

Kinase Inhibitors and Activators

ABSTRACT

Promega has recently made important additions to its line of highly pure, quality-tested products for cell signaling research. Here we examine questions commonly posed by investigators considering the utility of these products in a variety of experimental applications.

Cell Notes 1, 10-11

James J. Cali
Promega Corporation
Publication Date: 2001



Download Article

(55 KB)