

Technically Speaking: Anti-ACTIVE® Antibodies and MAPK Signaling Pathways

ABSTRACT

The mitogen-activated protein kinase (MAPK) superfamily of intracellular kinases is involved in many cell signaling and regulatory pathways. These kinases are activated through dual phosphorylation of the threonine and tyrosine residues in response to a multitude of extracellular stimuli.

1

Promega Notes 76, 23-24.

Michael Curtin
Promega Corporation
Publication Date: 2000



Download Article

(57 KB)