

## Identify P-glycoprotein Substrates and Inhibitors with the Rapid HTS Pgp-Glo™ Assay System

## **ABSTRACT**

The  $Pgp-Glo^{TM}$  Assay Systems are luminescent P-glycoprotein (Pgp) ATPase assays. The  $Pgp-Glo^{TM}$  Assay detects the effects of compounds on recombinant human Pgp in a cell membrane fraction. The assay couples the ATPase activity of Pgp to the ATP dependence of the light-generating reaction of firefly luciferase. The rate of ATP consumption by Pgp is inversely correlated with the light output of the luciferase reaction. The  $Pgp-Glo^{TM}$  Assay is easily amenable to high-throughput screening and can complement cell-based Pgp studies.

Promega Notes 96, 11-14.

**Dongping Ma and James J. Cali**Promega Corporation
Publication Date: 2007



**Download Article** 

(254 KB)