

ONE-Glo™ EX Luciferase Assay System

Promega Corporation
2800 Woods Hollow Rd., Madison, WI 53711



This product uses less hazardous components and generates less waste to support sustainable practices in your lab.

Materials Affected

- ONE-Glo™ EX Luciferase Assay System, 10ml (Cat.# E8110)
- ONE-Glo™ EX Luciferase Assay System, 100ml (Cat.# E8120)
- ONE-Glo™ EX Luciferase Assay System, 10 × 10ml (Cat.# E8130)
- ONE-Glo™ EX Luciferase Assay System, 10 × 100ml (Cat.# E8150)

Introduction

We are committed to reducing the environmental effects from our operations globally, and with the products that we provide. We are working to reduce greenhouse gas emissions, energy, water, waste and effects from product distribution as part of our Corporate Responsibility Program. To learn more, visit: promega.com/responsibility

This green sheet discloses the basis of our environmental claims and highlights how the ONE-Glo™ EX Luciferase Assay System formulation reduces its environmental effects.

Product Description

ONE-Glo™ EX Luciferase Assay System is a highly sensitive assay for the detection of firefly luciferase with long-lived luminescence. Once added to cells, this assay provides a bright, stable signal that can be measured for hours. The reconstituted reagent has increased stability allowing for room temperature or 4°C storage, reduced sensitivity to phenol red in the culture media and elimination of odor-causing thiol compounds, such as DTT.

Sustainability Features

The formulation of the ONE-Glo™ EX Reagent has the following environmental benefits:

- **Less Waste:** Increased stability of the reconstituted reagent at room temperature or 4°C eliminates the need to dispense into aliquots and refreeze, simplifying repeated use over long experiments and reducing waste.
- **Energy Conservation:** Product stability at room temperature or 4°C reduces requirements for ultra-cold storage and saves energy.
- **Less Hazardous:** No thiols like DTT or other odor-causing compounds are used in the reagent, making transportation and handling nonhazardous.