

References

1. Conti, E. *et al.* (1996) *Structure* **4**, 287–298.
2. Croch, S.P.M. *et al.* (1993) *J. Immunol. Meth.* **160**, 81–88.
3. Bliss, T.V. and Collingridge, G.L. (1993) *Nature* **361**, 31–39.
4. Choi, D.W. *et al.* (1992) *J. Neurosci.* **7**, 357–368.
5. Obrenovitch, T.P. and Urenjak, J. (1997) *Progress in Neurobiol.* **51**, 39–87.
6. Rothstein, J.D. (1995) *Advances in Neurobiol.* **68**, 7–20.
7. Itano, Y. and Nomura, Y. (1995) *Brain Res.* **704**, 240–245.
8. Fall, C.P. and Bennett, J., Jr. (1999) *J. Neurosci. Res.* **55**, 620–628.
9. Banker, G. and Goslin, K., eds. (1998) *Culturing Nerve Cells*, 2nd ed. The MIT Press, Cambridge, Massachusetts.
10. Ray, J. and Gage, F.H. (1999) *In: Modern Techniques in Neuroscience Research*. Springer-Verlag.
11. Brewere, G.J. *et al.* (1993) *J. Neurosci. Res.* **35**, 567–576.
12. Los, G.V. *et al.* (1995) *Biochem. J.* **311**, 225–232.

Ordering Information

Product	Size	Cat. #
CellTiter-Glo™ Luminescent Cell Viability Assay ^(a)	10 × 100ml	G7573
	100ml	G7572
	10 × 10ml	G7571
	10ml	G7570

^(a)Patent Pending.

CellTiter-Glo is a trademark of Promega Corporation. AlexaFluor is a trademark of Molecular Probes. Axiovert is a registered trademark of Carl Zeiss, Inc. Hibernate and Neurobasal are trademarks of Life Technologies. Image-Pro is a registered trademark of Media Cybernetics, Inc. MLX is a registered trademark of DYNEX Technologies. Spot is a trademark of Diagnostic Instruments, Inc.

PROMEGA SCIENTISTS TO PRESENT TALKS AT AMERICAN SOCIETY FOR CELL BIOLOGY

A Fast and Sensitive “Single Addition” Luminescent Cell Viability Assay

Talk presented by Richard Moravec, M.S.
 Research and Development, Promega Corporation
 Sunday, December 9, 2001, 11:00am

We have developed a fast and sensitive, single-addition luminescent cell viability assay, the CellTiter-Glo™ Luminescent Cell Viability Assay^(a), which is based on the detection of ATP. The assay is performed by adding a single reagent directly to cells in culture. After mixing and incubating 10 minutes at ambient temperature, luminescence is measured with a luminometer or CCD imaging device. The luminescent signal is directly proportional to the number of viable cells, consistent with previously published reports using ATP as an indicator of cell viability. The luminescent signal produced has an extended half-life typically greater than 5 hours. The assay sensitivity is sufficient to detect 4 cells/well in a 384 well format and has a linear range of 3–4 logs, depending on cell type and plate format. We will present data demonstrating the use of the CellTiter-Glo™ Assay for cytotoxicity testing in a 384 well format and its compatibility with a variety of cell lines, media and solvents used to deliver drugs.

The Apo-ONE™ Homogeneous Caspase-3/7 Assay: The “Number ONE” Solution for Apoptosis Detection

Talk presented by Michael Curtin, B.S.
 Cell Signaling and Neuroscience, Promega Corporation
 Sunday December 9, 2001, 12:00pm

Promega Corporation has developed a one-step, highly sensitive, simple and robust assay to measure caspase-3 and caspase-7 activity. The Apo-ONE™ Homogeneous Caspase-3/7 Assay^(b) uses a rhodamine 110-based substrate in a specially formulated lysis buffer. This buffer, which is added directly to cells in culture, allows the user to quickly and accurately measure the activities of caspase-3 and caspase-7. We will present data comparing the Apo-ONE™ Assay to other currently existing technologies. Particular regard will be given to the assay’s use with primary, suspension and adherent cell lines in both apoptotic induction and inhibition screens as well as in potency testing. Furthermore, the inherent flexibility of the assay will be demonstrated from data collected on the Biomek® (Beckman) 2000 automated module. These demonstrations will show how quickly and routinely the assay can be used as an early and accurate indicator of apoptotic processes.

^(a)Patent Pending.

^(b)This product is covered by U.S. Pat. Nos. 4,557,862 and 4,640,893 and is sold for research use only. All other uses, including but not limited to use as a clinical diagnostic or therapeutic, require a separate license. Please contact Promega Corporation for details relating to obtaining a license for such other use.

Apo-ONE and CellTiter-Glo are trademarks of Promega Corporation. Biomek is a registered trademark of Beckman Instruments, Inc.