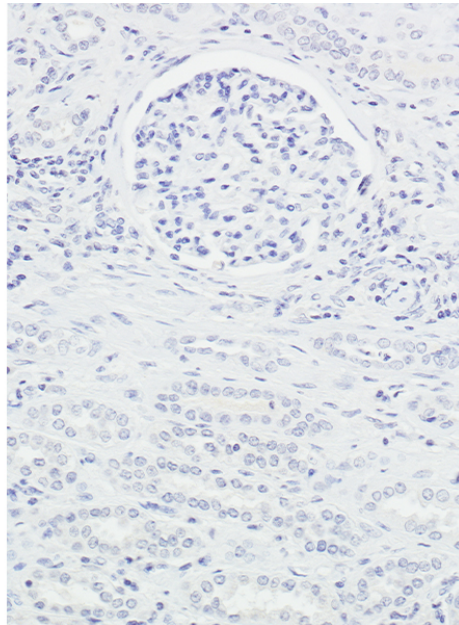
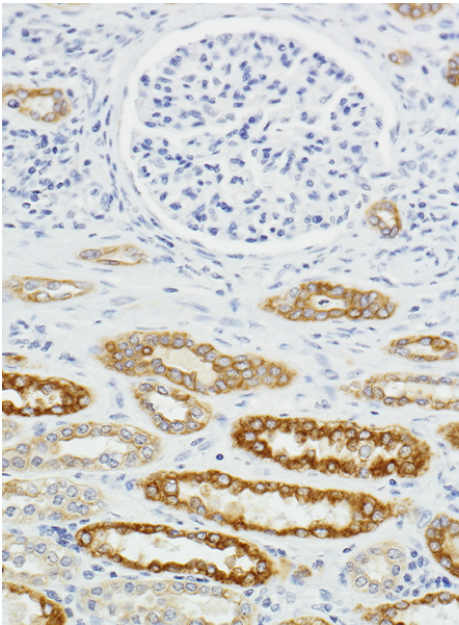




H.O.H.[™] (Human on Human) Immunodetection Kit

HOH-3000

Product Images



Short Description

The H.O.H. Immunodetection Kit is intended to detect human (or humanized) antibodies on frozen or paraffin embedded human tissue sections.

Key advantages of the H.O.H. Immunodetection Kit:

- Clear, crisp, specific staining of antigens of interest
- Procedures are simple and easy to follow
- No tedious calculations are required

The H.O.H. Immunodetection Kit contains enough stock reagents to produce about 5 ml of working solution, which is generally sufficient to stain approximately 50 - 100 tissue sections.

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Blotting Applications
Technology	Human on Human (H.O.H.) Systems
Detection Enzyme(s)	Peroxidase
Target Species	Human
Format	Concentrate
Blocking Action	Endogenous Human Ig

