

Example Experiment

Measuring Apoptosis in HeLa Cells Following Anisomycin Treatment



Published September 2012. Revised February 2020.
The most recent version of this document is posted at
[licor.com/bio/support](https://www.licor.com/bio/support).

Visit us on protocols.io! Explore an interactive
version of this protocol at bit.ly/2Qxb5al.

Table of Contents

	Page
I. Background	3
II. Introduction	3
III. Experimental Design	3
IV. Required Reagents	4
LI-COR Reagents	4
Additional Reagents	4
V. Prepare Cells	5
VI. Treat Cells	5
VII. Fix and Permeabilize Cells	6
Fix Cells	6
Permeabilize Cells	6
VIII. Block Cells	6
IX. Primary Antibody	7
Dilute Primary Antibody	7
Incubate with Primary Antibody	7
Wash	7
X. Secondary Antibody	7
Dilute Secondary Antibody	7
Incubate with Secondary Antibody	8
Wash	8
XI. Image	9
Suggested Scan Settings	9
XII. Experimental Results	10
XIII. References	11

I. Background

Accurate quantification of protein expression and/or post-translational modifications is important for advancing both basic and translational research. Given the flexibility, reproducibility, and higher throughput of the In-Cell Western™ Assay, it offers a convenient alternative to Western blotting and is a powerful platform for meaningful *in situ* analyses. The In-Cell Western microplate format can be used to analyze:

- Protein phosphorylation and signaling ([1](#) - [3](#))
- Off-target effects of drugs on signaling pathways ([4](#))
- Timing and kinetics of signaling events ([5](#), [6](#))
- Quantification of viral load ([7](#) - [11](#))
- Genotoxicity assays ([12](#), [13](#))
- Cell proliferation and apoptosis assays ([14](#))
- Bacterial-induced epithelial signaling ([15](#))
- Glycoprotein analysis ([16](#), [17](#))
- Library screening ([18](#) - [20](#))
- Screening of monoclonal antibody clones ([21](#))

II. Introduction

Understanding the processes leading to programmed cell death (apoptosis) is of utmost importance in the study of many diseases, including autoimmune diseases, neurological diseases, and cancer ([22](#)). For example, damaged cells that do not enter the apoptotic pathway may continue to proliferate and become cancerous ([23](#)). Due to the importance and complexity of the apoptosis signaling pathway, it is not surprising that this pathway plays an important role in numerous diseases. Consequently, understanding how new therapies impact this complex pathway is of great importance to the drug discovery field.

III. Experimental Design

In this experiment, we illustrate how the In-Cell Western Assay can be utilized to quantify the apoptotic response after drug treatment in a human cancer cell line. Anisomycin, a protein synthesis inhibitor, has been shown to trigger apoptosis in several cancer cell lines ([24](#) – [30](#)). Caspases (cysteinylnyl aspartate-specific proteases) are a family of important signaling molecules with essential roles that are both organ and subtype dependent ([31](#)). Caspase-3 has been implicated as an “effector” caspase associated with the initiation of apoptosis,

and its activity is a common marker used in cellular assays to quantify activators and inhibitors of the death cascade (32).

In the following example, caspase-3 activation is monitored using an antibody that detects endogenous levels of the large fragment (17/19 kDa) of activated caspase-3. This antibody does not recognize full length caspase-3 or other cleaved caspases. Cleaved caspase-3 levels are then normalized against CellTag™ 700 Stain (licor.com/CellTag). Therefore, we have investigated the effects of anisomycin on apoptosis as measured by caspase-3 activation in HeLa cells by In-Cell Western™.

IV. Required Reagents

LI-COR Reagents

- IRDye® 800CW Goat anti-Rabbit Secondary Antibody (LI-COR P/N 926-32211)
- CellTag™ 700 Stain (LI-COR P/N 926-41090)
- Intercept® (PBS) Blocking Buffer (licor.com/intercept)

Additional Reagents

- 1X PBS wash buffer
- HeLa cells (ATCC® CCL-2™)
- Tissue culture reagents (serum, DMEM, trypsin, 1X PBS)
- 20% Tween® 20
- Anisomycin (Sigma®, P/N A9789)
- 0.25% Trypsin/0.53 mM EDTA Solution (Sigma, P/N T-3924)
- 37% formaldehyde
- 10% Triton® X-100
- Black-sided 96-well or 384-well microplates with clear well bottoms
- Sealing film
- Cleaved Caspase-3 (Asp175) Antibody (Cell Signaling Technologies, P/N 9661)

