

NucView® References

[Click on a reference to link to the publication](#)

NucView® 405 in flow cytometry

Bohm, M., Wachtel, M., Marques, J. G., Streiff, N., Laubscher, D., Nanni, P., Mamchaoui, K., Santoro, R., and Schafer, B. W. (2016). [Helicase CHD4 is an epigenetic coregulator of PAX3-FOXO1 in alveolar rhabdomyosarcoma](#). *J Clin Invest* 126, 4237-4249. 10.1172/JCI85057 (Cell line: Rh4)

Grabocka, E., and Bar-Sagi, D. (2016). [Mutant KRAS Enhances Tumor Cell Fitness by Upregulating Stress Granules](#). *Cell* 167, 1803-1813 e1812. 10.1016/j.cell.2016.11.035 (Cell lines: DLD-1, NCI-H508)

NucView® 488 in flow cytometry

Amendola, R. S., Martin, A. C., Selistre-de-Araujo, H. S., Paula-Neto, H. A., Saldanha-Gama, R., and Barja-Fidalgo, C. (2015). [ADAM9 disintegrin domain activates human neutrophils through an autocrine circuit involving integrins and CXCR2](#). 10.1189/jlb.3A0914-455R (Cell type: primary human neutrophils)

Aronchik, I., Chen, T., Durkin, K. A., Horwitz, M. S., Preobrazhenskaya, M. N., Bjeldanes, L. F., and Firestone, G. L. (2011). [Target protein interactions of indole-3-carbinol and the highly potent derivative 1-Benzyl-I3C with the C-terminal domain of human elastase uncouples cell cycle arrest from apoptotic signaling](#). *Mol Carcinog*. DOI:10.1002/mc.20857 (Cell line: MDA-MB-231 (human breast adenocarcinoma))

Aronchik, I., Kundu, A., Quirrit, J. G., and Firestone, G. L. (2014). [The Anti-proliferative Response of Indole-3-carbinol in human melanoma cells is Triggered by an Interaction with NEDD4-1 and Disruption of Wild-type PTEN Degradation](#). *Mol Cancer Res*. 1541-7786. MCR-14-0018 (Cell line: G-361 (melanoma cells))

Ben Salem, I., Boussabbeh, M., Pires Da Silva, J., Guilbert, A., Bacha, H., Abid-Essefi, S., and Lemaire, C. (2017). [SIRT1 protects cardiac cells against apoptosis induced by zearalenone or its metabolites alpha- and beta-zearalenol through an autophagy-dependent pathway](#). *Toxicol Appl Pharmacol* 314, 82-90. 10.1016/j.taap.2016.11.012 (Cell line: H9c2)

Ben Salem, I., Prola, A., Boussabbeh, M., Guilbert, A., Bacha, H., Abid-Essefi, S., and Lemaire, C. (2015). [Crocic and Quercetin protect HCT116 and HEK293 cells from Zearalenone-induced apoptosis by reducing endoplasmic reticulum stress](#). *Cell Stress Chaperones* 20, 927-938. 10.1007/s12192-015-0613-0 (Cell lines: HCT116, HEK293)

Berenyi, E., Benko, I., Vamosi, G., Geresi, K., Tarkanyi, I., Szegedi, I., Lukacs, L., Juhasz, I., Kiss, C., Fesus, L., et al. (2011). [In vitro and in vivo activity of 4-thio-uridylylate against JY cells, a model for human acute lymphoid leukemia](#). *Biochem Biophys Res Commun* 410, 682-687. DOI:10.1016/j.bbrc.2011.06.056 (Cell line: JY (human lymphoblastic leukemia))

Boussabbeh, M., Ben Salem, I., Prola, A., Guilbert, A., Bacha, H., Abid-Essefi, S., and Lemaire, C. (2015). [Patulin induces apoptosis through ROS-mediated endoplasmic reticulum stress pathway](#). *Toxicol Sci* 144, 328-337. 10.1093/toxsci/kfu319 (Cell lines: HCT116, HEK293)

