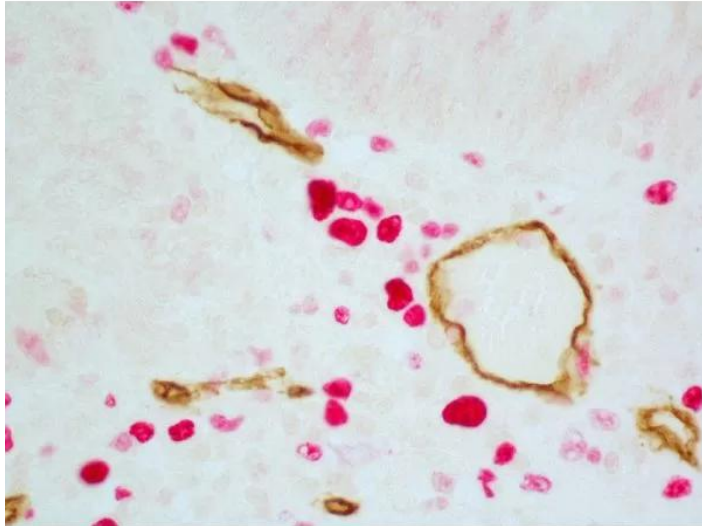




IMPRESS DUET DOUBLE STAINING HRP/AP POLYMER KIT (ANTI-RABBIT IGG-BROWN, ANTI-MOUSE IGG-MAGENTA)

SKU: MP-7714-15



DESCRIPTION

This ImmPRESS Duet Double Staining Polymer Kit enables fast, well-defined localization and visualization of two different target antigens on the same tissue section. Utilizing our unique ImmPRESS micropolymer chemistry, we independently conjugate highly-active horseradish peroxidase (HRP) and alkaline phosphatase (AP) enzyme polymers to anti-rabbit IgG and anti-mouse IgG affinity-purified, highly cross-adsorbed secondary antibodies.

Features:

Save time and streamline your double staining IHC workflow with the ImmPRESS Duet Double Staining Polymer Kit.

- Complete Kits for IHC Double Labeling
- Convenient Dual Specificity (anti-mouse IgG/rabbit IgG) ImmPRESS HRP/AP Polymer Reagent

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



- High Sensitivity and Specificity
- Non-Biotin Based Reagents
- Prediluted, Ready-to-use

Kit Contents:

- 15 ml BLOXALL® Blocking Solution
 - 15 ml 2.5% Normal Serum
 - 15 ml ImmPRESS Duet HRP/AP Polymer Reagent (mixture of ImmPRESS HRP Anti-Rabbit IgG and ImmPRESS AP Anti-Mouse IgG)
 - ImmPACT® DAB EqV Substrate (HRP, brown)
 - ImmPACT Vector® Red Substrate (AP, magenta)
- Note: The ImmPRESS Duet Double Staining Polymer Kit is intended for use on non-rodent tissue specimens.**

SPECIFICATIONS

Blocking Action	Endogenous Peroxidase, Endogenous Alkaline Phosphatase, Non-Specific Protein Blocking
Detection Enzyme	HRP/AP
Format	Ready-to-Use
Isolation Technology	Micropolymer Reagents
Unit Size	1 Kit
Applications	Immunohistochemistry / Immunocytochemistry
Target Species	Universal (Mouse/Rabbit)
Conjugate	Micropolymer HRP/Micropolymer AP
Platform	Micropolymer Reagents

TECHNICAL INFORMATION

Save time and streamline your double staining IHC workflow with the ImmPRESS Duet Double Staining Polymer Kit. These reagents are blended at optimized dilutions to provide a stable pre-diluted, ready-to-use HRP/AP formula that will detect mouse and rabbit primary antibodies, and facilitate a time-saving, simplified, one-step double label detection protocol. For maximum sensitivity and staining contrast between target antigens, ImmPACT DAB EqV (HRP, brown) and ImmPACT Vector® Red (AP, magenta) substrates are included to complete this easy to use

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



staining kit.

Reliable and reproducible results are obtained on tissue sections where the two target antigens do not over-lap (co-localize) in the same structure of the same cell, but rather are expressed in different cell compartments or different cell types.

Also included, BLOXALL Blocking Solution abolishes all endogenous peroxidase and alkaline phosphatase activity in a single 10 minute incubation. In addition, 2.5% Normal Serum blocks all non-specific binding of detection reagents to ensure crisp, specific staining.

