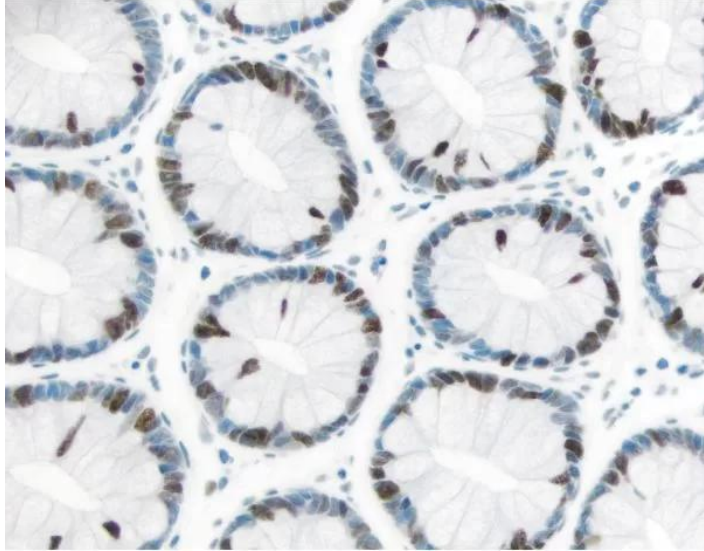




IMPRESS EXCEL AMPLIFIED POLYMER STAINING KIT, ANTI-RABBIT IGG, PEROXIDASE

SKU: MP-7601



DESCRIPTION

The ImmPRESS Excel Amplified HRP (peroxidase) Polymer Staining Kit is an enzymatic, amplification system that produces crisp, highly sensitive, specific staining with low background. It contains an amplifier antibody in a two-step method.

Features:

- Complete staining kits
- Precise, sharp antigen localization
- Convenient prediluted, ready-to-use reagents

Kit Contents:

- BLOXALL® Endogenous HRP/AP Blocking Solution (r.t.u.)
- 2.5% Normal serum for blocking non-specific binding (r.t.u.)

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



- Amplifier Antibody (either goat anti-mouse IgG or goat anti-rabbit IgG)
- ImmPRESS Excel Amplified HRP Polymer Reagent (anti-goat IgG)
- ImmPACT® DAB EqV Substrate/ImmPRESS Excel Staining Kit (15 ml) will stain approximately 75-150 sections and an ImmPRESS Excel Staining Kit (50 ml) will stain approximately 250-500 sections.

