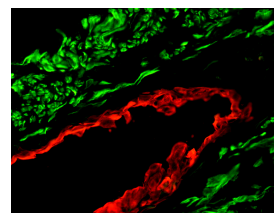
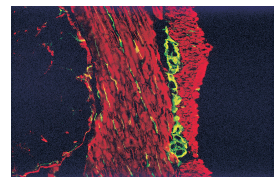
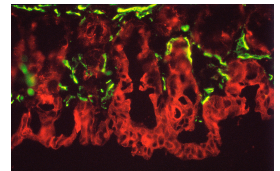
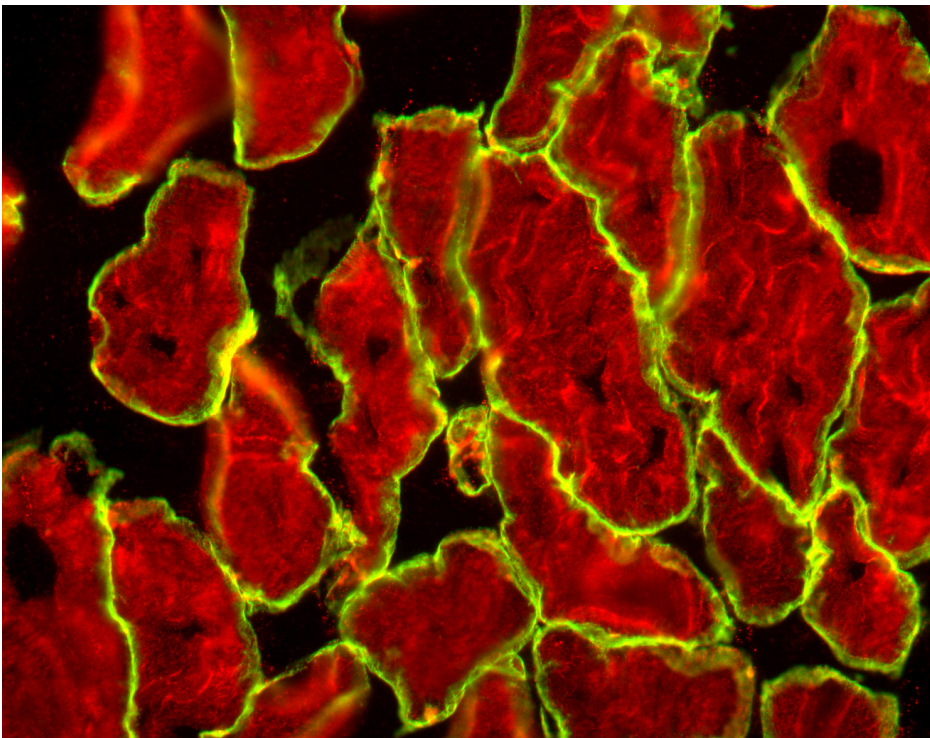


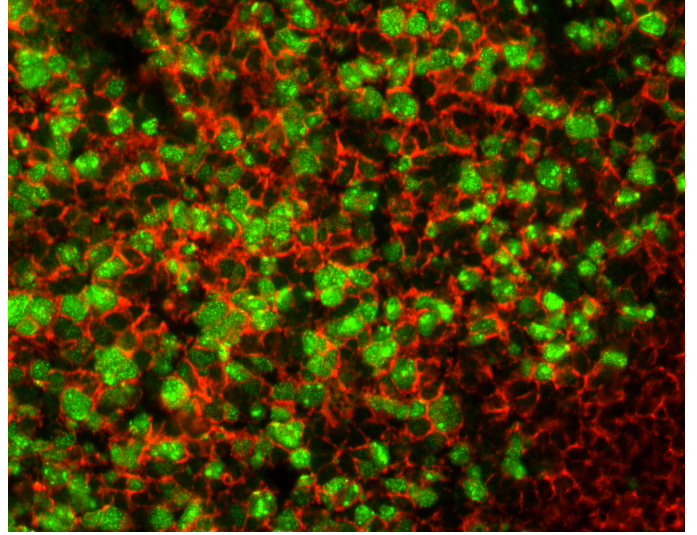
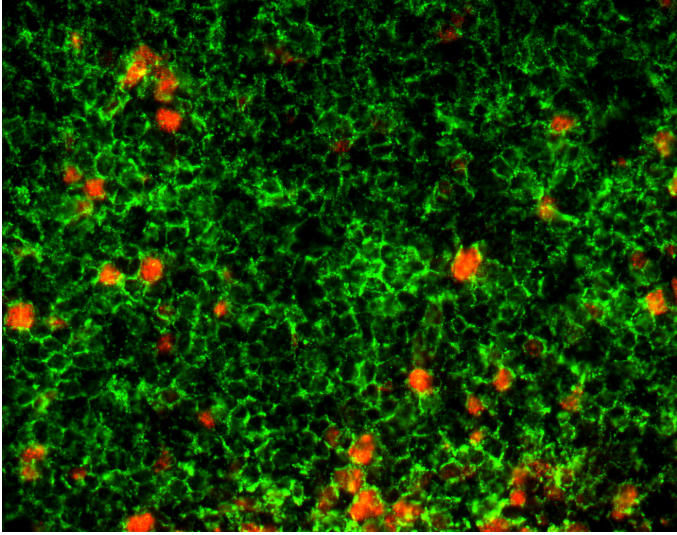


M.O.M.® (Mouse on Mouse) Immunodetection Kit, Basic

BMK-2202

Product Images





Short Description

The M.O.M. Basic Kit is 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. Kit 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.

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

- Significant reduction of endogenous mouse Ig staining when using mouse primary antibodies on mouse tissue
- Clear, crisp specific staining of antigens of interest
- Based on VECTASTAIN® ABC reagent protocols
- No tedious calculations or antibody prebinding steps required
- Contains enough stock reagents to produce about 25 ml of working solution, which is generally sufficient to stain approximately 250 tissue sections.

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence
Technology	ABC (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems
Target Species	Mouse
Format	Concentrate
Blocking Action	Endogenous Mouse Ig

