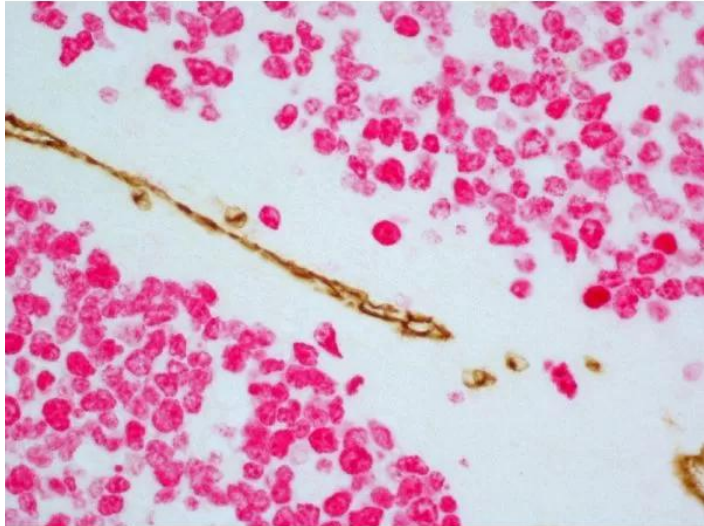


IMPRESS® DUET DOUBLE STAINING POLYMER KIT (HRP ANTI-MOUSE IGG-BROWN, AP ANTI-RABBIT IGG- MAGENTA)

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
- High sensitivity and Specificity
- Non-Biotin Based Reagents

For research use only. Not intended for animal or human therapeutic or diagnostic use.

- 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)

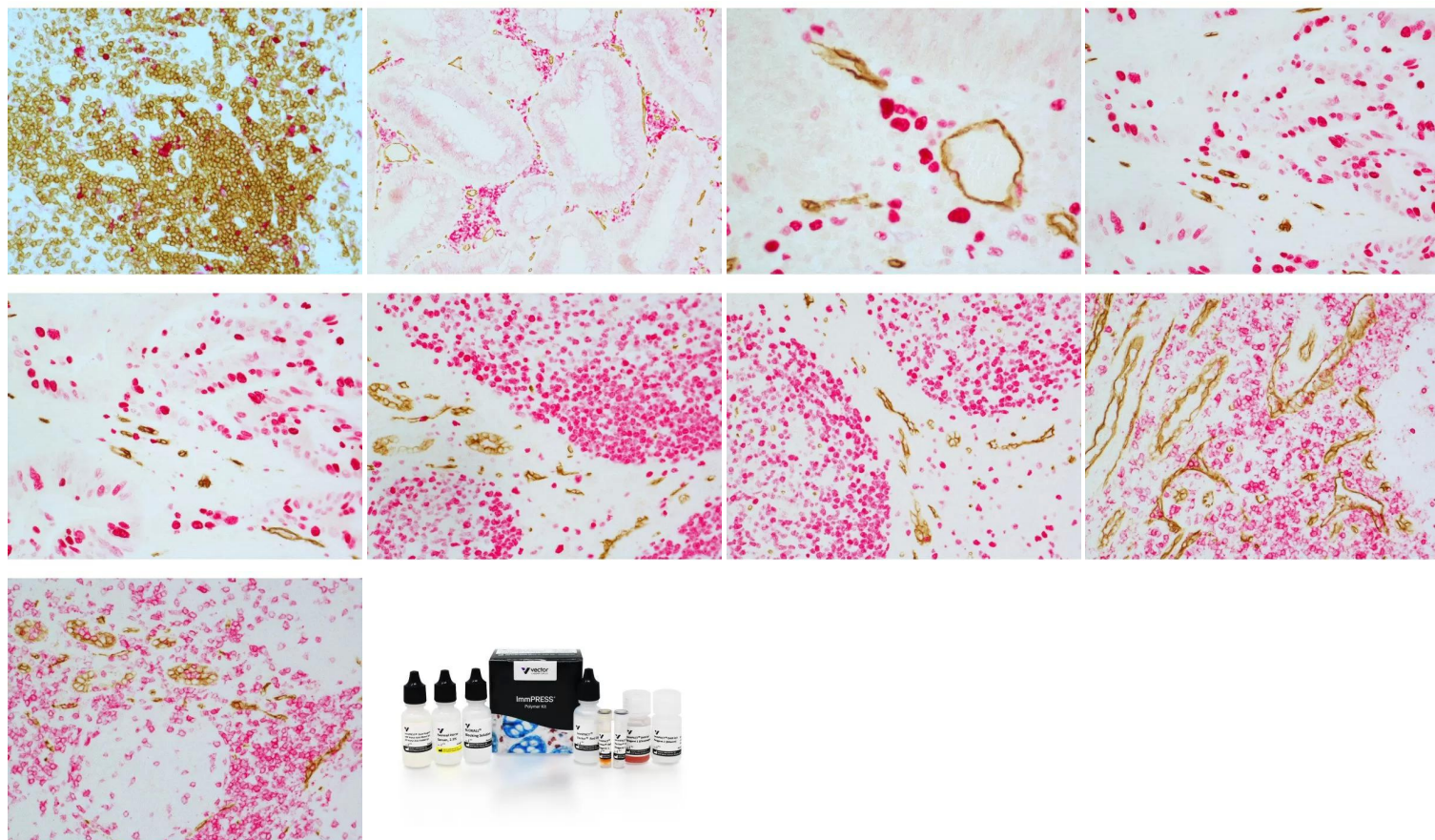
Note: The ImmPRESS Duet Double Staining Polymer Kit is intended for use on non-rodent tissue specimens.

Specifications

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

GALLERY IMAGES

For research use only. Not intended for animal or human therapeutic or diagnostic use.



For research use only. Not intended for animal or human therapeutic or diagnostic use.