SPECIFICATIONS

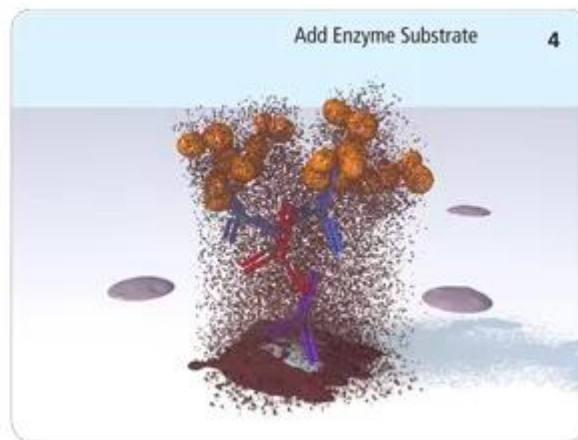
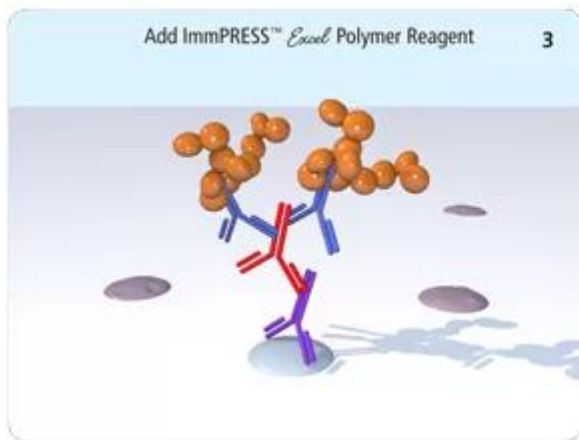
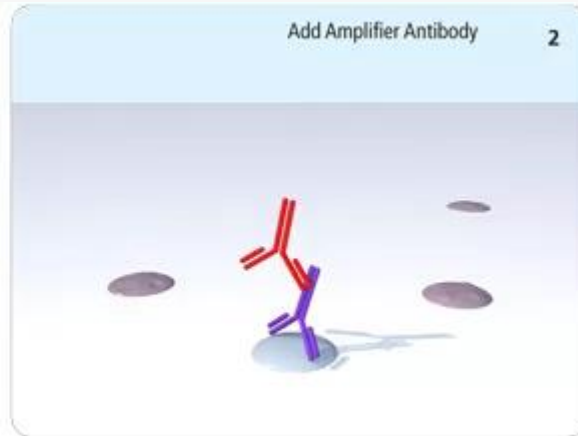
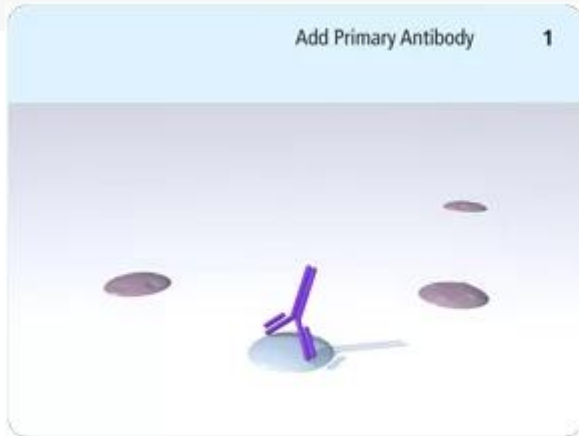
Detection Enzyme	Peroxidase
Format	Ready-to-Use
Isolation Technology	Amplified Micropolymer Systems
Unit Size	1 kit (15 or 50 ml)
Applications	ELISAs, Immunohistochemistry / Immunocytochemistry, In situ hybridization
Target Species	Rabbit
Conjugate	Micropolymer HRP
Platform	Amplified Micropolymer Systems
Source Species	Rabbit

TECHNICAL INFORMATION

The ImmPRESS Excel Amplified Staining System capitalizes on all the advantages of the ImmPRESS polymerized reporter enzyme staining system technology but offers additional sensitivity and convenience. This system employs a proprietary ready-to-use (RTU) Amplifier Antibody, followed by a ready-to-use (RTU) ImmPRESS Excel HRP Polymer Reagent. These reagents are affinity purified and extensively cross-adsorbed to ensure high sensitivity and low background. In addition, the ImmPRESS Excel Reagent is conjugated to horseradish peroxidase micropolymers, a novel approach that avoids the use of large dextrans or other macromolecules as a backbone. These unique micropolymers allow a higher density of enzymes per antibody to bind the target with minimal interference. The result is an increase in target accessibility, signal intensity, and significantly less background staining.

The included ImmPACT DAB EqV Substrate produces a crisp, dark brown reaction product with the excellent sensitivity and low background characteristic of the ImmPRESS /ImmPACT combination.

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



CITATIONS



Powered by Bioz © 2023 See more details on Bioz

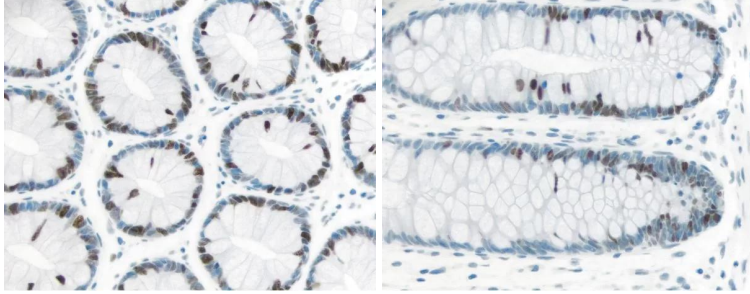
DOCUMENTS

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

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



GALLERY IMAGES



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