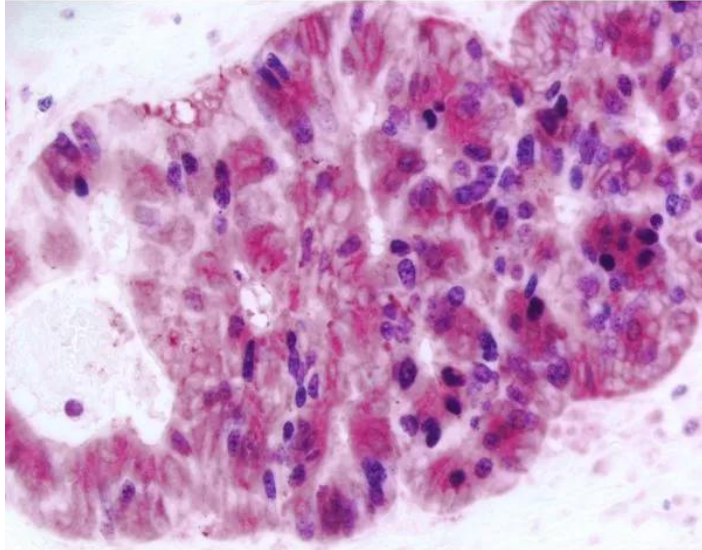




IMPRESS-AP HORSE ANTI-RABBIT IGG POLYMER DETECTION KIT, ALKALINE PHOSPHATASE

SKU: MP-5401



DESCRIPTION

The ImmPRESS polymerized reporter enzyme staining system uses novel conjugation and micropolymer chemistries to create a highly sensitive, ready-to-use, one-step detection system. This unique micropolymer of highly active enzyme is attached to our affinity purified secondary antibodies, producing reagents with outstanding sensitivity and low background.

Features:

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

Kit Contents:

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



- ImmPRESS-AP (alkaline phosphatase) Polymer Anti-Rabbit IgG Reagent (ready-to-use)
- 2.5% Normal Horse Serum for blocking (ready-to-use) The ImmPRESS (15 ml) Kit will stain approximately 75-150 sections and the ImmPRESS (50 ml) Kit will stain approximately 250-500 sections.

SPECIFICATIONS

Detection Enzyme	Alkaline Phosphatase
Format	Ready-to-Use
Isolation Technology	Micropolymer Reagents
Unit Size	15 ml, 50 ml
Applications	Immunohistochemistry / Immunocytochemistry, ELISAs, In situ hybridization
Target Species	Rabbit
Conjugate	Micropolymer AP
Source Species	Rabbit

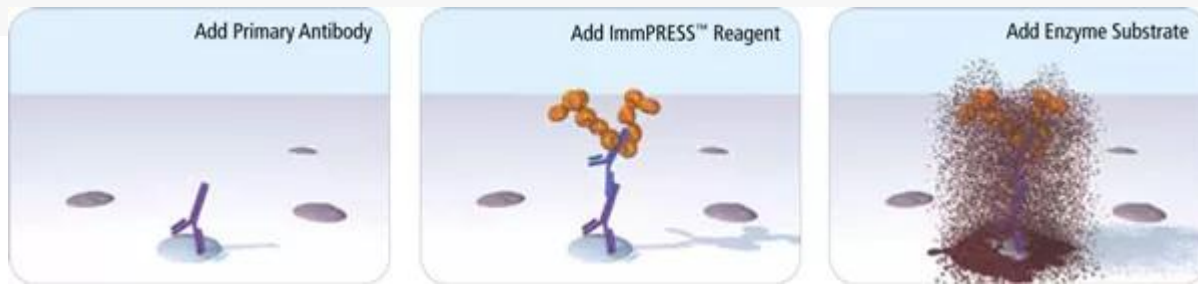
TECHNICAL INFORMATION

The ImmPRESS-AP Alkaline Phosphatase Polymer Reagent micropolymer technology limits steric interference and provides enhanced accessibility to the target avoiding the disadvantages of other polymer systems that use large dextrans or other macromolecules as backbones. The result is crisp, strong staining of antibody targets, especially nuclear and membrane antigens, and greater sensitivity than conventional alkaline phosphatase antibody conjugates and other alkaline phosphatase polymer systems for immunohistochemistry and other applications.