V. Prepare Cells

1. Allow HeLa cells to grow in a T75 flask using standard tissue culture procedures until ~80% confluency is achieved ($\sim 1.5 \times 10^7$ cells; DMEM, 10% FBS; Gibco®).
2. Remove growth media and wash cells with sterile 1X PBS (room temperature (RT)).
3. Add 5 mL Trypsin-EDTA (Sigma) and incubate 3-5 minutes at 37 °C to displace cells.
4. Neutralize displaced cells with culture media and pellet by centrifugation.
5. Remove supernatant and disrupt the cell pellet manually by hand-tapping the collection tube.

Note: To maintain cell integrity, do not pipet or vortex during pellet disruption.

6. Reconstitute cells in complete media so that 50,000 cells/mL is achieved.
7. Manually mix the cell suspension thoroughly.
8. Under sterile conditions, dispense 200 μ L of the cell suspension per well in a 96-well plate (10,000 cells plated per well).
9. Incubate cells at 37 °C with 5% CO₂ and monitor cell density until ~80% confluency is achieved.

VI. Treat Cells

1. Warm serum-free media (DMEM, Gibco) to 37 °C.
2. Remove cell culture media.
3. Add either serum-free media for resting cells (mock) or serum-free media containing dilution series (1:2) of Anisomycin ranging in concentration from 0.07 - 40 μ M. Add 100 μ L of resting or activation media per well.
4. Allow incubation at 37 °C with 5% CO₂ for 4 hours.

VII. Fix and Permeabilize Cells

Fix Cells

1. Prepare fresh Fixing Solution as follows:

1X PBS	45 mL
37% Formaldehyde	5 mL
<hr/>	
3.7% Formaldehyde	50 mL

2. When incubation period is complete, carefully remove activation media manually or by aspiration to avoid detaching the cells.
3. Using a multi-channel pipettor, add 150 μ L of fresh Fixing Solution (RT). Add the solution by pipetting down the sides of the wells carefully to avoid detaching the cells from the well bottom.
4. Allow incubation on bench top for 20 minutes at RT with no shaking.

Permeabilize Cells

1. Prepare Triton[®] Washing Solution as follows:

1X PBS	495 mL
10% Triton X-100	5 mL
<hr/>	
1X PBS + 0.1% Triton X-100	500 mL

2. Remove Fixing Solution to an appropriate waste container (contains formaldehyde).
3. Using a multi-channel pipettor, add 200 μ L Triton Washing Solution (RT). Add the solution down the sides of the wells carefully to avoid detaching the cells.
4. Allow plate to shake on a rotator for 5 minutes at RT.
5. Repeat washing steps 4 more times, removing wash manually each time.

Important: Do not allow cells to become dry during washing. Immediately add the next wash after manual disposal.

VIII. Block Cells

1. Using a multi-channel pipettor, add 150 μ L of Intercept[®] Blocking Buffer to each well. Add the solution by pipetting down the sides of the wells carefully to avoid detaching the cells.
2. Allow blocking for 1.5 hours at RT with moderate shaking on a plate shaker.

IX. Primary Antibody

Dilute Primary Antibody

1. Dilute the primary antibody in Intercept® Blocking Buffer. Cleaved Caspase-3; rabbit (1:100 dilution)
2. Mix the primary antibody solution thoroughly before adding to wells.

Incubate with Primary Antibody

1. Remove blocking buffer and add 50 µL of the desired primary antibody in Intercept Blocking Buffer to cover the bottom of each well.
2. Make sure to include control wells without primary antibody to serve as a source for background well intensity. Only add 50 µL of Intercept Blocking Buffer to control wells.
3. Securely seal plate with sealing film and incubate with primary antibody overnight with gentle shaking at 4° C.

Wash

1. Prepare Tween® Washing Solution as follows:

1X PBS	995 mL
20% Tween 20	5 mL
<hr/>	
1X PBS + 0.1% Tween 20	1000 mL

2. Remove primary antibody solution.
3. Using a multi-channel pipettor, add 200 µL Tween Washing Solution (RT). Add solution down the sides of the wells carefully to avoid detaching the cells from the well bottom.
4. Allow wash to shake on plate shaker for 5 minutes at RT.
5. Repeat washing steps 4 more times.

