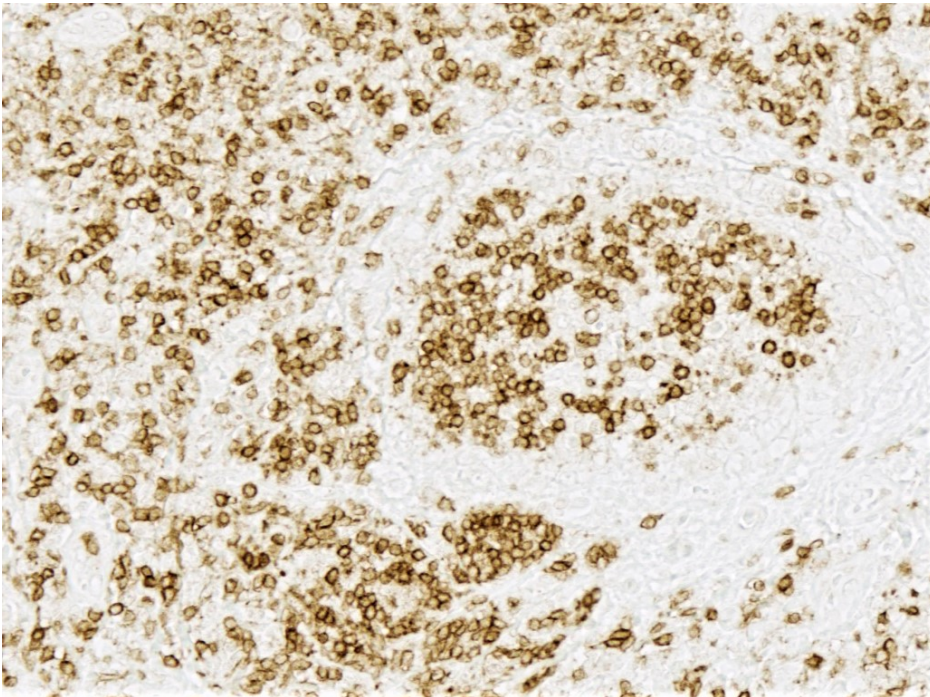




ImmPRESS® HRP Horse Anti-Rabbit IgG PLUS Polymer Kit, Peroxidase

MP-7801

[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 (peroxidase or alkaline phosphatase) is attached to our affinity purified secondary antibodies, producing reagents with outstanding sensitivity and low background. The inclusion of BLOXALL[®] endogenous enzyme quenching solution and ImmPACT[®] DAB EqV substrate in the PLUS kit provides a more complete detection system.

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	1 Kit
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, ELISAs
Technology	Micropolymer Reagents
Detection Enzyme(s)	Peroxidase
Conjugate	Micropolymer HRP
Target Species	Rabbit
Format	Ready-to-Use

