

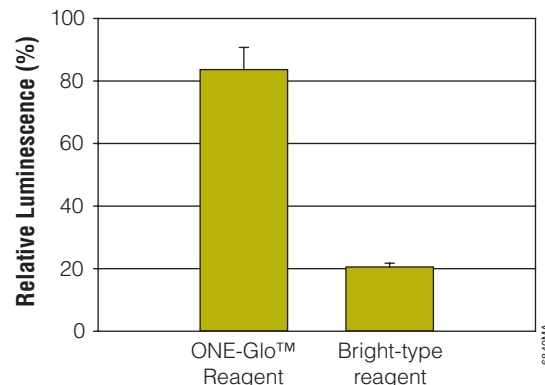
ONE-Glo™ Luciferase Assay System

Description

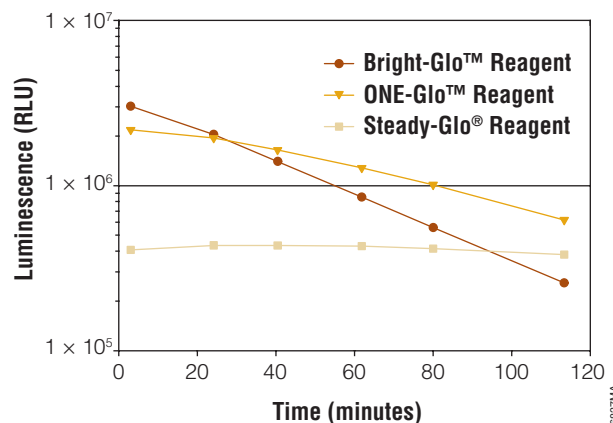
ONE-Glo™ Luciferase Assay Reagent provides a highly sensitive, robust, homogeneous assay for detection of firefly luciferase reporter gene expression in mammalian cells. Ideally suited for high- and ultrahigh-throughput applications, the ONE-Glo™ Assay contains a new luciferase substrate that is more stable, more tolerant to sample components, and has less odor than standard luciferase assay reagents. These features ensure that the ONE-Glo™ Reagent provides robust performance and also eliminates many of the handling inconveniences experienced using other reporter assays in a high-throughput setting.

Features

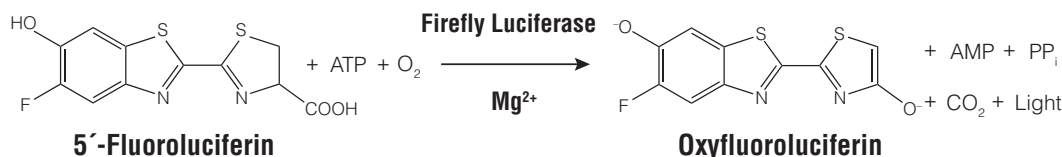
- **Simplify Your Assay Optimization:** The robust performance, reduced odor, improved storage, and larger sizes make assay optimization even easier and more efficient.
- **Store at Room Temperature or at 4°C:** The extended stability of the ONE-Glo™ Reagent at room temperature and 4°C makes it much more convenient for everyday use.
- **Improve Assay Precision:** Because the ONE-Glo™ Reagent is less sensitive to mixing and dispensing conditions, reproducibility is enhanced. The assay is ideal for use in high-density (384- and 1536-well) microplates.
- **Get a Brighter, Longer Signal:** Optimized for batch and continuous-process handling, the extended bright light output of the ONE-Glo™ Assay allows high sensitivity—especially when extended incubation is required prior to reading results.
- **Reduce Unwanted Effects from Sample Components:** The novel chemistry used in the ONE-Glo™ Assay is less sensitive to culture media, phenol red, and luciferase inhibitors than other luciferase assays.



ONE-Glo™ Reagent protects the luciferase reaction from inhibition by 10µM resveratrol, a known firefly luciferase inhibitor.

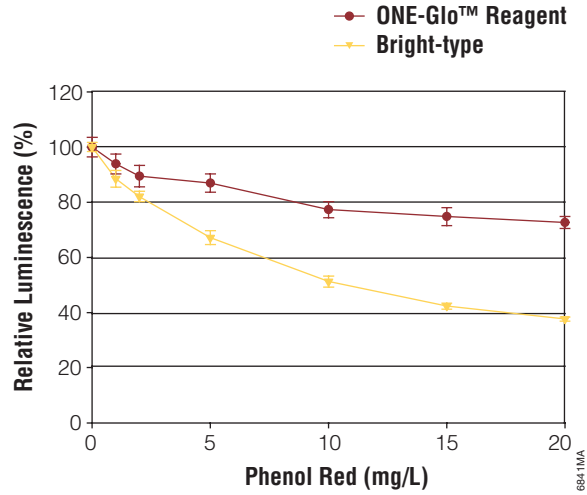


ONE-Glo™ Reagent generates bright and stable luminescence that is measurable for hours.

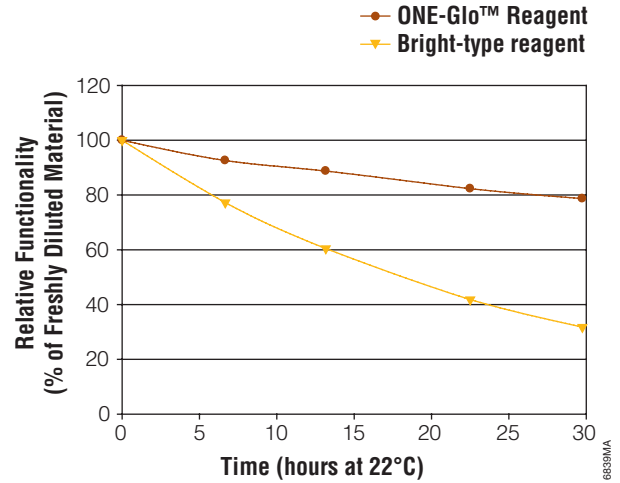


ONE-Glo™ uses a novel firefly luciferase substrate, 5'-fluoroluciferin.





ONE-Glo™ Reagent is more tolerant of phenol red than other luciferin-based reagents.



Reconstituted ONE-Glo™ Reagent can be used much longer than other Bright-type reagents.

Ordering Information

Product	Size	Cat. #
ONE-Glo™ Luciferase Assay System	10ml	E6110
	100ml	E6120
	1L	E6130

The method of recombinant expression of *Coleoptera* luciferase is covered by U.S. Pat. Nos. 5,583,024, 5,674,713 and 5,700,673.

Certain applications of this product may require licenses from others.

Products may be covered by pending or issued patents. Please visit www.promega.com for more information.



Promega Corporation • 2800 Woods Hollow Road • Madison, WI 53711-5399 USA • Telephone 608-274-4330 • Fax 608-277-2601

www.promega.com

©2008 Promega Corporation. All Rights Reserved.
Prices and specifications subject to change without prior notice.

Printed in USA 01/08
16143-DS-CR
Part #DS279