Boussabbeh, M., Prola, A., Ben Salem, I., Guilbert, A., Bacha, H., Lemaire, C., and Abis-Essefi, S. (2015). [Crocic and quercetin prevent PAT-induced apoptosis in mammalian cells: Involvement of ROS-mediated ER stress pathway](#). *Environ Toxicol*. 10.1002/tox.22185 (Cell lines: HCT116, HEK293)

Brodsky, I. E., and Medzhitov, R. (2008). [Reduced secretion of YopJ by Yersinia limits in vivo cell death but enhances bacterial virulence](#). *PLoS Pathog* 4, e1000067. DOI:10.1371/journal.ppat.1000067 (Primary mouse macrophages and dendritic cells)

- Cen, H., Mao, F., Aronchik, I., Fuentes, R. J., and Firestone, G. L. (2008). [DEVD-NucView488: a novel class of enzyme substrates for real-time detection of caspase-3 activity in live cells](#). *FASEB J* 22, 2243-2252. DOI:10.1096/fj.07-099234 (Cell lines: Jurkat (human T lymphocyte), HeLa (human cervical cancer))
- Cestari, I., Ansa-Addo, E., Deolindo, P., Inal, J. M., and Ramirez, M. I. (2012). [Trypanosoma cruzi immune evasion mediated by host cell-derived microvesicles](#). *J Immunol* 188, 1942-1952. 10.4049/jimmunol.1102053 (Cell line: THP-1 cells (human monocyte))
- Claudio, E., Saret, S., Wang, H., and Siebenlist, U. (2009). [Cell-autonomous role for NF-kappa B in immature bone marrow B cells](#). *J Immunol* 182, 3406-3413. DOI:10.4049/jimmunol.0803360 (Primary mouse immature B cells)
- Contreras, L., Calderon, R. I., Varela-Ramirez, A., Zhang, H. Y., Quan, Y., Das, U., Dimmock, J. R., Skouta, R., and Aguilera, R. J. (2018). [Induction of apoptosis via proteasome inhibition in leukemia/lymphoma cells by two potent piperidones](#). *Cell Oncol (Dordr)*. 10.1007/s13402-018-0397-1 (Cell lines: HL-60, Ramos)
- Cunha, B., Aguiar, T., Silva, M. M., Silva, R. J., Sousa, M. F., Pineda, E., Peixoto, C., Carrondo, M. J., Serra, M., and Alves, P. M. (2015). [Exploring continuous and integrated strategies for the up- and downstream processing of human mesenchymal stem cells](#). *J Biotechnol*. 10.1016/j.jbiotec.2015.02.023 (Cell lines: HFF, stem cells)
- Dahlawi, H., Jordan-Mahy, N., Clench, M., McDougall, G. J., and Le Maitre, C. L. (2013). [Polyphenols are responsible for the proapoptotic properties of pomegranate juice on leukemia cell lines](#). *Food Science & Nutrition* 1, 196-208. DOI:10.1002/fsn3.26 (Cell lines: CCRF-CEM, HL-60, and MOLT-3 (human leukemia) and THP-1 (human monocyte))
- de Sica, R. C., Rodrigues, C. J., Maria, D. A., and Cuce, L. C. (2016). [Study of 1550nm Erbium Glass Laser Fractional non-ablative treatment of photoaging: Comparative clinical effects, histopathology, electron microscopy and immunohistochemistry](#). *J Cosmet Laser Ther*, 1-31. 10.3109/14764172.2015.1114645 (Cell type: human skin biopsy cells)
- Dikovskaya, D., Cole, J. J., Mason, S. M., Nixon, C., Karim, S. A., McGarry, L., Clark, W., Hewitt, R. N., Sammons, M. A., Zhu, J., et al. (2015). [Mitotic Stress Is an Integral Part of the Oncogene-Induced Senescence Program that Promotes Multinucleation and Cell Cycle Arrest](#). *Cell Rep* 12, 1483-1496. 10.1016/j.celrep.2015.07.055 (Cell type: primary human fibroblast)
- Eicher, C., Dewerth, A., Kirchner, B., Warmann, S. W., Fuchs, J., and Armeanu-Ebinger, S. (2011). [Treatment effects of the multikinase inhibitor sorafenib on hepatoblastoma cell lines and xenografts in NMRI-Foxn1\(nu\) mice](#). *Liver Int*. 10.1111/j.1478-3231.2011.02729.x (human hepatoblastoma lines HepT1, HUH6)
- Feng, J. H., Nakagawa-Goto, K., Lee, K. H., and Shyr, L. F. (2016). [A Novel Plant Sesquiterpene Lactone Derivative, DETD-35, Suppresses BRAFV600E Mutant Melanoma Growth and Overcomes Acquired Vemurafenib Resistance in Mice](#). *Mol Cancer Ther* 15, 1163-1176. 10.1158/1535-7163.MCT-15-0973 (Cell line: A-375)
- Gallerne, C., Prola, A., and Lemaire, C. (2013). [Hsp90 inhibition by PU-H71 induces apoptosis through endoplasmic reticulum stress and mitochondrial pathway in cancer cells and overcomes the resistance conferred by Bcl-2](#). *Biochim Biophys Acta*. 10.1016/j.bbamcr.2013.02.014 (Cell line: HeLa)
- Gomez, M. C., Pope, C. E., Kutner, R. H., Ricks, D. M., Lyons, L. A., Ruhe, M., Dumas, C., Lyons, J., Lopez, M., Dresser, B. L., et al. (2008). [Nuclear transfer of sand cat cells into enucleated domestic cat oocytes is affected by cryopreservation of donor cells](#). *Cloning Stem Cells* 10, 469-483. 10.1089/clo.2008.0021 (Cell type: primary sand cat skin fibroblasts)
- Handa, P., Tupper, J. C., Jordan, K. C., and Harlan, J. M. (2011). [FLIP \(Flice-like inhibitory protein\) suppresses cytoplasmic double-stranded-RNA-induced apoptosis and NF-kappaB and IRF3-mediated signaling](#). *Cell Commun Signal* 9, 16. 10.1186/1478-811X-9-16 (mouse embryonic fibroblasts)

