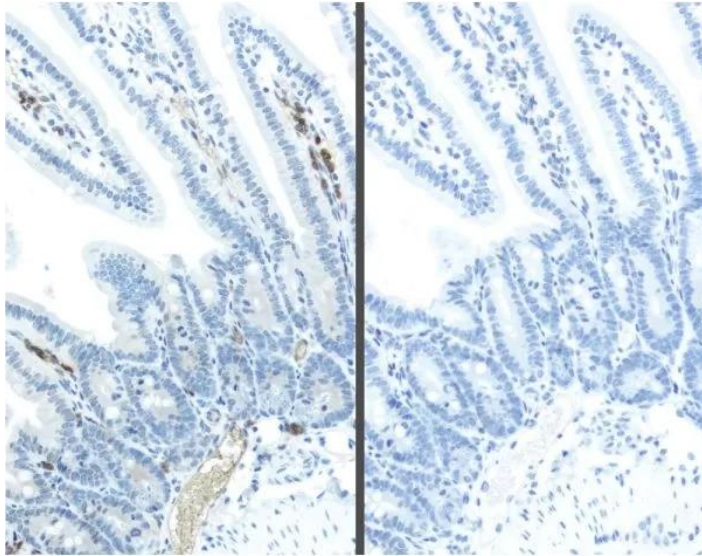




M.O.M. (MOUSE ON MOUSE) IMPRESS HRP (PEROXIDASE) POLYMER KIT

SKU: MP-2400



DESCRIPTION

The M.O.M. ImmPRESS Kit contains our proprietary M.O.M. Mouse Ig Blocking Reagent paired with a specialized, ready-to-use, one-step M.O.M. ImmPRESS Peroxidase Polymer reagent. It provides a significant reduction of endogenous mouse Ig staining when using mouse primary antibodies on mouse tissue.

Features:

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

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



SPECIFICATIONS

Blocking Action	Endogenous Mouse Ig
Detection Enzyme	Peroxidase
Isolation Technology	Micropolymer Reagents, Mouse on Mouse (M.O.M.) Systems
Unit Size	1 Kit
Applications	Immunohistochemistry / Immunocytochemistry
Target Species	Mouse
Conjugate	Micropolymer HRP
Platform	Micropolymer Reagents, Mouse on Mouse (M.O.M.) Systems

KIT COMPONENTS

- 0.6 ml concentrated stock solution of M.O.M. Mouse Ig Blocking Reagent
- 30 ml ready-to-use (R.T.U.) 2.5% Normal Horse Serum (NHS) for general protein blocking
- 15 ml R.T.U. M.O.M. ImmPRESS Anti-Mouse Ig Reagent (made in horse, ready-to-use)

TECHNICAL INFORMATION

For years, researchers attempting to localize mouse primary antibodies on mouse tissue have relied on the Vector[®] M.O.M. Kit to produce crisp, strong, specific staining of antibody targets. The Vector M.O.M. Kits are specifically designed to significantly reduce endogenous mouse Ig staining when using mouse primary antibodies on mouse tissue. The inability of an anti-mouse detection 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 Vector M.O.M. Kits. The result is clear, crisp staining of the antigens of interest. Tedious calculations and primary antibody pre-binding steps are not required.

The M.O.M. ImmPRESS HRP polymer reagent is affinity purified and specifically optimized for use with the M.O.M. Mouse Ig Blocking Reagent. The highly active peroxidase micropolymers of the ImmPRESS Detection System provide outstanding sensitivity, limit steric interference, and enhance accessibility. 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.



Key advantages of the ImmPRESS Polymer System:

- High sensitivity
- Low background
- Micropolymer technology
- Enhanced accessibility to nuclear and membrane antigens
- One-step detection
- Ready-to-use in a convenient dropper bottle
- Shorter assay time
- Simplified multiple labeling

The Vector M.O.M. ImmPRESS Kit (Cat. No. MP-2400) contains sufficient reagents to stain approximately 150 mouse sections.

CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

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

GALLERY IMAGES

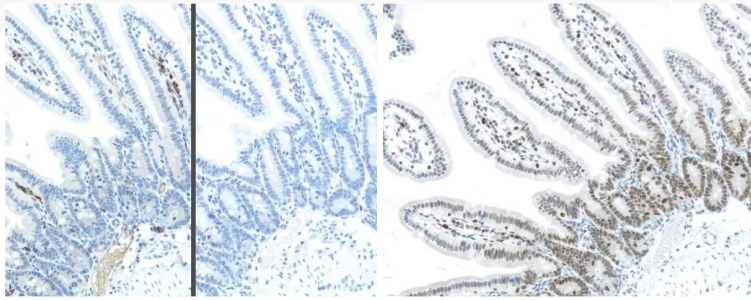
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



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. \(Mouse on Mouse\) ImmPRESS HRP \(Peroxidase\) Polymer Kit](https://vectorlabs.com/products/mouse-on-mouse-immpress-hrp-polymer-kit/)

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