

Assessing Viability in Primary Cultures Using a Luminescent ATP-Based Assay:...

Assessing Viability in Primary Cultures Using a Luminescent ATP-Based Assay: Applications for the CellTiter-Glo® Assay

ABSTRACT

Here we highlight previously unpublished data and several peer-reviewed publications in which primary cell cultures were studied using the CellTiter-Glo® Assay. The sensitivity of the CellTiter-Glo® Assay allows researchers to conserve sample when working with extremely precious primary cell cultures.

Cell Notes **19**, 14–15.

Publication Date: 2007



Download Article

(98 KB)