Jangamreddy, J. R., Ghavami, S., Grabarek, J., Kratz, G., Wiechec, E., Fredriksson, B. A., Rao Pariti, R. K., Cieslar-Pobuda, A., Panigrahi, S., and Los, M. J. (2013). [Salinomycin induces activation of autophagy, mitophagy and affects mitochondrial polarity: differences between primary and cancer cells](#). *Biochim Biophys Acta* 1833, 2057-2069. 10.1016/j.bbamcr.2013.04.011

Jeong, J. C., Kim, M. S., Kim, T. H., and Kim, Y. K. (2009). [Kaempferol induces cell death through ERK and Akt-dependent down-regulation of XIAP and survivin in human glioma cells](#). *Neurochem Res* 34, 991-1001. 10.1007/s11064-008-9868-5 (Cell type: A172 human glioma cell line)

Kato, M., Ospelt, C., Gay, R. E., Gay, S., and Klein, K. (2013). [Dual role of autophagy in stress-induced cell death in rheumatoid arthritis synovial fibroblasts](#). *Arthritis Rheum.* 10.1002/art.38190 (Cell type: human synovial fibroblasts)

Knapp, D., Kannan, N., Pellacani, D., and Eaves, C. J. (2017). [Mass Cytometric Analysis Reveals Viable Activated Caspase-3\(+\) Luminal Progenitors in the Normal Adult Human Mammary Gland](#). *Cell Rep* 21, 1116-1126. 10.1016/j.celrep.2017.09.096 (Cells: mammary luminal progenitor cells)

Kopec, A. K., Spada, A. P., Contreras, P. C., Mackman, N., and Luyendyk, J. P. (2018). [Caspase Inhibition Reduces Hepatic Tissue Factor-Driven Coagulation In Vitro and In Vivo](#). *Toxicol Sci* 162, 396-405. 10.1093/toxsci/kfx268 (Cells: hepatocytes)

Li, K., Li, Y., Wu, W., Gordon, W. R., Chang, D. W., Lu, M., Scoggin, S., Fu, T., Vien, L., Histen, G., et al. (2008). [Modulation of Notch signaling by antibodies specific for the extracellular negative regulatory region of NOTCH3](#). *J Biol Chem* 283, 8046-8054. 10.1074/jbc.M800170200 (Cell types: 293-T)

