

GloMax[®] 20/20 Luminometer Method for ENLITEN[®] ATP Assay System Bioluminescence...

GloMax[®] 20/20 Luminometer Method for ENLITEN[®] ATP Assay System Bioluminescence Detection Kit for ATP

ABSTRACT

The GloMax[®] 20/20 Luminometer combined with the ENLITEN[®] ATP Assay System Bioluminescence Detection Kit for ATP provides a sensitive, rapid method for measuring adenosine 5'-triphosphate (ATP). A reliable method for ATP detection is useful for studying enzymes that produce or degrade ATP. ATP detection also provides an indirect measurement of microbes, food residue, or other biological material.



Download Article

(121 KB)
