



High-Throughput Screening Using a Universal Luminescent Kinase Assay

ABSTRACT

Phosphotransferases represent a major group of cellular enzymes that use ATP as a substrate and are implicated in a wide variety of cellular functions. The homogeneous Kinase-Glo® Luminescent Kinase Assay can monitor the activity of these enzymes, which have become important drug discovery targets. This non-radioactive assay is robust and easily amenable to high-throughput screening applications.

Cell Notes **10**, 20–23.

Said Goueli¹, Brad Larson¹, Kevin Hsiao¹, Tracy Worzella¹, Aoife Gallagher², and Eric Matthews³

¹Promega Corporation, ²Deerac Fluidics, ³BMG LABTECH

Publication Date: 2004



Download Article

(176 KB)
