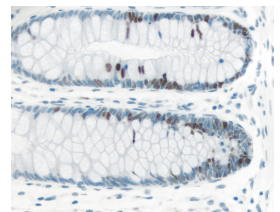
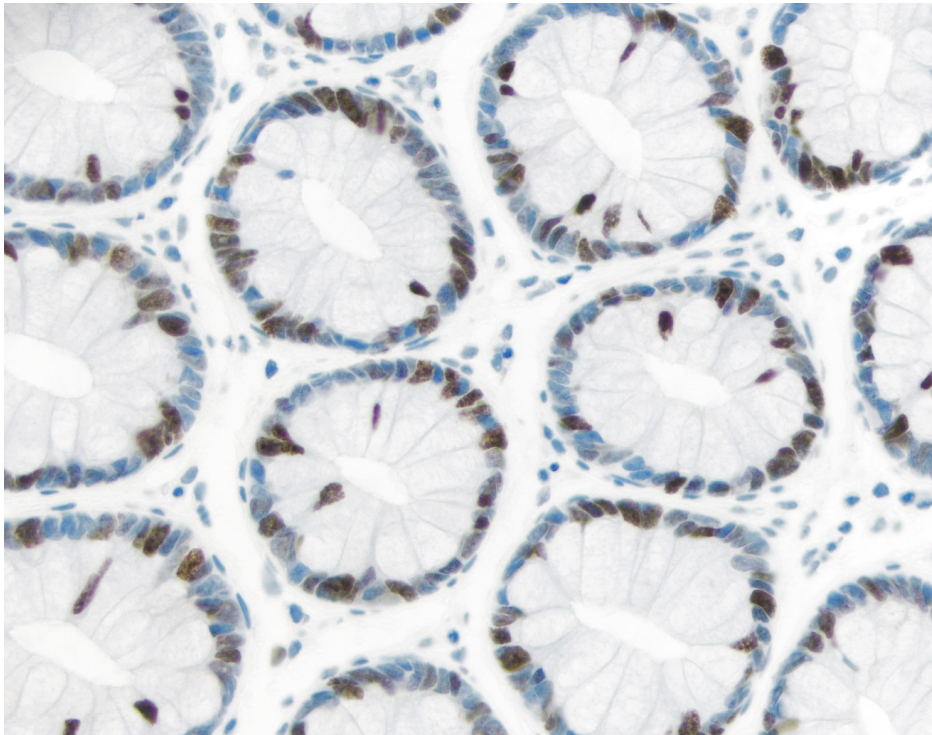




# ImmPRESS® Excel Amplified Polymer Staining Kit, Anti-Rabbit IgG, Peroxidase

## MP-7601

[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* 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.

## 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	Rabbit
Format	Ready-to-Use

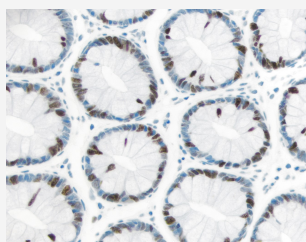
### ImmPRESS® Excel Amplified Polymer Staining Kit, Anti-Rabbit IgG, Peroxidase

#### 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)
- 20 ml ImmPACT® DAB *EqV* Substrate



### ImmPRESS® Excel Amplified Polymer Staining Kit, Anti-Rabbit 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)
- 20 ml ImmPACT® DAB *EqV* Substrate

