

In Vitro Toxicology and Cellular Fate Determination Using Promega Cell-Based Assays

ABSTRACT

This article describes options for measuring cellular fate by using reagents and systems developed to detect specific cellular markers and events. Here we show that a combination of Promega cell-based assays provides cytotoxicity researchers the tools needed to differentiate cell fates such as apoptosis and necrosis.

1

Cell Notes 4, 2-5.

Randy Hoffman
Promega Corporation
Publication Date: 2002



Download Article

(110 KB)