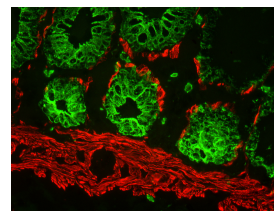
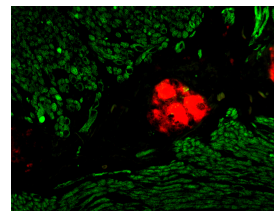
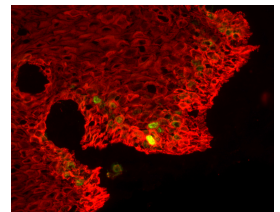
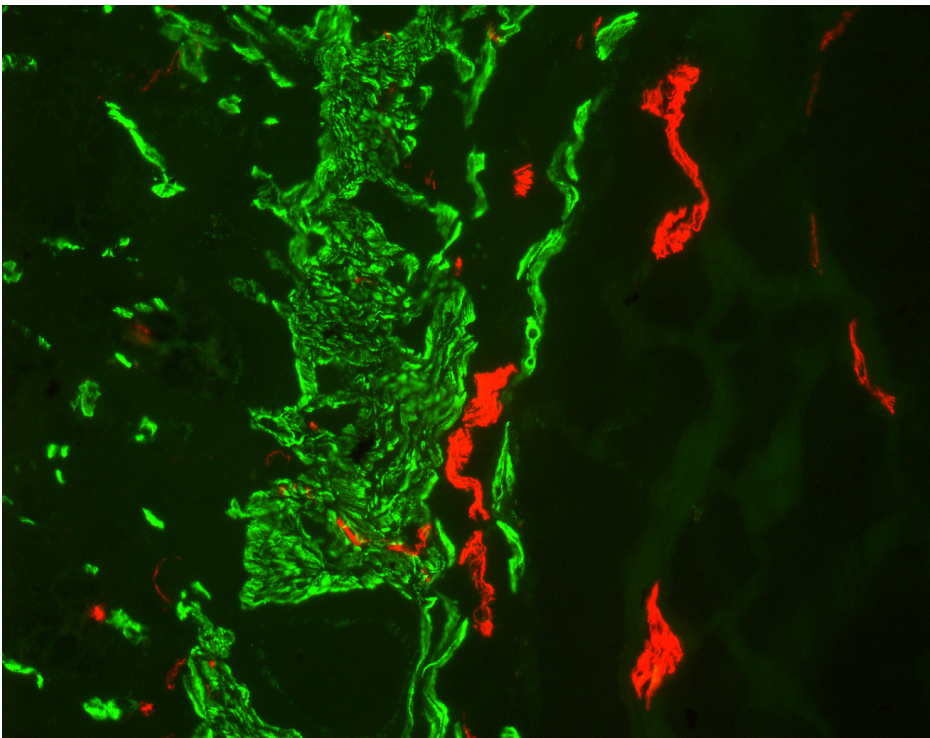


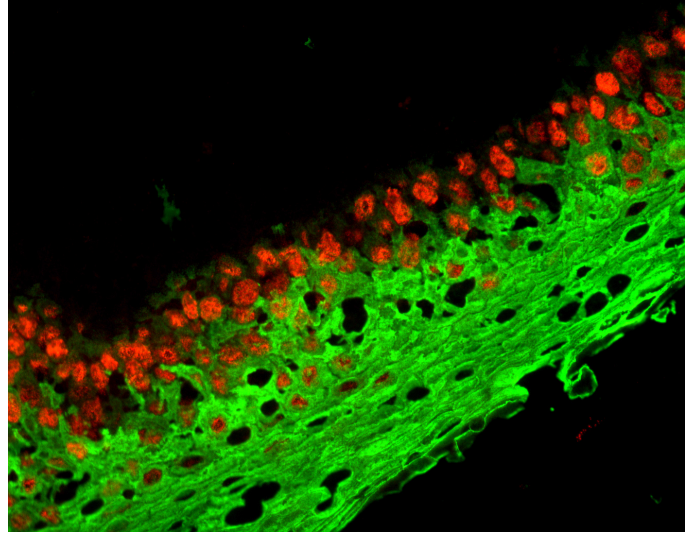
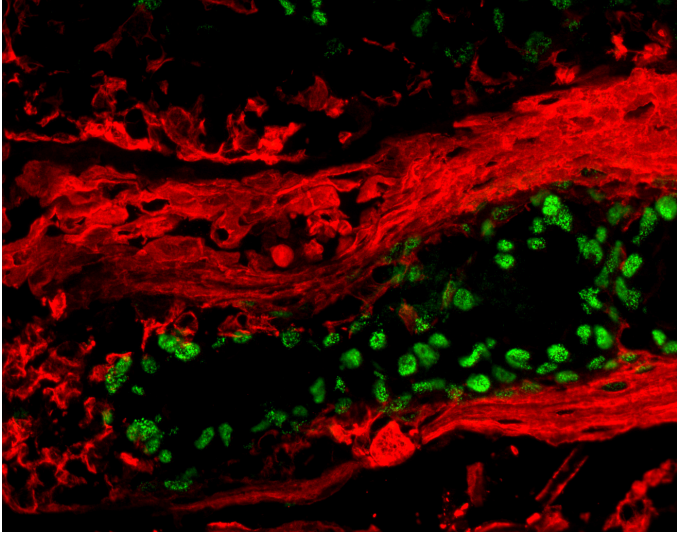


M.O.M.® (Mouse on Mouse) Immunodetection Kit, Fluorescein

FMK-2201

Product Images





Short Description

The Vector M.O.M. Fluorescein Kit is specifically designed to reduce endogenous mouse IgG staining when using mouse primary antibodies on mouse tissue. It contains a proprietary Mouse IgG Blocking Reagent, a M.O.M. Biotinylated Anti-Mouse Ig Reagent, and Fluorescein Avidin DCS and is based on a simple and easy to follow avidin/biotin protocol.

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

- Clear, crisp specific staining of antigens of interest
- Based on VECTASTAIN® ABC reagent protocols
- Procedures are simple and easy to follow
- No tedious calculations or antibody prebinding steps required

Kit Contents:

- 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
- 0.4 ml Fluorescein Avidin DCS

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.

Additional Information

Unit Size	1 kit
Applications	Immunofluorescence
Technology	ABC (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems
Conjugate	Fluorescein
Color of Fluorescence	Green
Target Species	Mouse
Format	Concentrate
Blocking Action	Endogenous Mouse Ig

