

The Bioluminescence Advantage

ABSTRACT

Although scientists may draw upon a range of photon-emitting chemistries to serve different experimental purposes, increasingly bioluminescence is becoming a preferred choice for many types of analyses. This article outlines the underlying reasons for the advantage bioluminescence brings to bioanalytical methodologies.

Promega Notes 96, 3–5.

Keith V. Wood

Promega Corporation

Publication Date: 2007



Download Article

(234 KB)
