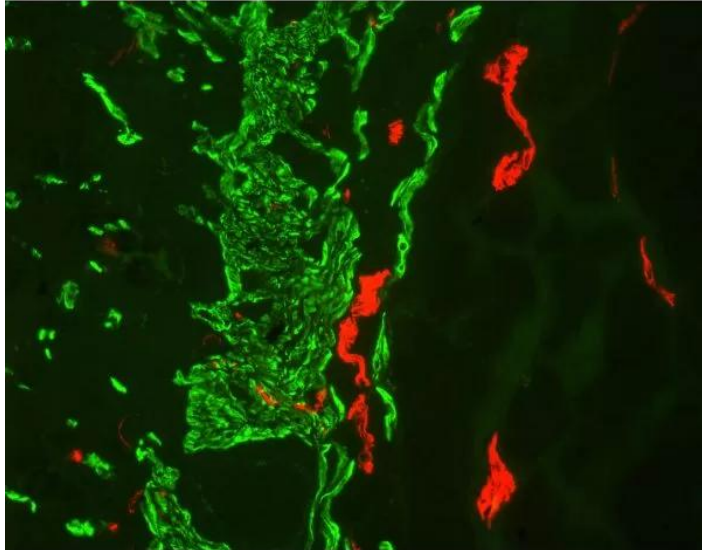




M.O.M. (MOUSE ON MOUSE) IMMUNODETECTION KIT, FLUORESC EIN

SKU: FMK-2201



DESCRIPTION

The Vector M.O.M. Fluorescein Kit is specifically designed to reduce endogenous mouse IgG staining when using mouse primary antibodies on mouse tissue. It contains a proprietary Mouse IgG Blocking Reagent, a M.O.M. Biotinylated Anti-Mouse Ig Reagent, and Fluorescein Avidin DCS and is based on a simple and easy to follow avidin/biotin protocol.

Key advantages of the Vector M.O.M. Immunodetection Kits:

- Clear, crisp specific staining of antigens of interest
- Based on VECTASTAIN® ABC reagent protocols
- Procedures are simple and easy to follow
- No tedious calculations or antibody prebinding steps required

Kit Contents:

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



- 6 ml of M.O.M. Protein Concentrate
- 1 ml M.O.M. Mouse Ig Blocking Reagent
- 0.1 ml M.O.M. Biotinylated Anti-Mouse IgG Reagent
- 0.4 ml Fluorescein Avidin DCSThe Vector® M.O.M. Immunodetection Kit contains enough stock reagents to produce about 25 ml of working solution which is generally sufficient to stain approximately 250 tissue sections.

SPECIFICATIONS

Blocking Action	Endogenous Mouse Ig
Color of Fluorescence	Green
Format	Concentrate
Isolation Technology	ABC (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems
Unit Size	1 Kit
Applications	Immunofluorescence
Target Species	Mouse
Conjugate	Fluorescein
Platform	ABC (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems
Source Species	Mouse

TECHNICAL INFORMATION

The inability of an anti-mouse secondary antibody to distinguish between a primary antibody produced in mice and the endogenous mouse immunoglobulins present in mouse tissue results in high background staining which obscures specific staining. This problem can be essentially eliminated by using a Vector M.O.M. Kit. The result is clear, crisp specific staining of the antigens of interest. Excellent staining results for a once difficult application have now become routine with the Vector M.O.M. System.

The Vector M.O.M. Immunodetection Kits are useful in any mouse-on-mouse application, such as studies in genetically engineered mice, including transgenic and knock-out models, as well as mouse xenografts.

The M.O.M. Kits can be also used in multiple antigen staining protocols to localize two different mouse primary antibodies on the same mouse tissue sections using fluorescent or enzyme-

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



based detection systems. Detailed information for multiple antigen labeling can be found in our brochure, *Discovery Through Color: A Guide to Multiple Antigen Labeling*, available on our website.

Four Vector M.O.M. Kits are available. The three kits based on avidin/biotin technology use the same proprietary M.O.M. Mouse Ig Blocking Reagent and the M.O.M. Biotinylated Anti-Mouse Ig Reagent, but offer a choice of an enzyme-based or fluorescent-based visualization method. M.O.M. Mouse Ig Blocking Reagent and M.O.M. Biotinylated Anti-Mouse Ig Reagent are also available separately.

The Vector M.O.M. ImmPRESS® Kit contains our proprietary M.O.M. Mouse Ig Blocking Reagent paired with a specialized, ready-to-use, one-step, non-biotin M.O.M. ImmPRESS™ Peroxidase Polymer reagent.

CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

- [M.O.M. \(Mouse on Mouse\) Troubleshooting Guide](#)
- [User Guide](#)
- [Protocol for Multiple Immunofluorescent Staining Using Mouse Antibodies](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES

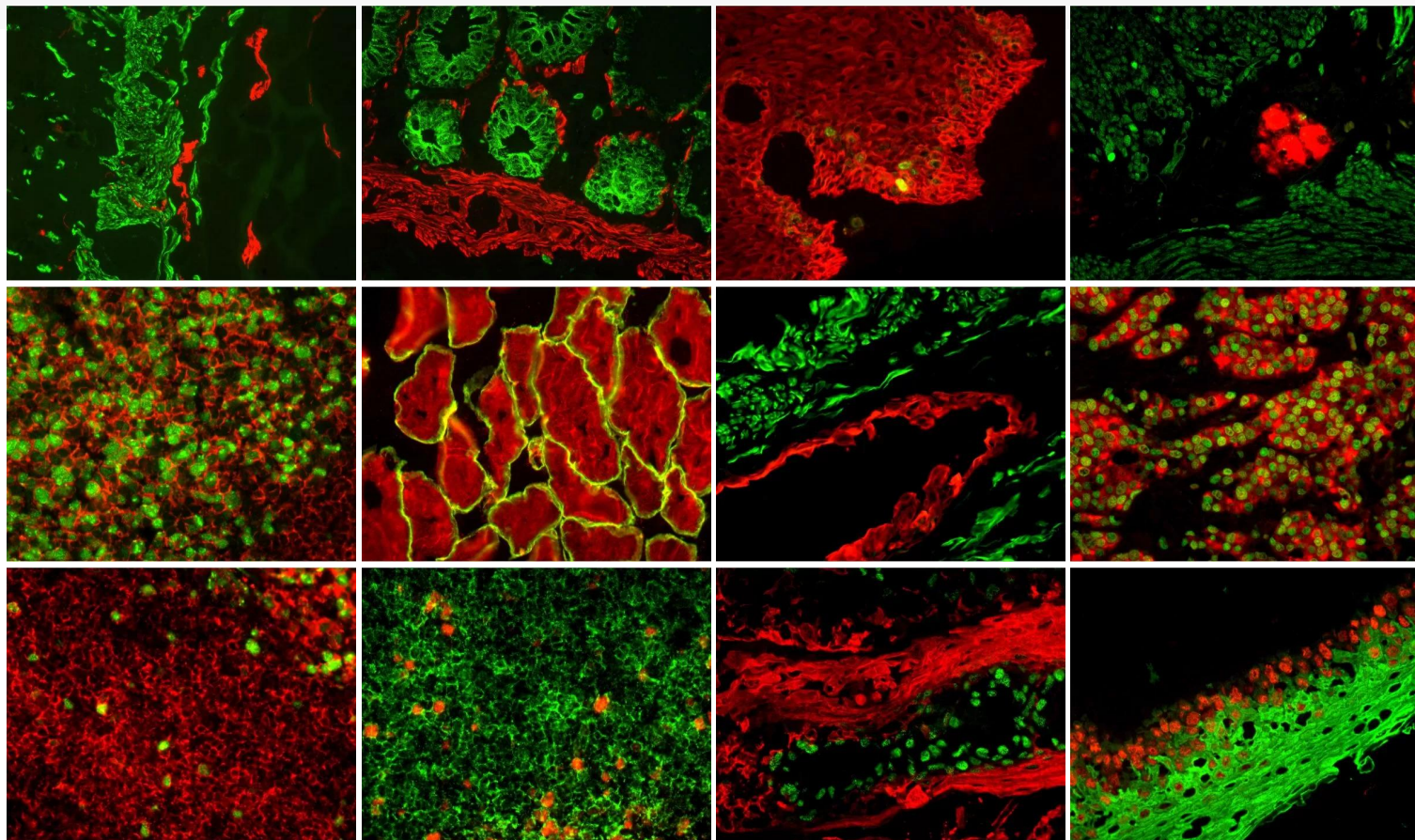
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



www.vectorlabs.com

Email: customerservice@vectorlabs.com

Telephone: [\(650\) 697-3600](tel:(650)697-3600)



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

[M.O.M. \(Mouse on Mouse\) Immunodetection Kit, Fluorescein](https://vectorlabs.com/products/mouse-on-mouse-fluorescein-kit/)

<https://vectorlabs.com/products/mouse-on-mouse-fluorescein-kit/>