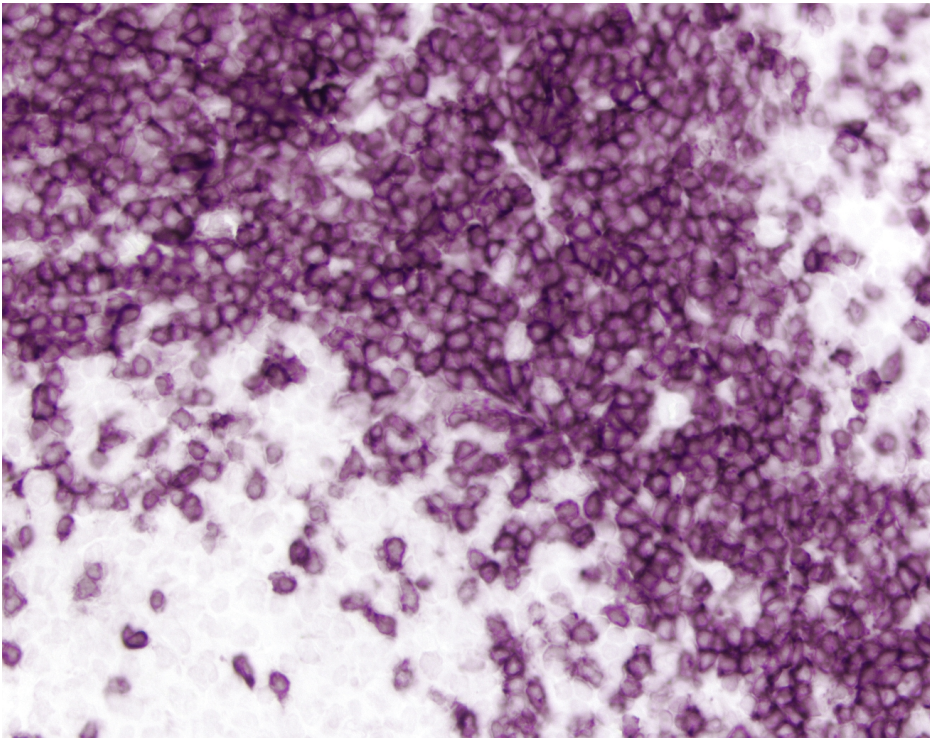




# ImmPRESS® HRP Goat Anti-Rat IgG Polymer Detection Kit, Peroxidase

## MP-7404-50

[Product Images](#)



## Short Description

---

The ImmPRESS polymerized reporter enzyme staining system uses novel conjugation and micropolymer chemistries to create a highly sensitive, ready-to-use, one-step detection system. This unique micropolymer of highly active enzyme is attached to our affinity purified secondary antibodies, producing reagents with outstanding sensitivity and low background.

### Features:

- Enhanced accessibility to nuclear and membrane antigens
- Ready-to-use in a convenient dropper bottle
- Shorter assay time
- Simplified multiple labeling
- Blocking solution included

The ImmPRESS (50 ml) Kit will stain approximately 250-500 sections.

## Additional Information

---

Unit Size	50 ml
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, ELISAs
Technology	Micropolymer Reagents
Detection Enzyme(s)	Peroxidase
Conjugate	Micropolymer HRP
Target Species	Rat
Format	Ready-to-Use

