

# Dual-Glo<sup>®</sup> Luciferase Assay System: A Homogeneous Dual-Reporter System

---

## ABSTRACT

*Promega has developed a new, homogeneous reporter assay, the Dual-Glo<sup>®</sup> Luciferase Assay System, for monitoring both firefly and Renilla luciferases expressed in mammalian cells in 96- and 384-well plates. The new assay system integrates cell lysis and luminescence chemistry, and the luminescence kinetics have been extended over several hours. The homogeneous assay format allows dual-reporter assays to be more easily and rapidly performed by reducing sample processing requirements and eliminating the need for reagent injectors in luminometers.*

*Cell Notes 4, 14–15.*

**Erika Hawkins, Braeden Butler, Keith Wood, Michael O'Grady, Laurie Orr, and Michael Beck**

*Promega Corporation*

Publication Date: 2002



## Download Article

(106 KB)

---