X. Secondary Antibody

Dilute Secondary Antibody

1. Dilute the fluorescently-labeled secondary antibody in Intercept® Blocking Buffer as specified below. The recommended antibody dilution range is 1:200 to 1:1,200. To lower background, directly add CellTag™ 700 Stain to the antibody dilution to obtain a 1:500

dilution of CellTag 700 Stain.
IRDye® 800CW Goat anti-Rabbit (1:800 dilution)
CellTag™ 700 Stain (1:500 dilution)

Note: Minimize exposure of the antibody vials to light.

2. Thoroughly mix the antibody-stain solution then add 50 µL of the solution to each well.

Incubate with Secondary Antibody

1. Incubate for 60 minutes with gentle shaking at RT. Protect plate from light during incubation.

Wash

1. Remove secondary antibody solution.
2. Using a multi-channel pipettor, add 200 µL of Tween Washing Solution at RT (see "Wash" on the previous page). Add solution down the sides of the wells carefully to avoid detaching the cells from the well bottom.
3. Allow wash to incubate on a plate shaker for 5 minutes at RT.
4. Repeat washing steps 4 more times. Protect plate from light during washing.

XI. Image

1. After final wash, remove wash solution completely from wells. Turn the plate upside down and tap or blot gently on paper towels to remove traces of wash buffer. For best results, scan plate immediately; plates may also be stored at 4 °C for several weeks (sealed and protected from light).
2. Before plate scanning, clean the bottom plate surface and the Odyssey® Imager scanning bed (if applicable) with moist, lint-free tissue to avoid any obstructions during scanning.
3. Scan plate with detection in both 700 and 800 nm channels.

Suggested Scan Settings

All settings may require adjustment for optimal data quality. Higher resolutions or scan qualities can be used, but the scan time will increase.

Instrument	Resolution	Scan Quality	Intensity Setting (700 nm)	Intensity Setting (800 nm)
Odyssey Classic	169 µM	lowest	5	5
Odyssey CLx	169 µM	lowest	Auto Mode	Auto Mode
Odyssey Sa	200 µM	lowest	7	7
Aerius™ Imager	200 µM	lowest	7	7

XII. Experimental Results

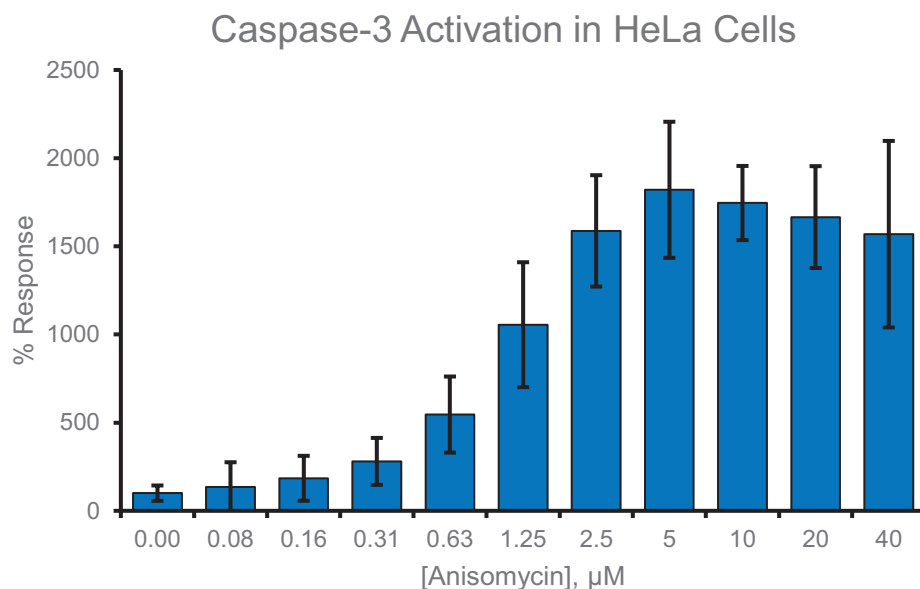


Figure 1. Induction of apoptosis in HeLa cells was achieved with increasing concentrations of anisomycin. An increase in cleaved caspase-3, a cleaved by-product indicative of apoptosis, is illustrated in the graph. Error bars represent the standard deviation from four technical replicates. The ultimate result of apoptosis induction is cell death. The reduction in cell number per well is taken into account when normalizing with CellTag™ 700 Stain. In an assay such as this, normalization is very important. See www.licor.com/ICWnormalization for more information on normalization options for an In-Cell Western™ Assay.

