

## Bright-Glo™ and Steady-Glo® Luciferase Assay Systems: Assay Reagents for Academic and...

Bright-Glo™ and Steady-Glo® Luciferase Assay Systems: Assay Reagents for Academic and Industrial Applications

## **ABSTRACT**

The Bright-Glo™ and Steady-Glo® Luciferase Assay Systems have been designed to meet the needs of virtually any research application of the firefly luciferase reporter gene. These reagents have been optimized for homogeneous or nonhomogeneous assay formats, manual or automated sample processing, small-scale analysis to ultrahigh-throughput screening. Assays may be performed in a single step by adding the reagent into the culture medium containing cells expressing luciferase. The reagents provide exceptional assay precision with the highest sensitivity available. Characteristics of each reagent are further described to assist researchers in choosing the best reagent for their experiments.

Promega Notes 75, 3-6.

Erika Hawkins, Braeden Butler and Keith V. Wood Promega Corporation
Publication Date: 2000



**Download Article** 

(261 KB)