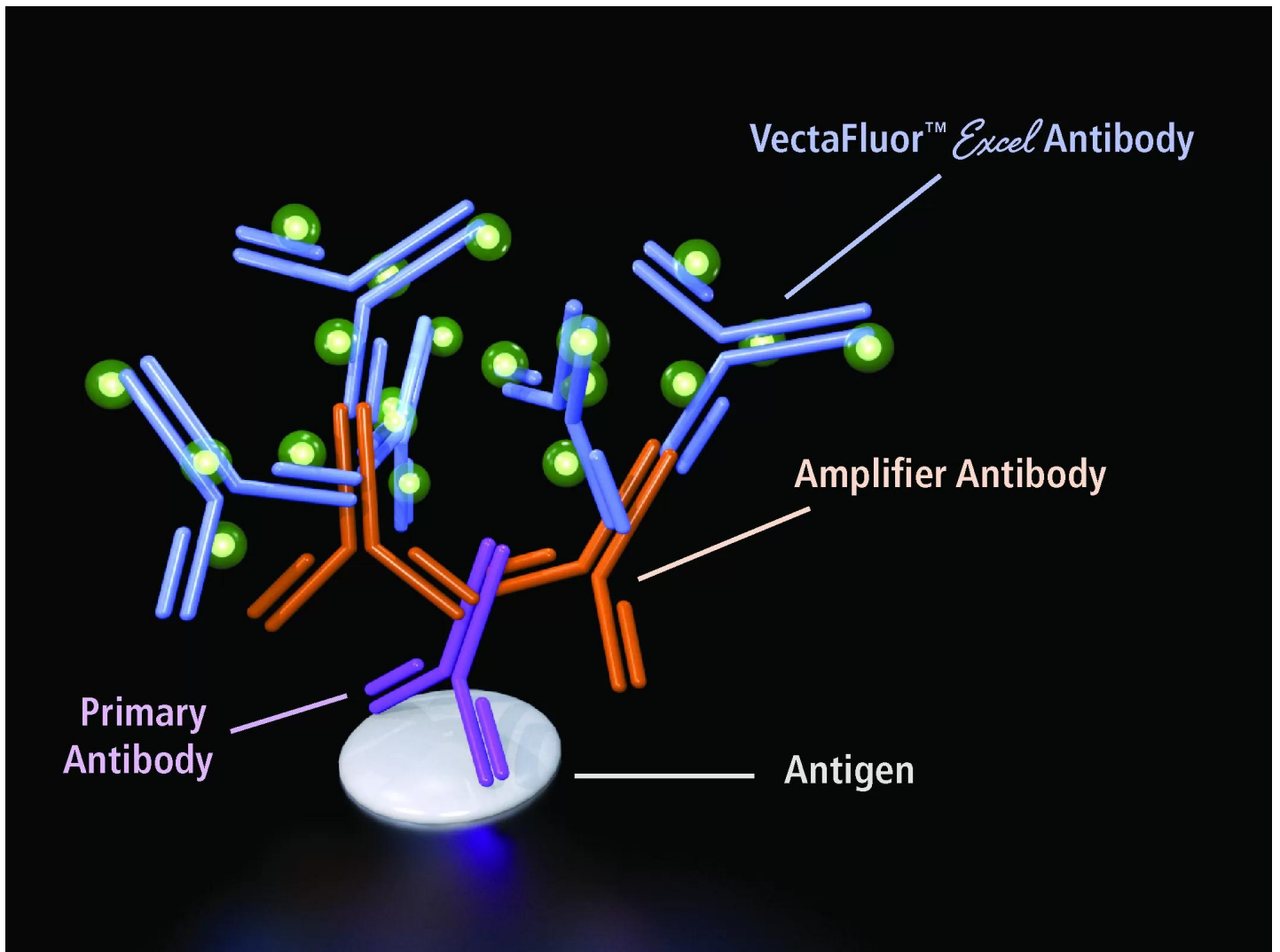
- 15 ml 2.5% Normal Horse Serum (ready-to-use) for blocking
- 15 ml Amplifier Antibody (Goat anti-rabbit IgG; ready-to-use)
- 15 ml VectaFluor™ DyLight™ labeled Horse Anti-Goat IgG (ready-to-use)

- Excitation maximum: 493 nm
- Emission maximum: 518 nm

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



- Color: Green



CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

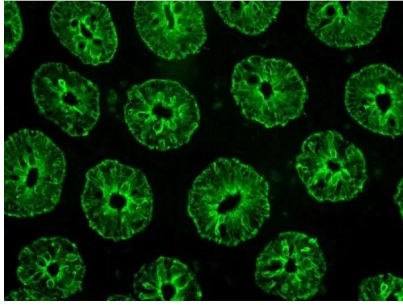
- [User Guide](#)

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.