

GloMax[®] 96 Microplate Luminometer Method for the Caspase-Glo[®] 3/7 Assay

ABSTRACT

The GloMax[®] 96 Microplate Luminometer in combination with the Caspase-Glo[®] Assay provides a convenient procedure for measuring caspase-3 or -7 activities. In the Caspase-Glo[®] Assay a proluminescent caspase-3/7 substrate contains the signature DEVD peptide sequence. Cell lysis occurs following the addition of Caspase-Glo[®] 3/7 Reagent. Lysed cells release caspase-3 or -7, which cleave the DEVD substrate from the aminoluciferin. Finally, the oxidation of the luciferin by luciferase produces light, which is measured by the GloMax[®] 96 plate reader. The amount of light produced is proportional to the amount of caspase-3 or -7 activity.



Download Article

(113 KB)
