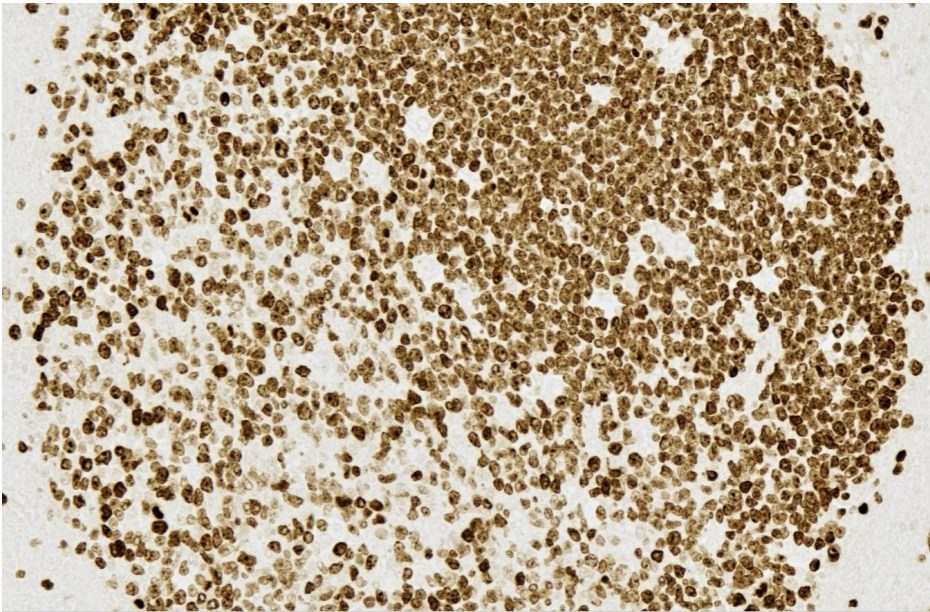




# ImmPRESS® HRP Horse Anti-Mouse IgG PLUS Polymer Kit, Peroxidase

## MP-7802

[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) is attached to our affinity purified secondary antibodies, producing reagents with outstanding sensitivity and low background. The inclusion of BLOXALL<sup>®</sup> endogenous enzyme quenching solution and ImmPACT<sup>®</sup> 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
Format	Ready-to-Use