Li, Y., and Shively, J. E. (2013). [CEACAM1 Regulates Fas-mediated Apoptosis in Jurkat T-cells via its Interaction with beta-catenin](#). *Exp Cell Res.* 10.1016/j.yexcr.2013.02.020 (Cell line: Jurkat)

Macsik, L. L., Somogyi, I., Opper, B., Bovari-Biri, J., Pollak, E., Molnar, L., Nemeth, P., and Engelmann, P. (2015). [Induction of apoptosis-like cell death by coelomocyte extracts from Eisenia andrei earthworms](#). *Mol Immunol* 67, 213-222. 10.1016/j.molimm.2015.05.015 (Cell line: SP2/O-AG14)

Oldham, E. D., Nunes, L. M., Varela-Ramirez, A., Rankin, S. E., Knutson, B. L., Aguilera, R. J., and Lehmler, H. J. (2015). [Cytotoxic activity of triazole-containing alkyl beta-D-glucopyranosides on a human T-cell leukemia cell line](#). *Chem Cent J* 9, 3. 10.1186/s13065-014-0072-1 (Cell line: Jurkat)

Park, J. Y., Kim, M. J., Kim, Y. K., and Woo, J. S. (2011). [Ceramide induces apoptosis via caspase-dependent and caspase-independent pathways in mesenchymal stem cells derived from human adipose tissue](#). *Arch Toxicol* 85, 1057-1065. 10.1007/s00204-011-0645-x (Cell type: adipose mesenchymal stem cells)

Pedroza, D. A., De Leon, F., Varela-Ramirez, A., Lema, C., Aguilera, R. J., and Mito, S. (2014). [The cytotoxic effect of 2-acylated-1,4-naphthoquinones on leukemia/lymphoma cells](#). *Bioorg Med Chem* 22, 842-847. 10.1016/j.bmc.2013.12.007 (CEM T cell line)

Prola, A., Pires Da Silva, J., Guilbert, A., Lecru, L., Piquereau, J., Ribeiro, M., Mateo, P., Gressette, M., Fortin, D., Boursier, C., et al. (2017). [SIRT1 protects the heart from ER stress-induced cell death through eIF2alpha deacetylation](#). *Cell Death Differ* 24, 343-356. 10.1038/cdd.2016.138 (Cell line: H9c2)

Robles-Escajeda, E., Das, U., Ortega, N. M., Parra, K., Francia, G., Dimmock, J. R., Varela-Ramirez, A., and Aguilera, R. J. (2016). [A novel curcumin-like dienone induces apoptosis in triple-negative breast cancer cells](#). *Cell Oncol (Dordr)* 39, 265-277. 10.1007/s13402-016-0272-x (Cell line: MDA-MB-231)

Robles-Escajeda, E., Lerma, D., Nyakeriga, A. M., Ross, J. A., Kirken, R. A., Aguilera, R. J., and Varela-Ramirez, A. (2013). [Searching in mother nature for anti-cancer activity: anti-proliferative and pro-apoptotic effect elicited by green barley on leukemia/lymphoma cells](#). *PLoS One* 8, e73508. 10.1371/journal.pone.0073508 (NALM-6 cells)

Shulga, N., and Pastorino, J. G. (2015). [Fructose sensitizes pancreatic beta cells to TNFalpha-induced necroptosis](#). Biol Open. 10.1242/bio.014712 (Cell line: INS-IE (mouse pancreatic islet))

Stewart, S. G., Braun, C. J., Ng, S. L., Polomska, M. E., Karimi, M., and Abraham, L. J. (2010). [New thalidomide analogues derived through Sonogashira or Suzuki reactions and their TNF expression inhibition profiles](#). Bioorg Med Chem 18, 650-662. 10.1016/j.bmc.2009.12.001 (Cell line: Jurkat)

