

Six Luminescent Substrates for Cytochrome P450 Assays

ABSTRACT

Luminogenic CYP assays use prosubstrates for the light-generating reaction of firefly luciferase. CYPs convert the prosubstrates to an active luciferin, which produces light in a second reaction with luciferase. The amount of light generated is proportional to the amount of luciferin produced by the CYP and, therefore, to CYP enzyme activity.

Cell Notes 23, 19–20.

Promega Corporation

Publication Date: 2009



Download Article

(65 KB)
