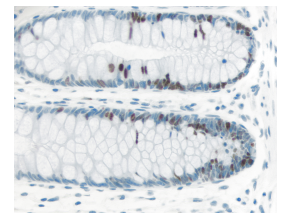
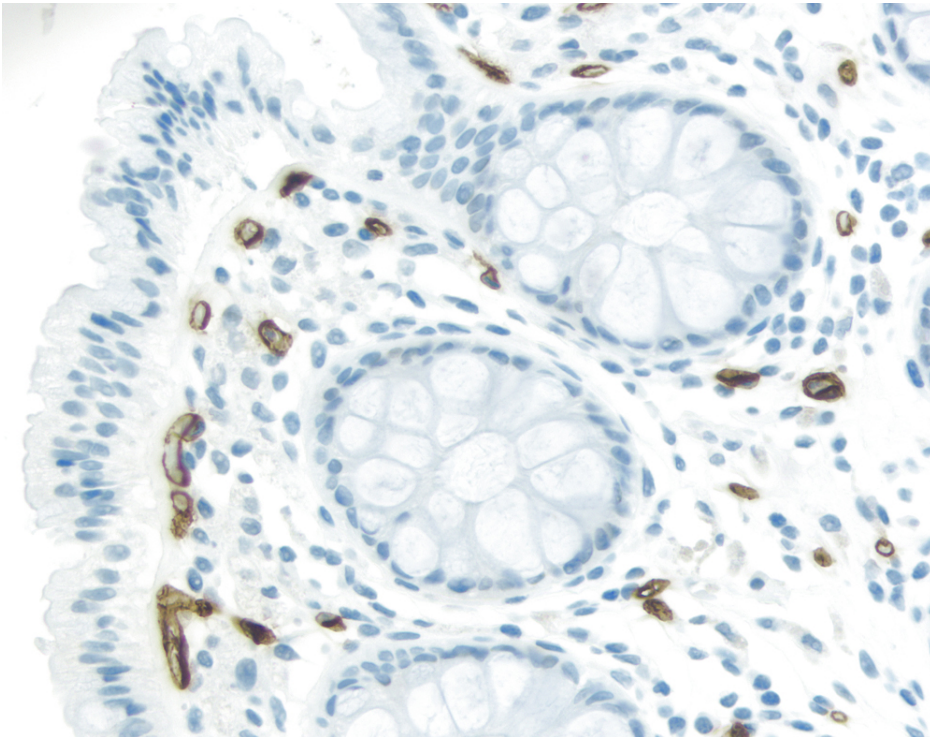




ImmPRESS® Excel Amplified Polymer Staining Kit, Anti-Mouse IgG, Peroxidase

MP-7602

[Product Images](#)



Short 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.)
- Amplifier Antibody (either goat anti-mouse IgG or goat anti-rabbit IgG)
- ImmPRESS Excel Amplified HRP Polymer Reagent (anti-goat IgG)
- ImmPACT[®] DAB EqV Reagent 1 (Chromogen)
- ImmPACT DAB EqV Reagent 2 (Buffer)

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.

Additional Information

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

ImmPRESS® Excel Amplified Polymer Staining Kit, Anti-Mouse IgG, Peroxidase

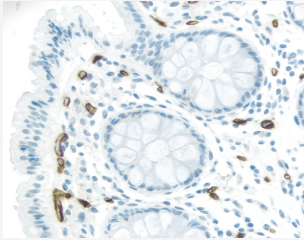
Features:

- Complete staining kits
- Highest sensitivity
- Precise, sharp antigen localization
- Low background
- Amplified polymer technology
- 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.)
- Amplifier Antibody (either goat anti-mouse IgG or goat anti-rabbit IgG)
- ImmPRESS Excel Amplified HRP Polymer Reagent (anti-goat IgG)
- ImmPACT® DAB EqV Reagent 1 (Chromogen)
- ImmPACT DAB EqV Reagent 2 (Buffer)

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.



ImmPRESS® Excel Amplified Polymer Staining Kit, Anti-Mouse IgG, Peroxidase

Features:

- Complete staining kits
- Highest sensitivity
- Precise, sharp antigen localization
- Low background
- Amplified polymer technology
- Convenient prediluted, ready-to-use reagents

Kit Contents:

- 15 ml BLOXALL® Endogenous HRP/AP Blocking Solution (r.t.u.)
- 15 ml 2.5% Normal serum for blocking non-specific binding (r.t.u.)
- 15 ml Amplifier Antibody (either goat anti-mouse IgG or goat anti-rabbit IgG)
- 15 ml ImmPRESS Excel Amplified HRP Polymer Reagent (anti-goat IgG)
- 10 ml ImmPACT® DAB EqV Reagent 1 (Chromogen)
- 10 ml ImmPACT DAB EqV Reagent 2 (Buffer)

