

# M.O.M.<sup>®</sup> (Mouse on Mouse) ImmPRESS<sup>®</sup> HRP Polymer Kit

Peroxidase

**Cat. No.** MP-2400

**Storage** Store reagents in original bottles at 2–8 °C. Do not freeze.

**Description** The Vector<sup>®</sup> M.O.M. (Mouse on Mouse) Immunodetection Kits are specifically designed to localize mouse primary antibodies on mouse tissues.

The M.O.M. Immunodetection Kit can be used with normal and genetically engineered mouse models, including transgenic, xenograft, knock out and other mutant strains.

## Kit Components

Product Name	Volume
M.O.M. Mouse IgG Blocking Reagent	0.6 ml
M.O.M. Normal Horse Serum, 2.5% (ready to use)	30 ml
M.O.M. ImmPRESS Horse Anti-Mouse IgG Polymer Reagent (ready to use)	15 ml

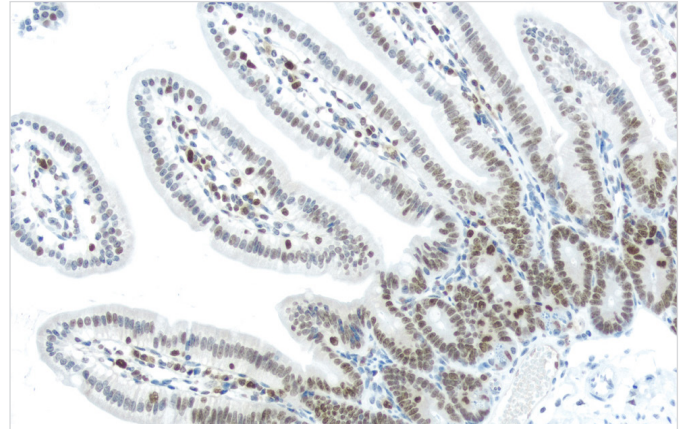
The M.O.M. ImmPRESS Polymer Kit contains enough reagents to stain approximately 75–150 mouse sections.

## Preparation of M.O.M. Working Solutions

- M.O.M. Mouse IgG Blocking Reagent: add 2 drops (~90 µl) of M.O.M. Blocking Reagent stock solution to 2.5 ml of PBS or TBS. †
- † PBS: 10 mM sodium phosphate, 0.15 M NaCl, pH 7.4–7.8  
TBS: 50 mM TRIS, 0.15 M NaCl, pH 7.5–7.8

## Staining Procedure

1. For paraffin sections, deparaffinize and hydrate tissue sections through xylenes or other clearing agents and graded alcohol series.  
For frozen sections or cell preparations, fix with acetone or an appropriate fixative for the antigen under study. Air dry.  
Rinse for 5 minutes in tap water.
2. If antigen unmasking is required, perform this procedure using a Vector<sup>®</sup> Antigen Unmasking Solution, Citrate-based, pH 6.0 (H-3300) or Tris-based, pH 9.0 (H-3301).
3. Block endogenous enzyme activity, if necessary, by incubating sections with BLOXALL<sup>®</sup> Blocking Solution (SP-6000) for 10 minutes.



Mouse Intestine - Single Label: Ki67 (m) detected with M.O.M. ImmPRESS Polymer Kit and Vector DAB Substrate Kit (brown). Counterstained with Hematoxylin (blue).

4. Wash 2 x 2 minutes in PBS or TBS.
  5. Incubate for 1 hour in working solution of prepared M.O.M. Mouse IgG Blocking Reagent.
  6. Wash 2 x 2 minutes in PBS or TBS.\*
  7. Incubate for 5 minutes in M.O.M. Normal Horse Serum, 2.5% (NHS)\*.
  8. Tip off excess NHS. Dilute primary antibody in NHS. Incubate section in diluted primary antibody for 30 minutes.\*
  9. Wash for 2 x 2 minutes in PBS or TBS.\*
  10. Apply the M.O.M. ImmPRESS Reagent and incubate for 10 minutes.\*
  11. Wash for 2 x 2 minutes in PBS or TBS.
  12. Prepare and apply peroxidase substrate solution (not included in kit) according to substrate kit instructions.
- \* It is recommended that the exact times described in steps 6–10 be used in the staining protocol. Longer incubation may result in an increase in background staining.

Detailed product listings, specifications, protocols and additional support information such as our Troubleshooting Guide: Mouse Antibodies on Mouse Tissue are available on our website: [vectorlabs.com](http://vectorlabs.com)