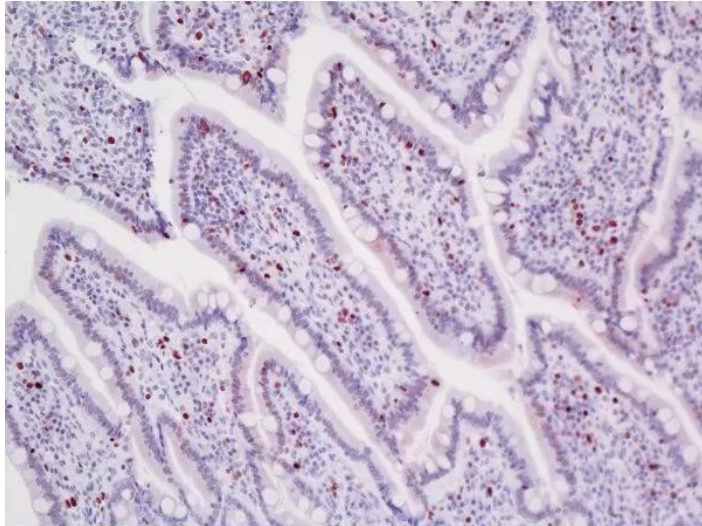




IMPRESS-VR HORSE ANTI-RABBIT IGG POLYMER DETECTION KIT, PEROXIDASE, VETERINARY REAGENT

SKU: MP-6401-15



DESCRIPTION

ImmPRESS VR Polymer Detection Reagents are cross-adsorbed to ensure minimal cross-reactivity against endogenous tissue elements in animal species commonly used for diagnostics and research-based animal model systems. ImmPRESS VR Polymer HRP Anti-Rabbit IgG is designed to be used on the following tissues: bovine, goat, sheep, swine, horse, cat, dog, rat and mouse (as well as human).

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.

Features:

- High sensitivity
- Low background

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



- One-step detection
- Micropolymer technology
- Enhanced accessibility to nuclear and membrane antigens
- Ready-to-use in a convenient dropper bottle
- Shorter assay time
- Simplified multiple labeling

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	Rabbit
Conjugate	Micropolymer HRP

KIT COMPONENTS

- ImmPRESS VR Polymer HRP Anti-Rabbit 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

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.

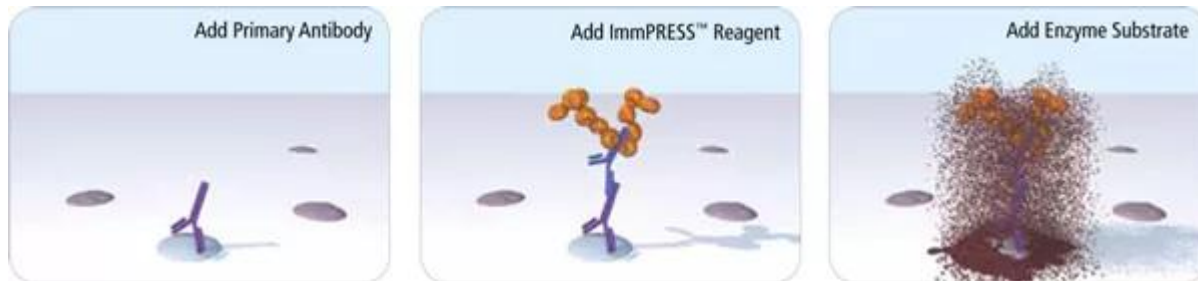
Staining Procedure

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

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



wash, the appropriate ImmPRESS Reagent is added to the sections and incubated for 30 minutes. Sections are again rinsed and the slides are developed with the peroxidase substrate of choice.

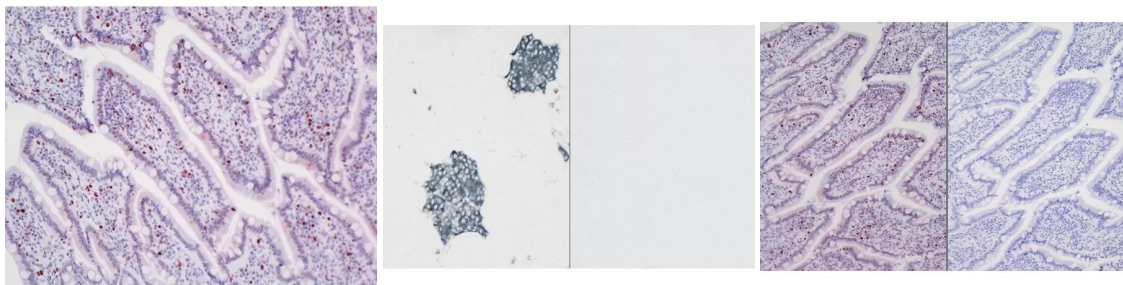


Product review on Biocompare.com.

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.