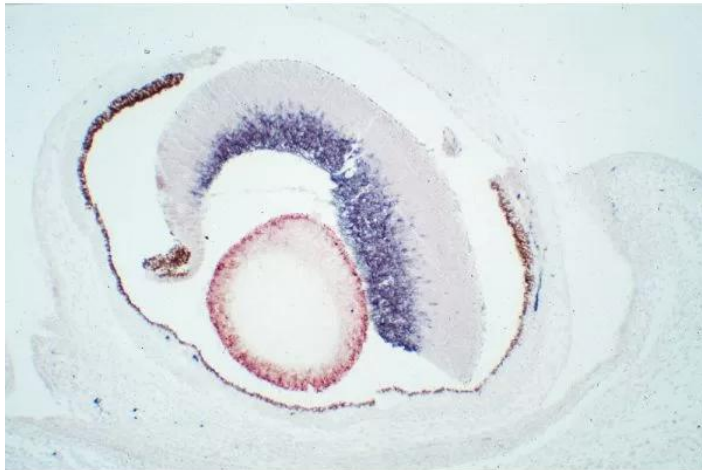




M.O.M. (MOUSE ON MOUSE) ELITE® IMMUNODETECTION KIT, PEROXIDASE

SKU: PK-2200



DESCRIPTION

The M.O.M. (Mouse on Mouse) Elite Peroxidase Kit significantly reduces endogenous mouse Ig staining when using mouse primary antibodies on mouse tissue. It is based on VECTASTAIN® ABC reagent protocols.

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

- Clear, crisp specific staining of antigens of interest
 - Procedures are simple and easy to follow
 - No tedious calculations or antibody prebinding steps required
- The 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.

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



SPECIFICATIONS

Blocking Action	Endogenous Mouse Ig
Detection Enzyme	Peroxidase
Format	Concentrate
Isolation Technology	ABC Elite (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems
Unit Size	1 Kit
Applications	Immunohistochemistry / Immunocytochemistry, Blotting Applications
Target Species	Mouse
Platform	ABC Elite (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems

KIT COMPONENTS

- 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
- VECTASTAIN Elite ABC Reagents A(1 ml) and B (1 ml)

TECHNICAL INFORMATION

The Vector M.O.M. Immunodetection Kits are specifically designed to reduce endogenous mouse IgG staining when using mouse primary antibodies on mouse tissue. 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. Elite Kit. The result is clear, crisp specific staining of the antigens of interest.

The Vector M.O.M. Elite Kit contains a proprietary Mouse IgG Blocking Reagent and is based on the simple and easy to follow VECTASTAIN® ABC reagent protocols; no tedious calculations or antibody pre-binding steps are required. Excellent staining results for a once difficult application have now become routine with the Vector M.O.M. System.

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



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-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.

Choose from a wide variety of peroxidase substrates available for chromogenic development.

Product review on Biocompare.com.

CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

- [User Guide](#)
- [M.O.M. \(Mouse on Mouse\) Troubleshooting Guide](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

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

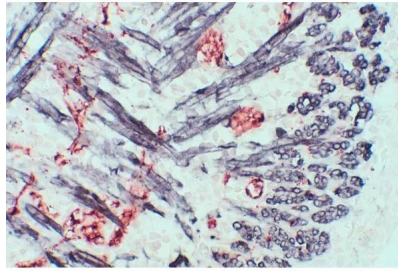
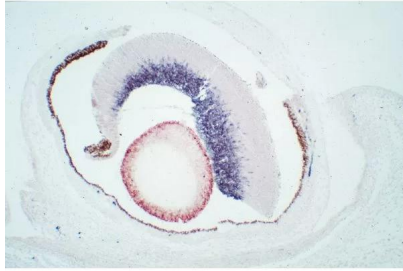


www.vectorlabs.com

Email: customerservice@vectorlabs.com

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

GALLERY IMAGES



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

[M.O.M. \(Mouse on Mouse\) Elite® Immunodetection Kit, Peroxidase](https://vectorlabs.com/products/mouse-on-mouse-elite-hrp-kit/)

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