



ImmPRESS® HRP Universal (Horse Anti-Mouse/Rabbit IgG) PLUS Polymer Kit, Peroxidase

MP-7800

[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. The inclusion of BLOXALL[®] endogenous enzyme quenching solution and ImmPACT[®] DAB EqV substrate in the PLUS kit provides added convenience with optimized, matched reagents in one product.

Features:

- High sensitivity and low background
- Enhanced accessibility to nuclear and membrane antigens
- Ready-to-use in a convenient dropper bottle
- Shorter assay time
- Simplified multiple labeling
- Designed for use with both rabbit and mouse primary antibodies

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

Note: This polymer detection kit should not be used to stain tissues from rodents or rabbits due to its reactivity with IgG present in the tissues of these species.

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	Universal (Mouse/Rabbit)
Format	Ready-to-Use