XIII. References

1. Chen, H., Kovar, J., Sissons, S., Cox, K., Matter, W., Chadwell, F., Luan, P., Vlahos, C. J., Schutz-Geschwender, A., and Olive, D. M. (2005) A cell-based immunocytochemical assay for monitoring kinase signaling pathways and drug efficacy. *Analytical biochemistry* 338, 136-142
2. Aguilar, H. N., Zielnik, B., Tracey, C. N., and Mitchell, B. F. (2010) Quantification of rapid Myosin regulatory light chain phosphorylation using high-throughput in-cell Western assays: comparison to Western immunoblots. *PLoS One* 5, e9965
3. Wong, S. K. (2004) A 384-well cell-based phospho-ERK assay for dopamine D2 and D3 receptors. *Analytical biochemistry* 333, 265-272
4. Kumar, N., Afeyan, R., Kim, H. D., and Lauffenburger, D. A. (2008) Multipathway model enables prediction of kinase inhibitor cross-talk effects on migration of Her2-overexpressing mammary epithelial cells. *Mol Pharmacol* 73, 1668-1678
5. Hannoush, R. N. (2008) Kinetics of Wnt-driven beta-catenin stabilization revealed by quantitative and temporal imaging. *PLoS One* 3, e3498
6. Chen, W. W., Schoeberl, B., Jasper, P. J., Niepel, M., Nielsen, U. B., Lauffenburger, D. A., and Sorger, P. K. (2009) Input-output behavior of ErbB signaling pathways as revealed by a mass action model trained against dynamic data. *Mol Syst Biol* 5, 239
7. Counihan, N. A., Daniel, L. M., Chojnacki, J., and Anderson, D. A. (2006) Infrared fluorescent immunofocus assay (IR-FIFA) for the quantitation of non-cytopathic and minimally cytopathic viruses. *J Virol Methods* 133, 62-69
8. Lin, Y. C., Li, J., Irwin, C. R., Jenkins, H., DeLange, L., and Evans, D. H. (2008) Vaccinia virus DNA ligase recruits cellular topoisomerase II to sites of viral replication and assembly. *J Virol* 82, 5922-5932
9. Weldon, S. K., Mischnick, S. L., Urlacher, T. M., and Ambroz, K. L. (2010) Quantitation of virus using laser-based scanning of near-infrared fluorophores replaces manual plate reading in a virus titration assay. *J Virol Methods* 168, 57-62
10. Lopez, T., Silva-Ayala, D., Lopez, S., and Arias, C. F. (2012) Methods suitable for high-throughput screening of siRNAs and other chemical compounds with the potential to inhibit rotavirus replication. *J Virol Methods* 179, 242-249
11. Wan, Y., Zhou, Z., Yang, Y., Wang, J., and Hung, T. (2010) Application of an In-Cell Western assay for measurement of influenza A virus replication. *J Virol Methods* 169, 359-364

