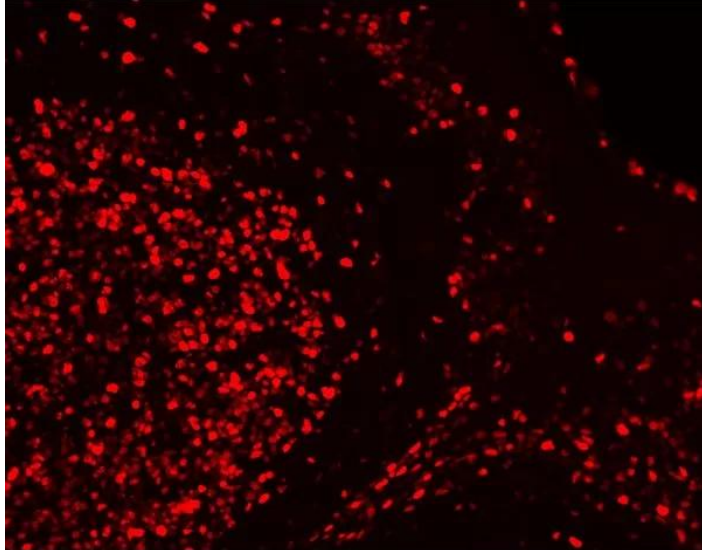




VECTAFLUOR EXCEL AMPLIFIED ANTI-RABBIT IGG, DYLIGHT 594 ANTIBODY KIT

SKU: DK-1594



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 594 labeled anti-goat IgG antibody for detection
- Excitation: 592 nm
- Emission: 617 nm
- Color: Red

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



SPECIFICATIONS

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

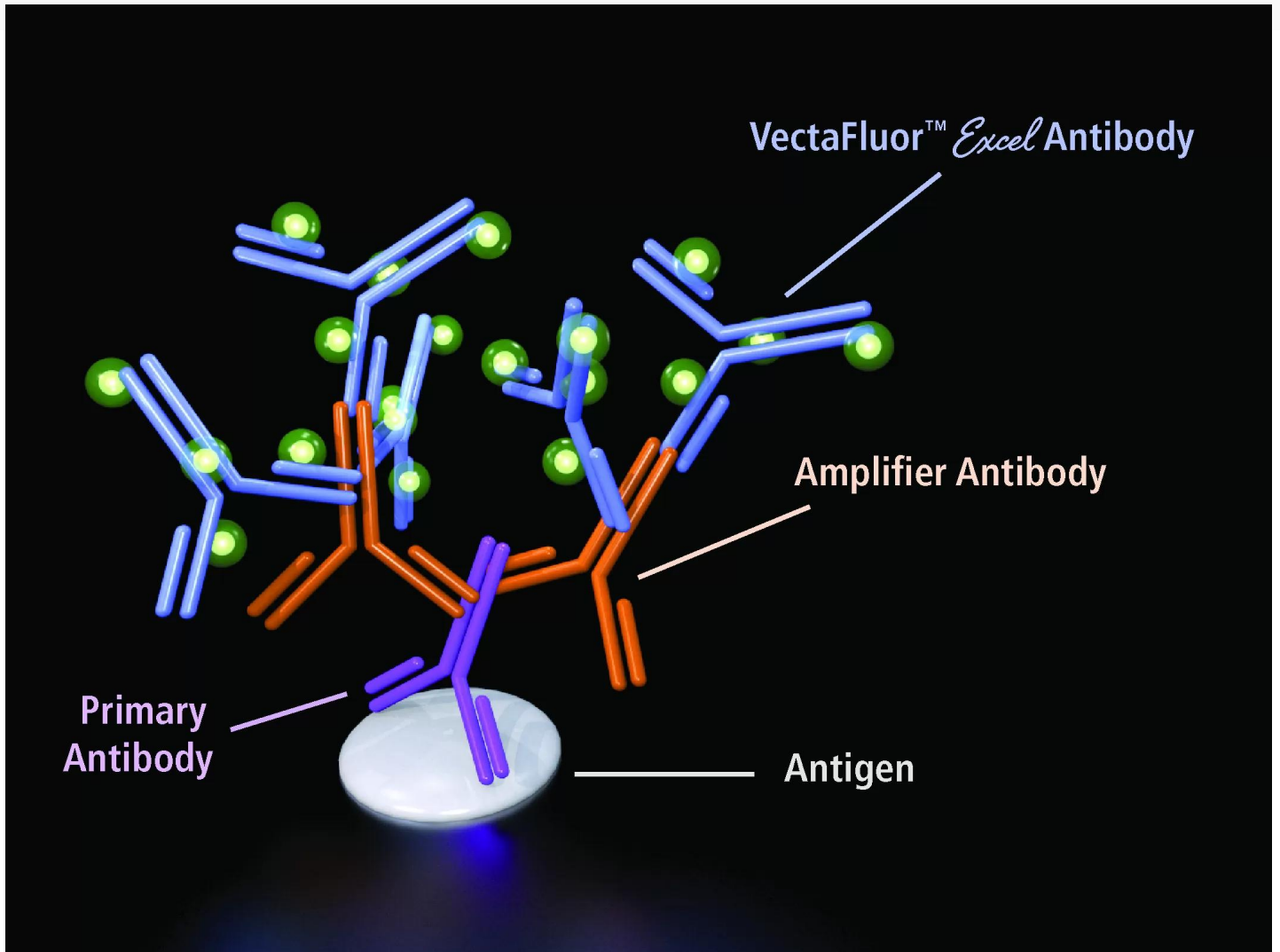
All the reagents in the VectaFluor Excel Kit are affinity purified and extensively cross-adsorbed to ensure high sensitivity and low background. In addition, the VectaFluor Reagents are conjugated to the DyLight dyes using technology that ensures maximum labeling without compromising antibody affinity 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)

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



CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

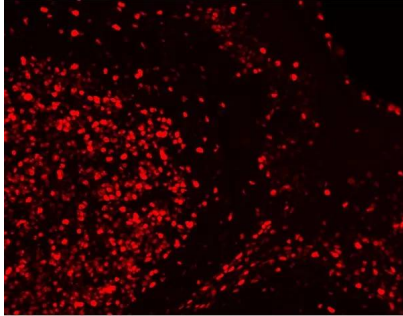
- [User Guide](#)
- [Safety Data Sheet](#)
- [Download CoA](#)

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



- [Datasheet](#)

GALLERY IMAGES



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