

Selecting Cell-Based Assays for Drug Discovery Screening

ABSTRACT

Promega produces a complete portfolio of homogeneous "add-mix-measure" assays. Choosing a cell-based screening method requires understanding the endpoint measured, the correlation with cell viability and the limitations of the assay chemistries. Here we provide recommendations for characterizing a model assay system, describe some of the factors to consider when choosing cell-based assays for automated systems and show examples of multiplexing cell-based assays.

Cell Notes **13**, 16–21.

Terry Riss, Rich Moravec and Andrew Niles

Promega Corporation

Publication Date: 2005



Download Article

(282 KB)
