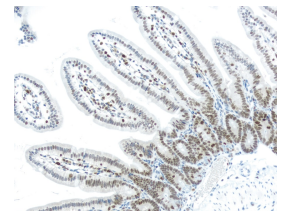
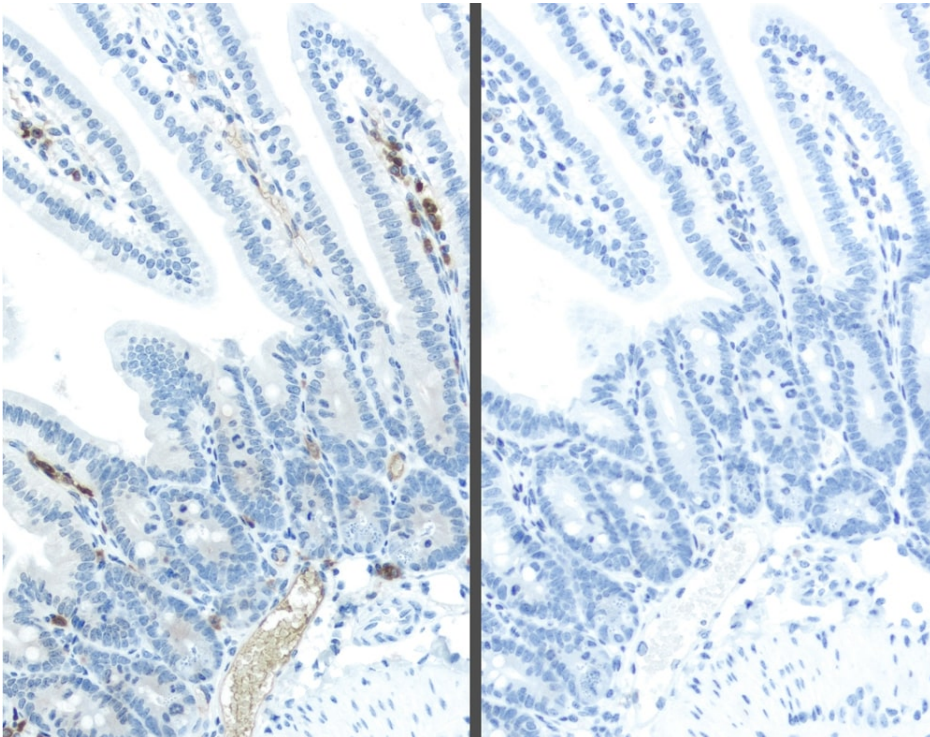




# M.O.M.<sup>®</sup> (Mouse on Mouse) ImmPRESS<sup>®</sup> HRP (Peroxidase) Polymer Kit

## MP-2400

Product Images



## Short 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

## Additional Information

---

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