Su, N. W., Wu, S. H., Chi, C. W., Liu, C. J., Tsai, T. H., and Chen, Y. J. (2017). [Metronomic Cordycepin Therapy Prolongs Survival of Oral Cancer-Bearing Mice and Inhibits Epithelial-Mesenchymal Transition](#). Molecules 22. 10.3390/molecules22040629 (Cell lines: SAS, OECM-1)

Sztiller-Sikorska, M., Jakubowska, J., Wozniak, M., Stasiak, M., and Czyz, M. (2009). [A non-apoptotic function of caspase-3 in pharmacologically-induced differentiation of K562 cells](#). Br J Pharmacol 157, 1451-1462. 10.1111/j.1476-5381.2009.00333.x (Cell types: K562 cells)

Takata, N., Ohshima, Y., Suzuki-Karasaki, M., Yoshida, Y., Tokuhashi, Y., and Suzuki-Karasaki, Y. (2017). [Mitochondrial Ca²⁺ removal amplifies TRAIL cytotoxicity toward apoptosis-resistant tumor cells via promotion of multiple cell death modalities](#). Int J Oncol 51, 193-203. 10.3892/ijo.2017.4020 (Cell lines: HOS, A-375)

Tribulatti, M. V., Mucci, J., Cattaneo, V., Agüero, F., Gilmartin, T., Head, S. R., and Campetella, O. (2007). [Galectin-8 induces apoptosis in the CD4\(high\)CD8\(high\) thymocyte subpopulation](#). Glycobiology 17, 1404-1412. 10.1093/glycob/cwm104 (Cell type: mouse thymocytes)

Valente, M., Baey, C., Louche, P., Dutertre, C. A., Vimeux, L., Maranon, C., Hosmalin, A., and Feuillet, V. (2014). [Apoptotic cell capture by DCs induces unexpectedly robust autologous CD4 T-cell responses](#). Eur J Immunol. 10.1002/eji.201344191 (Cell type: primary human T cells)

Wang, H., Liu, J., Wang, F., Chen, M., Xiao, Z., Ouyang, R., Fei, A., Shen, Y., and Pan, S. (2013). [The role of charged multivesicular body protein 5 in programmed cell death in leukemic cells](#). Acta Biochim Biophys Sin (Shanghai) 45, 383-390. 10.1093/abbs/gmt028 (Cell line: U937)

Zhang, X., Paun, A., Claudio, E., Wang, H., and Siebenlist, U. (2013). [The tumor promoter and NF-kappaB modulator Bcl-3 regulates splenic B cell development](#). J Immunol 191, 5984-5992. 10.4049/jimmunol.1300611 (Cell type: primary mouse B cells)

NucView® 530 in microscopy

Bueno, O., Estevez Gallego, J., Martins, S., Prota, A. E., Gago, F., Gomez-SanJuan, A., Camarasa, M. J., Barasoain, I., Steinmetz, M. O., Diaz, J. F., et al. (2018). [High-affinity ligands of the colchicine domain in tubulin based on a structure-guided design](#). Sci Rep 8, 4242. 10.1038/s41598-018-22382-x (Cell line: MDA-MB-231)

Manas, A., Wang, S., Nelson, A., Li, J., Zhao, Y., Zhang, H., Davis, A., Xie, B., Maltsev, N., and Xiang, J. (2017). [The functional domains for Bax2 aggregate-mediated caspase 8-dependent cell death](#). Exp Cell Res 359, 342-355. 10.1016/j.yexcr.2017.08.016 (Cell line: HCT116)

Yao, Q., Zhang, H., Zhao, Y., Ye, Z., Lee, Y. J., and Xiang, J. (2018). [The C-terminus of Ubl4A is critical for pro-death activity and association with the Arp2/3 complex](#). Biochem Biophys Res Commun 503, 3192-3197. 10.1016/j.bbrc.2018.08.123 (Cell type: MEFs)

NucView® 488 in microscopy

Adhikari, S. (2008). [Role of mitogen-activated protein kinases in crambene-induced pancreatic acinar cell apoptosis](#). International Journal of Integrative Biology 3 (Cell type: Primary mouse pancreatic acinar cells)

