

Detecting Toxicological Responses in Cells with the Bioluminescent GSH-Glo™ Glutathione...

Detecting Toxicological Responses in Cells with the Bioluminescent GSH-Glo™ Glutathione Assay

ABSTRACT

Here we present a new bioluminescent assay for determining the concentration of the reduced form of glutathione in cells or tissue lysates. The bioluminescent format reduces test compound interference and uses a protocol easily amenable to high-throughput screening applications.

Cell Notes **20**, 15–17.

Nancy Murphy, John Schultz, Wenhui Zhou and Keith V. Wood

Promega Corporation

Publication Date: 2008



Download Article

(269 KB)
