

BacTiter-Glo™ Assay for Antimicrobial Drug Discovery and General Microbiology

ABSTRACT

The BacTiter-Glo™ Microbial Cell Viability Assay determines the number of viable cells in culture by measuring ATP. This simple assay can provide results in as little as five minutes, and its excellent sensitivity allows you to detect microbial growth sooner than with conventional O.D. measurements. It is easily adaptable to automated and multiwell formats, and automated reagent injectors are not required.

Promega Notes 89, 25-27.

Frank Fan, Braeden Butler, Terry Riss and Keith Wood Promega Corporation

Publication Date: 2005



Download Article

(211 KB)