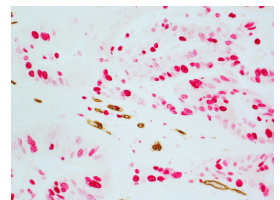
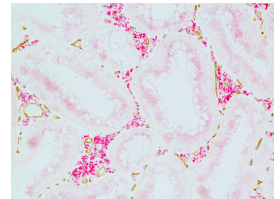
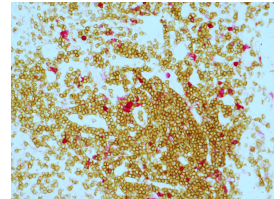
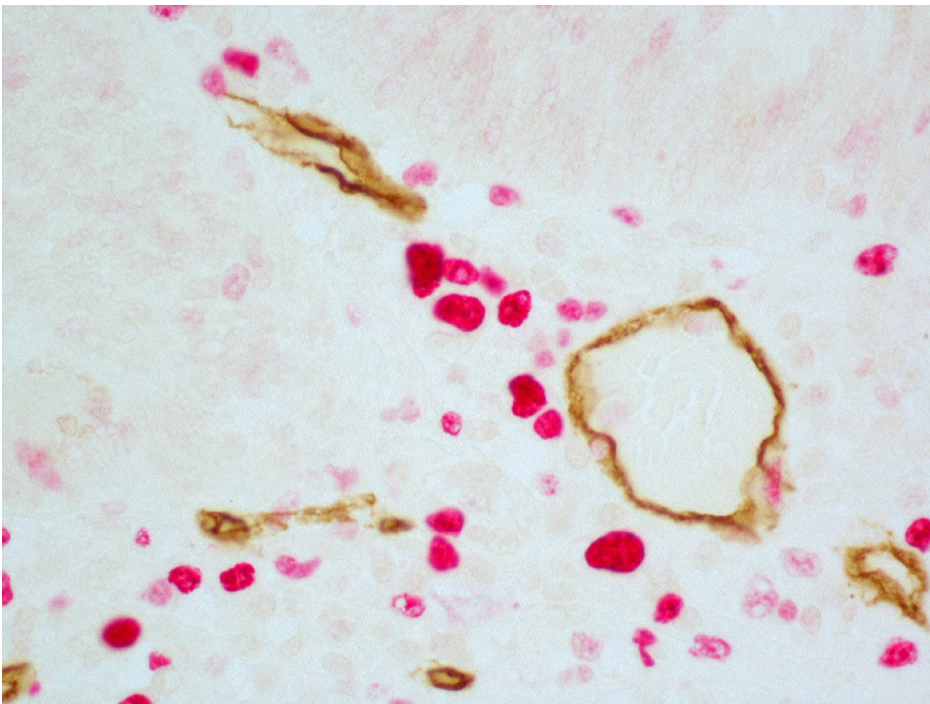


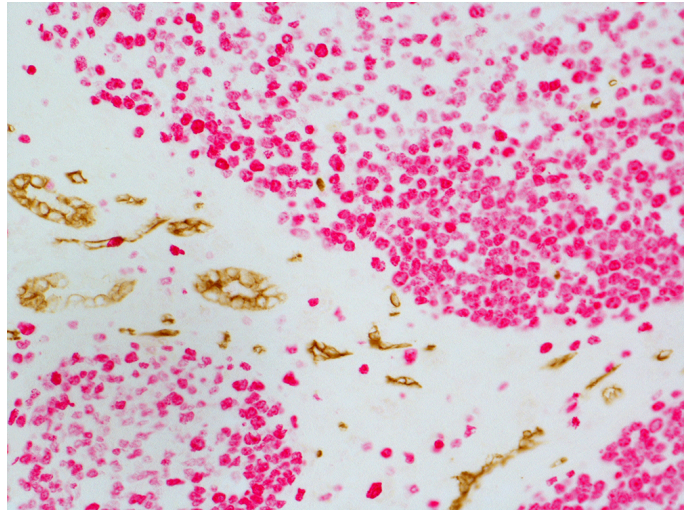
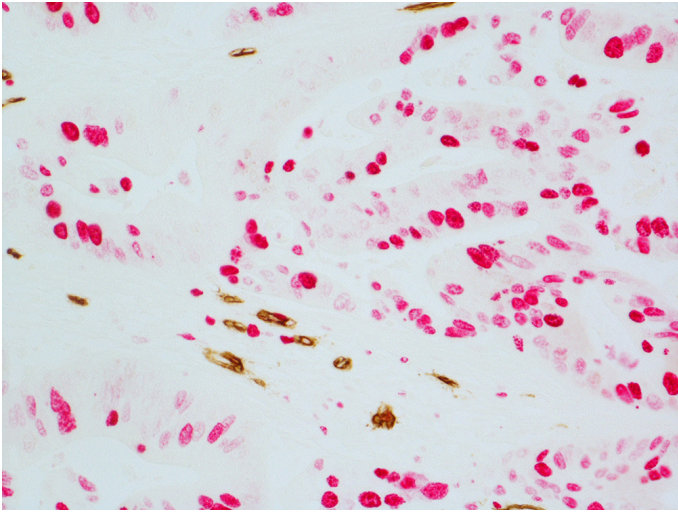


ImmPRESS® Duet Double Staining Polymer Kit (HRP – Anti-Rabbit IgG-brown, AP – Anti-Mouse IgG-magenta)

MP-7714

[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-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
- 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.

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

