

# Use of pluripotent stem cell-derived cardiomyocytes to understand mechanisms of...

Use of pluripotent stem cell-derived cardiomyocytes to understand mechanisms of cardiotoxic compounds

---

## ABSTRACT

*Human pluripotent stem cell technology has the potential to overcome inefficiencies in early drug screening and clinical testing by allowing researchers to use human cells that reflect a variety of genetic backgrounds in assays that require minimal sample and test compound. Here we provide information about current practices and brief examples of the use of pluripotent stem cells in cytotoxicity testing using assays to detect mechanisms of cytotoxicity, such as apoptosis.*

*Cell Notes* **23**, 10–13.

**Jia-Quang He, Junyi Ma, and Blake Anson**

*Cellular Dynamics International, Inc.*

Publication Date: 2009



## Download Article

(153 KB)

---

## HOW TO CITE THIS ARTICLE

Scientific Style and Format, 7th edition, 2006

He, J., Ma, J. and Anson, B. Use of pluripotent stem cell-derived cardiomyocytes to understand mechanisms of cardiotoxic compounds. [Internet] 2009. [cited: year, month, date]. Available from: <http://www.promega.com/resources/pubhub/cellnotes/use-of-pluripotent-stem-cell-derived-cardiomyocytes-to-understand-mechanisms-of-cardiotoxic-compounds/>

American Medical Association, Manual of Style, 10th edition, 2007

He, J., Ma, J. and Anson, B. Use of pluripotent stem cell-derived cardiomyocytes to understand mechanisms of cardiotoxic compounds. Promega Corporation Web site. <http://www.promega.com/resources/pubhub/cellnotes/use-of-pluripotent-stem-cell-derived-cardiomyocytes-to-understand-mechanisms-of-cardiotoxic-compounds/> Updated 2009. Accessed Month Day, Year.