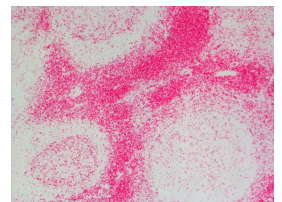
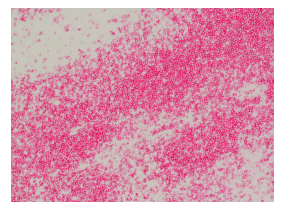
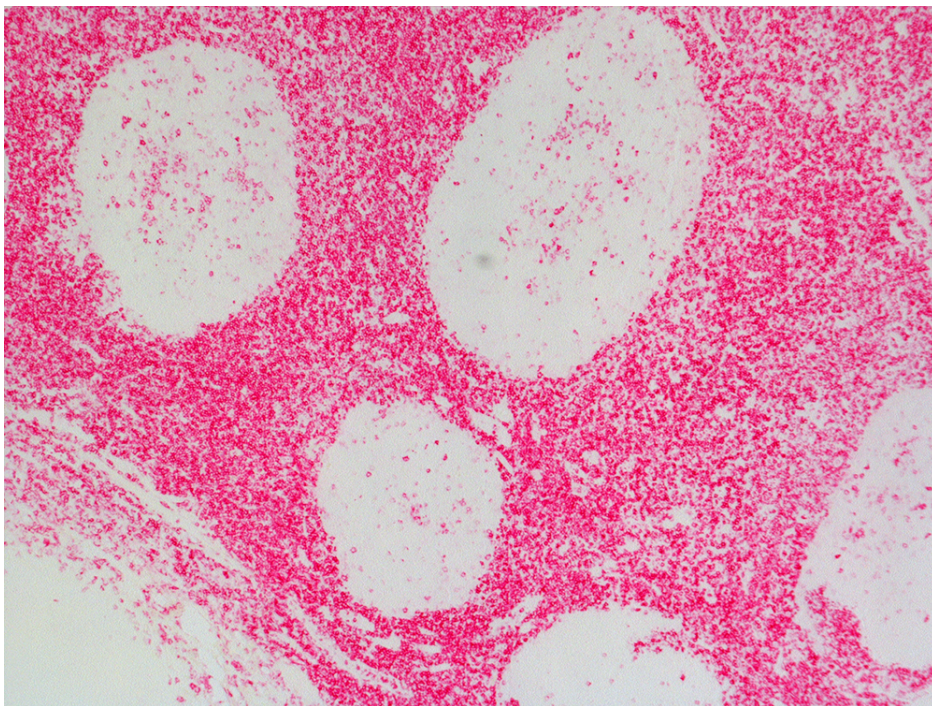




# ImmPRESS®-AP Goat Anti-Rat IgG Polymer Detection Kit, Alkaline Phosphatase

## MP-5404-15

[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, biotin-free detection system. This unique micropolymer of highly active enzyme (alkaline phosphatase) 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

The ImmPRESS (15 ml) Kit will stain approximately 75-150 sections.

## Additional Information

Unit Size	15 ml
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, ELISAs
Technology	Micropolymer Reagents
Detection Enzyme(s)	Alkaline Phosphatase
Conjugate	Micropolymer AP
Target Species	Rat
Format	Ready-to-Use