- Aftab, O., Nazir, M., Fryknas, M., Hammerling, U., Larsson, R., and Gustafsson, M. G. (2014). [Label free high throughput screening for apoptosis inducing chemicals using time-lapse microscopy signal processing](#). *Apoptosis*. 10.1007/s10495-014-1009-9 (Cell line: HCT116 (human colorectal carcinoma))
- Alinezhad, S., Vaananen, R. M., Mattsson, J., Li, Y., Tallgren, T., Tong Ochoa, N., Bjartell, A., Akerfelt, M., Taimen, P., Bostrom, P. J., et al. (2016). [Validation of Novel Biomarkers for Prostate Cancer Progression by the Combination of Bioinformatics, Clinical and Functional Studies](#). *PLoS One* 11, e0155901. 10.1371/journal.pone.0155901 (Cell type: VCaP spheroids)
- Angeloni, M. B., Silva, N. M., Castro, A. S., Gomes, A. O., Silva, D. A., Mineo, J. R., and Ferro, E. A. (2009). [Apoptosis and S phase of the cell cycle in BeWo trophoblastic and HeLa cells are differentially modulated by Toxoplasma gondii strain types](#). *Placenta* 30, 785-791. DOI:10.1016/j.placenta.2009.07.002 (Cell lines: BeWo (human trophoblast), HeLa (human cervical cancer))
- Angelova, P. R., Baev, A. Y., Berezhnov, A. V., and Abramov, A. Y. (2016). [Role of inorganic polyphosphate in mammalian cells: from signal transduction and mitochondrial metabolism to cell death](#). *Biochem Soc Trans* 44, 40-45. 10.1042/BST20150223 (Cell type: primary rat astrocyte)
- Antczak, C., Takagi, T., Ramirez, C. N., Radu, C., and Djaballah, H. (2009). [Live-cell imaging of caspase activation for high-content screening](#). *J Biomol Screen* 14, 956-969. DOI:10.1177/1087057109343207 (Cell line: HeLa (human cervical cancer))
- Aronchik, I., Appleton, B. A., Basham, S. E., Crawford, K., Del Rosario, M., Doyle, L. V., Estacio, W. F., Lan, J., Lindvall, M. K., Luu, C. A., et al. (2014). [Novel Potent and Selective Inhibitors of p90 Ribosomal S6 Kinase Reveal the Heterogeneity of RSK Function in MAPK Driven Cancers](#). *Mol Cancer Res*. 10.1158/1541-7786.MCR-13-0595 (Cell lines: MDA-MB-231, H358)
- Arsic, N., Ho-Pun-Cheung, A., Evelyne, C., Assenat, E., Jarlier, M., Anguille, C., Colard, M., Pezet, M., Roux, P., and Gadea, G. (2017). [The p53 isoform delta133p53ss regulates cancer cell apoptosis in a RhoB-dependent manner](#). *PLoS One* 12, e0172125. 10.1371/journal.pone.0172125 (Cell lines: CCL-227, CCL-228 (colon cancer))
- Awasthi, B. P., Kathuria, M., Pant, G., Kumari, N., and Mitra, K. (2016). [Plumbagin, a plant-derived naphthoquinone metabolite induces mitochondria mediated apoptosis-like cell death in Leishmania donovani: an ultrastructural and physiological study](#). *Apoptosis* 21, 941-953. 10.1007/s10495-016-1259-9 (Cell type: Leishmania)
- Baar, M. P., Brandt, R. M. C., Putavet, D. A., Klein, J. D. D., Derks, K. W. J., Bourgeois, B. R. M., Stryeck, S., Rijksen, Y., van Willigenburg, H., Feijtel, D. A., et al. (2017). [Targeted Apoptosis of Senescent Cells Restores Tissue Homeostasis in Response to Chemotoxicity and Aging](#). *Cell* 169, 132-147 e116. 10.1016/j.cell.2017.02.031 (Cell line: IMR-90)
- Balez, R., Steiner, N., Engel, M., Munoz, S. S., Lum, J. S., Wu, Y., Wang, D., Vallotton, P., Sachdev, P., O'Connor, M., et al. (2016). [Neuroprotective effects of apigenin against inflammation, neuronal excitability and apoptosis in an induced pluripotent stem cell model of Alzheimer's disease](#). *Sci Rep* 6, 31450. 10.1038/srep31450 (Cell type: primary neurons)
- Balijepalli, M. K., Tandra, S., and Pichika, M. R. (2010). [Antiproliferative activity and induction of apoptosis in estrogen receptor-positive and negative human breast carcinoma cell lines by Gmelina asiatica roots](#). *Pharmacognosy Res* 2, 113-119. DOI:10.4103/0974-8490.62949 (Cell lines: MCF-7, MDA-MB-231 (human breast adenocarcinoma))
- Balzarini, J., Thomas, J., Liekens, S., Noppen, S., Dehaen, W., and Romagnoli, R. (2014). [2-aminothiophene-3-carboxylic acid ester derivatives as novel highly selective cytostatic agents](#). *Invest New Drugs* 32, 200-210. 10.1007/s10637-013-9981-4 (Cell line: HeLa, PC-3)
- Benetti, L., and Roizman, B. (2007). [In transduced cells, the US3 protein kinase of herpes simplex virus 1 precludes activation and induction of apoptosis by transfected procaspase 3](#). *J Virol* 81, 10242-10248. DOI:10.1128/JVI.00820-07 (Cell line: U2OS (human osteosarcoma))