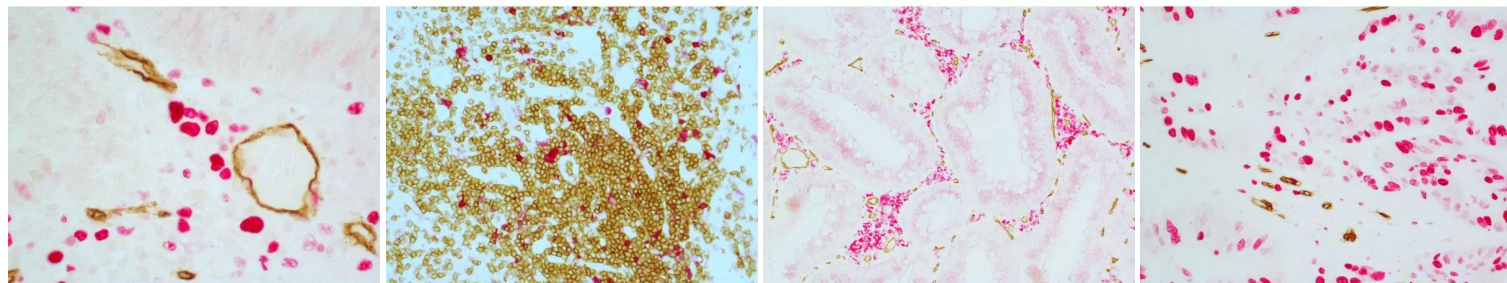
All reagents are supplied in stabilized, ready-to-use (R.T.U.) solutions for maximum convenience, and consistency between assays. The ImmPRESS Duet Double Staining Polymer Kits offer a complete and simple solution for visualizing target antigens in cell and tissue preparations and contains sufficient reagents to stain approximately 150 sections.

Product review on Biocompare.com.

DOCUMENTS

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

GALLERY IMAGES



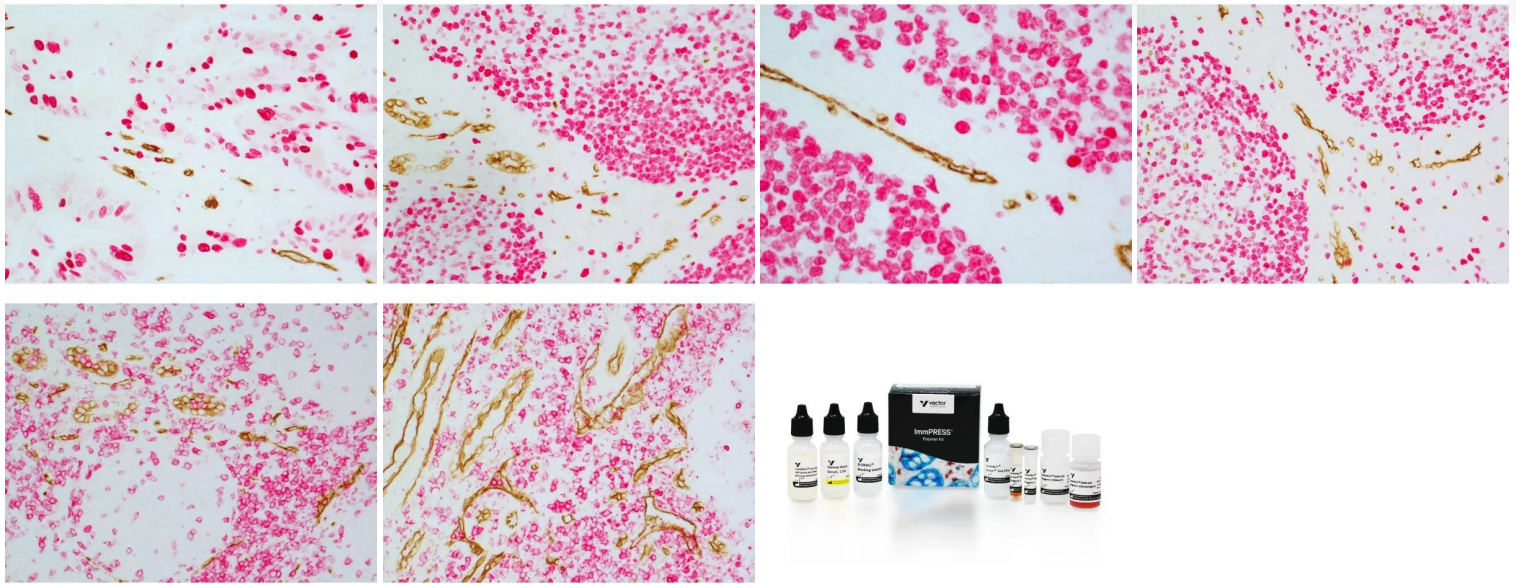
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



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 Duet Double Staining HRP/AP Polymer Kit \(Anti-Rabbit IgG-brown, Anti-Mouse IgG-magenta\)](https://vectorlabs.com/products/immpress-duet-staining-kit-rb-br-ms-red/)

<https://vectorlabs.com/products/immpress-duet-staining-kit-rb-br-ms-red/>