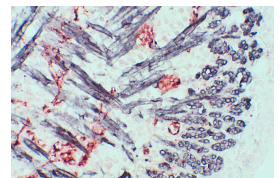
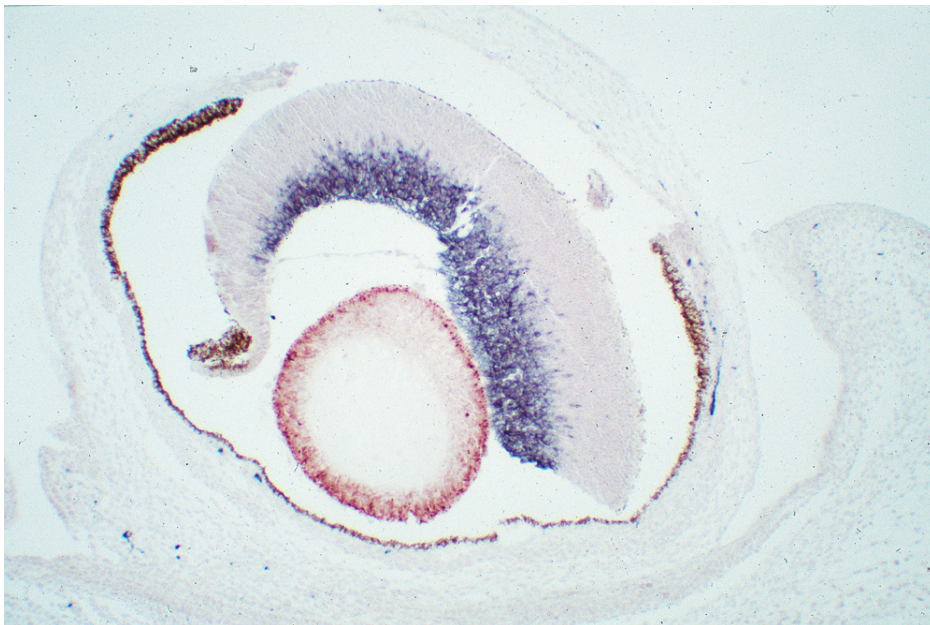




M.O.M.® (Mouse on Mouse) Elite® Immunodetection Kit, Peroxidase

PK-2200

Product Images



Short Description

The M.O.M. (Mouse on Mouse) Elite Peroxidase Kit significantly reduces endogenous mouse Ig staining when using mouse primary antibodies on mouse tissue. It is based on VECTASTAIN® ABC reagent protocols.

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

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

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	Immunohistochemistry / Immunocytochemistry, Blotting Applications
Technology	ABC Elite (Avidin/Biotin) Systems, Mouse on Mouse (M.O.M.) Systems
Detection Enzyme(s)	Peroxidase
Target Species	Mouse
Format	Concentrate
Blocking Action	Endogenous Mouse Ig