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

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



Consider Species Cross-Reactivity

When choosing the optimal detection system for your application, it is important to consider not only the species of the primary antibody but also the species of the tissue. If the species of the primary antibody and the species of the tissue are closely related (e.g. rat and mouse), the secondary antibody may bind to endogenous IgG in the tissue section leading to background. The following options minimize background staining in these instances:

- Use a secondary antibody specifically adsorbed to remove cross-reacting antibodies of closely-related species (e.g. ImmPRESS Anti-Mouse IgG, Rat Adsorbed).
- Use the M.O.M. ImmPRESS Kit (Cat. No. MP-2400) for applications of mouse primary antibodies on mouse tissue.

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](#)

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

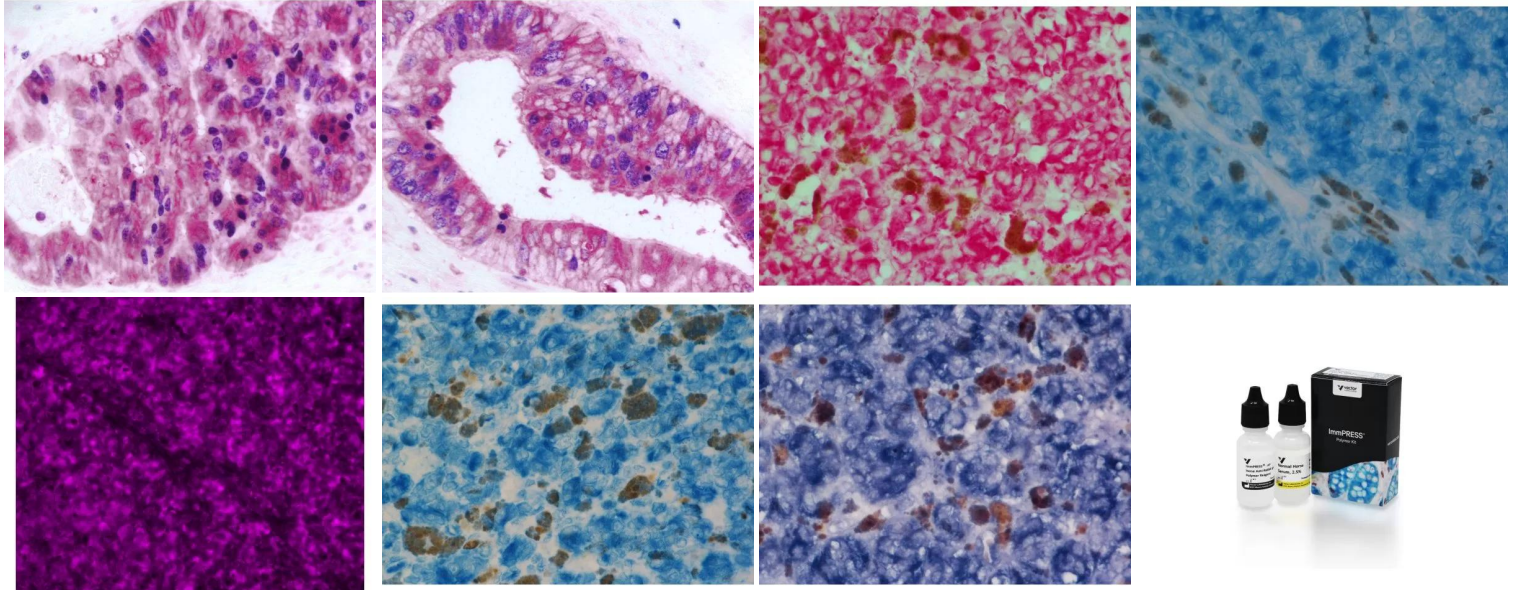


www.vectorlabs.com

Email: customerservice@vectorlabs.com

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

GALLERY IMAGES



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

[ImmPRESS-AP Horse Anti-Rabbit IgG Polymer Detection Kit, Alkaline Phosphatase](https://vectorlabs.com/products/immpress-anti-rabbit-igg-ap-kit/)

<https://vectorlabs.com/products/immpress-anti-rabbit-igg-ap-kit/>