

Demonstration of Immunohistochemical Staining Using Promega Anti-ACTIVE® and Apoptosis...

Demonstration of Immunohistochemical Staining Using Promega Anti-ACTIVE® and Apoptosis Antibodies

ABSTRACT

This article illustrates tissue staining with Promega polyclonal antibodies in both mouse and rat systems. We show immunohistochemical data for Anti-ACTIVE® MAPK, Anti-ACTIVE® JNK, Anti-ACTIVE® p38 and Anti-ACTIVE® CaM KII pAbs. In addition, we present work using polyclonal antibodies for the study of apoptosis (Anti-PARP p85 Fragment pAb, Anti-ACTIVE® Caspase-3 pAb, and Anti-pS473 Akt pAb). Protocols for immunohistochemical data are provided.

Promega Notes **80**, 20–23.

Randy Hoffman¹, Martha O'Brien¹, and Bruce W. Jarvis²

¹Promega Corporation, ²School of Veterinary Medicine, University of Wisconsin, Madison, WI

Publication Date: 2002



Download Article

(246 KB)