12. Jamin, E. L., Riu, A., Douki, T., Debrauwer, L., Cravedi, J. P., Zalko, D., and Audebert, M. (2013) Combined genotoxic effects of a polycyclic aromatic hydrocarbon (B(a)P) and an heterocyclic amine (PhIP) in relation to colorectal carcinogenesis. *PLoS One* 8, e58591
13. Khoury, L., Zalko, D., and Audebert, M. (2013) Validation of high-throughput genotoxicity assay screening using gammaH2AX in-cell western assay on HepG2 cells. *Environ Mol Mutagen* 54, 737-746
14. Godin-Heymann, N., Ulkus, L., Brannigan, B. W., McDermott, U., Lamb, J., Maheswaran, S., Settleman, J., and Haber, D. A. (2008) The T790M "gatekeeper" mutation in EGFR mediates resistance to low concentrations of an irreversible EGFR inhibitor. *Mol Cancer Ther* 7, 874-879
15. Du, Y., Danjo, K., Robinson, P. A., and Crabtree, J. E. (2007) In-Cell Western analysis of Helicobacter pylori-induced phosphorylation of extracellular-signal related kinase via the transactivation of the epidermal growth factor receptor. *Microbes Infect* 9, 838-846
16. McInerney, M. P., Pan, Y., Short, J. L., and Nicolazzo, J. A. (2017) Development and Validation of an In-Cell Western for Quantifying P-Glycoprotein Expression in Human Brain Microvascular Endothelial (hCMEC/D3) Cells. *J Pharm Sci* 106, 2614-2624
17. Urlacher T, Xing K, Cheung L et al (2013) Glycoprotein applications using near-infrared detection. Poster presentation, Experimental Biology
18. Guo, K., Shelat, A. A., Guy, R. K., and Kastan, M. B. (2014) Development of a cell-based, high-throughput screening assay for ATM kinase inhibitors. *J Biomol Screen* 19, 538-546
19. Hoffman, G. R., Moerke, N. J., Hsia, M., Shamu, C. E., and Blenis, J. (2010) A high-throughput, cell-based screening method for siRNA and small molecule inhibitors of mTORC1 signaling using the In Cell Western technique. *Assay Drug Dev Technol* 8, 186-199
20. Schnaiter, S., Furst, B., Neu, J., Waczek, F., Orfi, L., Keri, G., Huber, L. A., and Wunderlich, W. (2014) Screening for MAPK modulators using an in-cell western assay. *Methods Mol Biol* 1120, 121-129
21. Daftarian, M. P., Vosoughi, A., and Lemmon, V. (2014) Gene-based vaccination and screening methods to develop monoclonal antibodies. *Methods Mol Biol* 1121, 337-346
22. Elmore, S. (2007) Apoptosis: a review of programmed cell death. *Toxicol Pathol* 35, 495-516
23. Hanahan, D., and Weinberg, R. A. (2000) The hallmarks of cancer. *Cell* 100, 57-70
24. Curtin, J. F., and Cotter, T. G. (2002) Anisomycin activates JNK and sensitises DU 145 prostate carcinoma cells to Fas mediated apoptosis. *Br J Cancer* 87, 1188-1194
25. Torocsik, B., and Szeberenyi, J. (2000) Anisomycin affects both pro- and antiapoptotic mechanisms in PC12 cells. *Biochem Biophys Res Commun* 278, 550-556

26. Kochi, S. K., and Collier, R. J. (1993) DNA fragmentation and cytolysis in U937 cells treated with diphtheria toxin or other inhibitors of protein synthesis. *Exp Cell Res* 208, 296-302
27. Kim, M., Lee, S. J., Shin, S., Park, K. S., Park, S. Y., and Lee, C. H. (2018) Novel natural killer cell-mediated cancer immunotherapeutic activity of anisomycin against hepatocellular carcinoma cells. *Sci Rep* 8, 10668
28. Cao, C., Yu, H., Wu, F., Qi, H., and He, J. (2017) Antibiotic anisomycin induces cell cycle arrest and apoptosis through inhibiting mitochondrial biogenesis in osteosarcoma. *J Bioenerg Biomembr* 49, 437-443
29. Chen, Q., Liu, X., Xu, L., Wang, Y., Wang, S., Li, Q., Huang, Y., and Liu, T. (2016) Long non-coding RNA BACE1-AS is a novel target for anisomycin-mediated suppression of ovarian cancer stem cell proliferation and invasion. *Oncol Rep* 35, 1916-1924
30. Liu, Y., Ge, J., Li, Q., Gu, L., Guo, X., Ma, Z. G., and Zhu, Y. P. (2013) Anisomycin induces apoptosis of glucocorticoid resistant acute lymphoblastic leukemia CEM-C1 cells via activation of mitogen-activated protein kinases p38 and JNK. *Neoplasma* 60, 101-110
31. Shalini, S., Dorstyn, L., Dawar, S., and Kumar, S. (2015) Old, new and emerging functions of caspases. *Cell Death Differ* 22, 526-539
32. Porter, A. G., and Janicke, R. U. (1999) Emerging roles of caspase-3 in apoptosis. *Cell Death Differ* 6, 99-104

© 2020 LI-COR, Inc. LI-COR, In-Cell Western, Aeries, Odyssey, Intercept, and IRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH

Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.

St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com