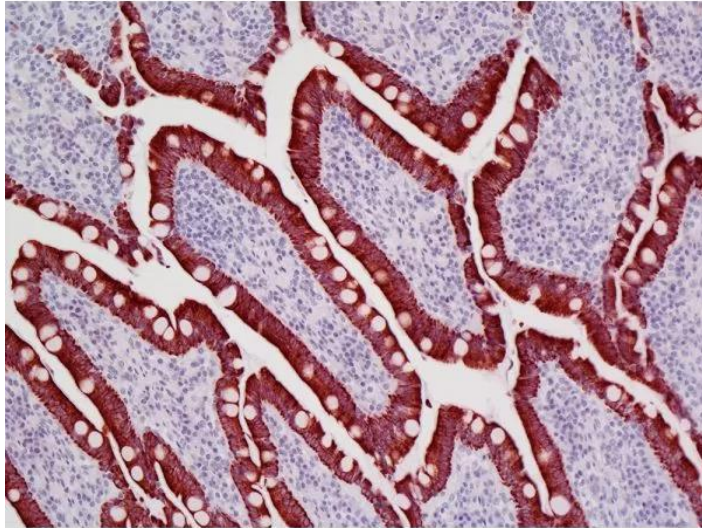




IMPRESS-VR HORSE ANTI-MOUSE IGG POLYMER DETECTION KIT, PEROXIDASE, VETERINARY REAGENT

SKU: MP-6402-15



DESCRIPTION

The ImmPRESS Polymerized Reporter Enzyme Staining System is an enzymatic, one-step detection kit. ImmPRESS Polymer Reagents consist of unique micropolymers of highly active peroxidase enzyme attached to highly cross-adsorbed, affinity purified secondary antibodies. The ImmPRESS Polymer Reagents produce outstanding immunohistochemistry and immunocytochemistry results due to increased target accessibility, binding specificity, and signal intensity along with low background staining.

ImmPRESS **VR** Polymer Detection Reagents are additionally cross-adsorbed to ensure minimal cross-reactivity against endogenous tissue elements in animal species commonly used for diagnostics and research-based animal model systems.

Features:

- Enhanced accessibility to nuclear and membrane antigens

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



- Ready-to-use in a convenient dropper bottle
- Shorter assay time
- Simplified multiple labeling
- Minimal cross-reactivity with bovine, goat, sheep, pig, horse, cat, dog, and rabbit (as well as human). For mouse and rat tissues, separate kits are recommended (for rat, MP-7422; for mouse, MP-2400).

SPECIFICATIONS

Detection Enzyme	Peroxidase
Format	Ready-to-Use
Isolation Technology	Micropolymer Reagents
Unit Size	15 ml
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, ELISAs
Target Species	Mouse
Conjugate	Micropolymer HRP
Platform	Micropolymer Reagents

KIT COMPONENTS

- ImmPRESS VR Polymer Anti-Mouse IgG Reagent (made in horse, ready-to-use)
- 2.5% Normal Horse Serum for blocking (ready-to-use)
- The ImmPRESS (15 ml) Kit will stain approximately 75-150 sections.

TECHNICAL INFORMATION

For mouse and rat tissues, separate kits are recommended (for rat, MP-7422; for mouse, MP-2400).

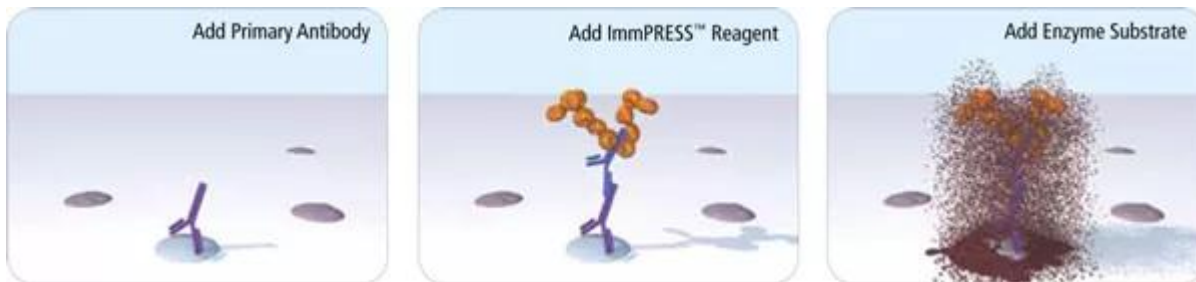
Staining Procedure

The staining procedure is simple as shown in the diagram below. Following a blocking step with the diluted normal horse serum, sections are incubated with primary antibody. After a brief wash, the appropriate ImmPRESS Reagent is added to the sections and incubated for 30

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



minutes. Sections are again rinsed and the slides are developed with the peroxidase substrate of choice.



Product review on [Biocompare.com](https://www.biocompare.com).

CITATIONS



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

- [ImmPRESS Polymer & Automated Staining App. Note](#)
- [ImmPRESS Troubleshooting Guide](#)
- [User Guide](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES

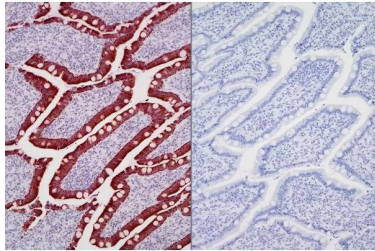
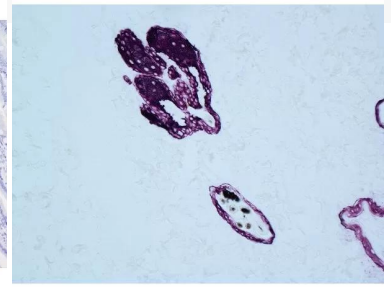
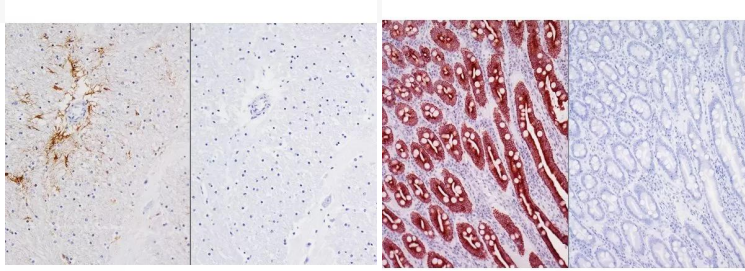
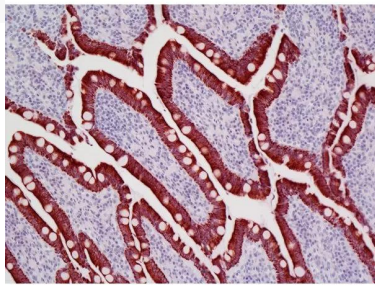
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



www.vectorlabs.com

Email: customerservice@vectorlabs.com

Telephone: [\(650\) 697-3600](tel:(650)697-3600)



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

[ImmPRESS-VR Horse Anti-Mouse IgG Polymer Detection Kit, Peroxidase, Veterinary Reagent](https://vectorlabs.com/products/immpress-hrp-vr-horse-anti-mouse-igg-kit/)

<https://vectorlabs.com/products/immpress-hrp-vr-horse-anti-mouse-igg-kit/>