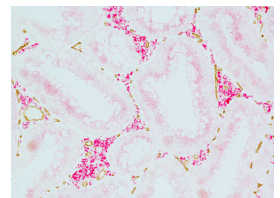
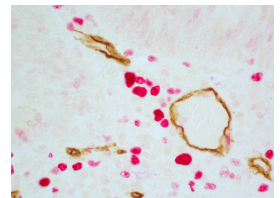
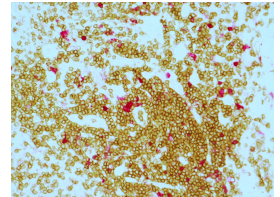
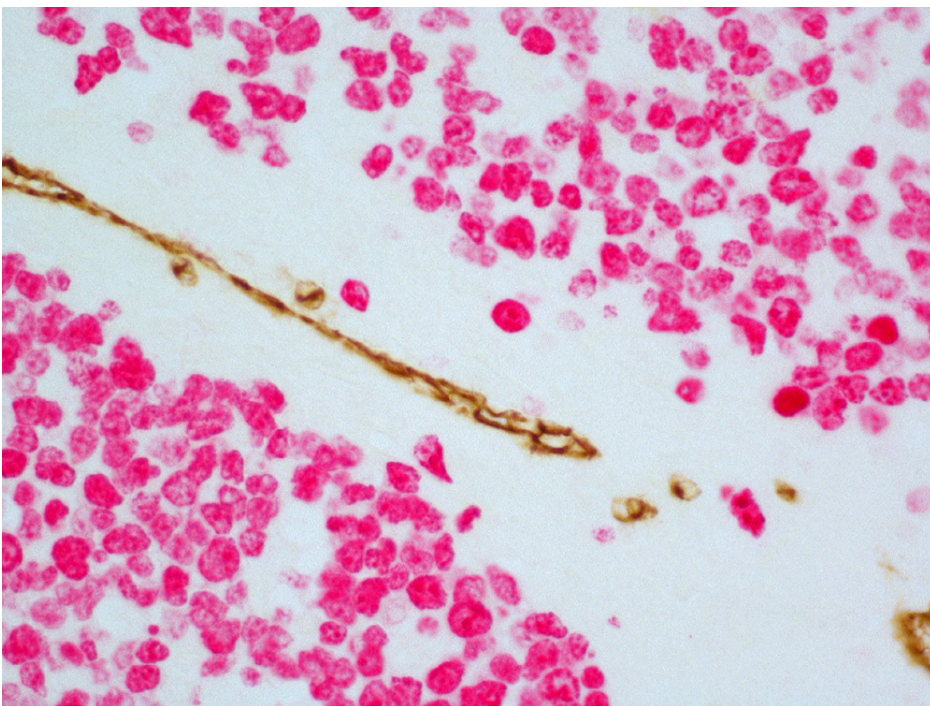


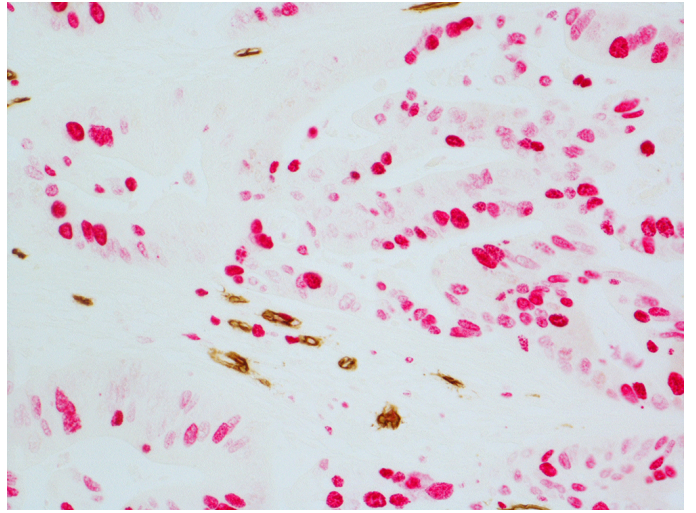
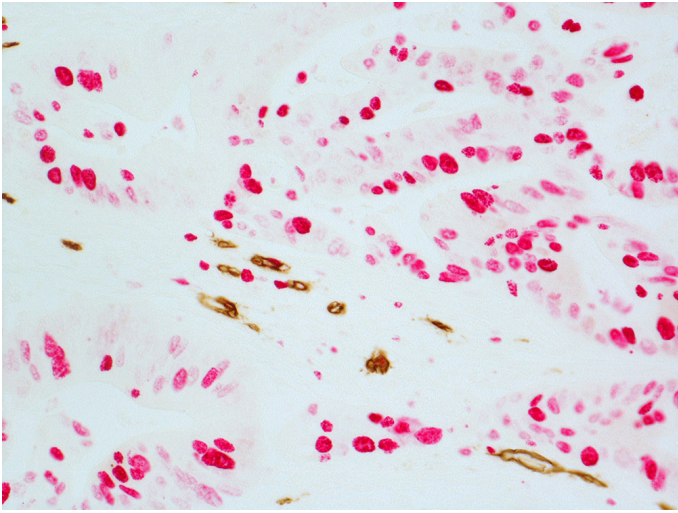


# ImmPRESS® Duet Double Staining Polymer Kit (HRP Anti-Mouse IgG-brown, AP Anti-Rabbit IgG-magenta)

## MP-7724

[Product Images](#)





## Short 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
- High sensitivity and Specificity
- Non-Biotin Based Reagents
- Prediluted, Ready-to-use

### Kit Contents:

- 15 ml BLOXALL<sup>®</sup> 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<sup>®</sup> DAB EqV Substrate (HRP, brown)
- ImmPACT Vector<sup>®</sup> Red Substrate (AP, magenta)

**Note: The ImmPRESS Duet Double Staining Polymer Kit is intended for use on non-rodent tissue specimens.**

## Additional Information

---

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

