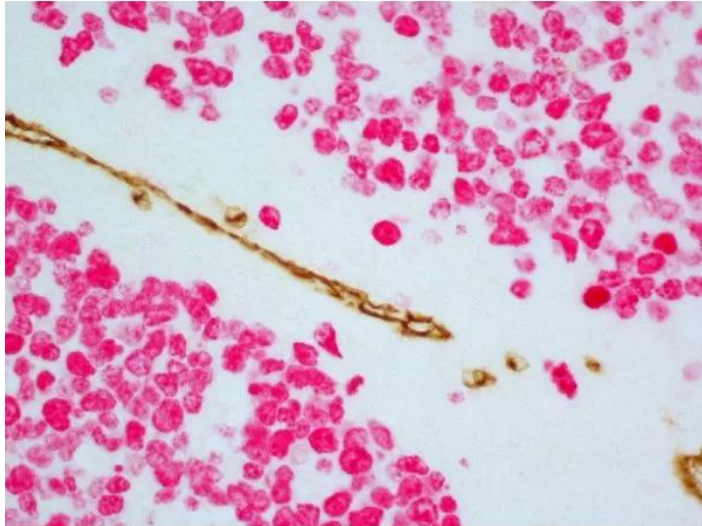




# **IMPRESS DUET DOUBLE STAINING HRP/AP POLYMER KIT (ANTI-MOUSE IGG-BROWN, ANTI-RABBIT IGG-MAGENTA)**

**SKU:** MP-7724-15



---

## **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-mouse IgG and anti-rabbit 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-Mouse IgG and ImmPRESS AP Anti-Rabbit 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

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

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



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 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](http://Biocompare.com).

## DOCUMENTS

- [Safety Data Sheet](#)
- [User Guide](#)
- [ImmPRESS Polymer & Automated Staining App. Note](#)
- [Download CoA](#)
- [Datasheet](#)
- [Datasheet2](#)

## GALLERY IMAGES

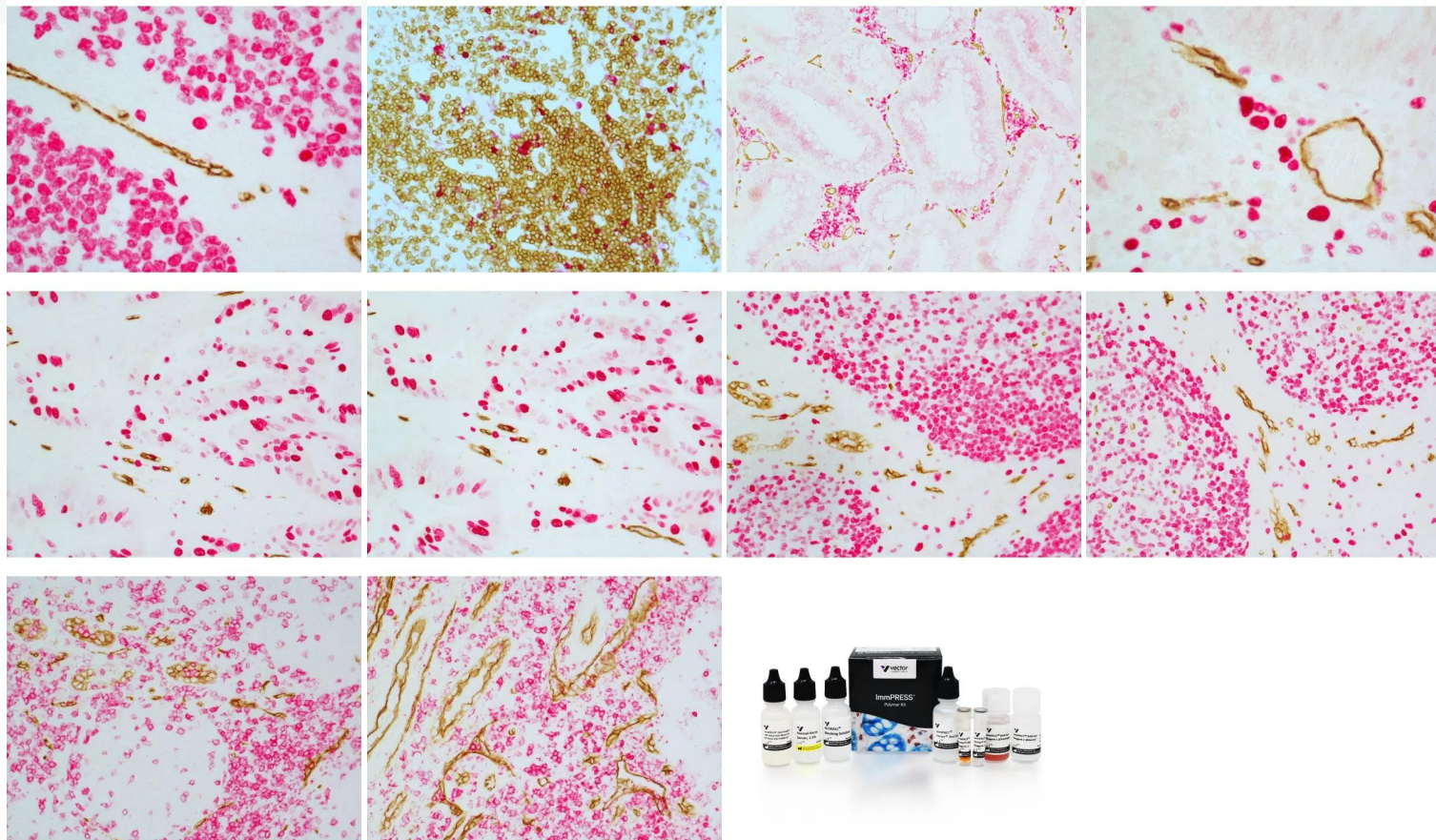
**For research use only. Not intended for therapeutic or diagnostic use in animals or humans.**



[www.vectorlabs.com](http://www.vectorlabs.com)

Email: [customerservice@vectorlabs.com](mailto: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 \(Anti-Mouse IgG-brown, Anti-Rabbit IgG-magenta\)](https://vectorlabs.com/products/immpress-duet-double-staining-kit-red/)

<https://vectorlabs.com/products/immpress-duet-double-staining-kit-red/>