- Bjork, J. K., Akerfelt, M., Joutsen, J., Puustinen, M. C., Cheng, F., Sistonen, L., and Nees, M. (2015). [Heat-shock factor 2 is a suppressor of prostate cancer invasion](#). *Oncogene*. 10.1038/onc.2015.241 (Cell line: PC-3; 3D culture)
- Bjorkman, M., Ostling, P., Harma, V., Virtanen, J., Mpindi, J. P., Rantala, J., Mirtti, T., Vesterinen, T., Lundin, M., Sankila, A., et al. (2011). [Systematic knockdown of epigenetic enzymes identifies a novel histone demethylase PHF8 overexpressed in prostate cancer with an impact on cell proliferation, migration and invasion](#). *Oncogene*. DOI:10.1038/onc.2011.512 (Cell line: PC-3 (human prostate cancer))
- Bosch, M., and Franklin-Tong, V. E. (2007). [Temporal and spatial activation of caspase-like enzymes induced by self-incompatibility in Papaver pollen](#). *Proc Natl Acad Sci U S A* 104, 18327-18332. DOI:10.1073/pnas.0705826104 (Papaver rhoeas (field poppy) pollen tube)
- Brassescio, M. S., Pezuk, J. A., de Oliveira, J. C., Valera, E. T., de Oliveira, H. F., Scrideli, C. A., Umezawa, K., and Tone, L. G. (2013). [Activator Protein-1 Inhibition by 3-\[\[Dodecylthiocarbonyl\]Methyl\]-Glutamaride Impairs Invasion and Radiosensitizes Osteosarcoma Cells In Vitro](#). *Cancer Biother Radiopharm*. DOI:10.1089/cbr.2012.1305 (Cell line: HOS (human osteosarcoma))
- Brassescio, M. S., Pezuk, J. A., Morales, A. G., de Oliveira, J. C., Roberto, G. M., da Silva, G. N., Francisco de Oliveira, H., Scrideli, C. A., and Tone, L. G. (2013). [In vitro targeting of Polo-like kinase 1 in bladder carcinoma: comparative effects of four potent inhibitors](#). *Cancer Biol Ther* 14, 648-657. 10.4161/cbt.25087 (Cell lines: RT4, 5637, and T24)
- Cai, Y., Geutjes, E. J., de Lint, K., Roepman, P., Bruurs, L., Yu, L. R., Wang, W., van Blijswijk, J., Mohammad, H., de Rink, I., et al. (2013). [The NuRD complex cooperates with DNMTs to maintain silencing of key colorectal tumor suppressor genes](#). *Oncogene*. 10.1038/onc.2013.178 (RKO cells; CHD4 KD + DAC treatment)
- Cen, H., Mao, F., Aronchik, I., Fuentes, R. J., and Firestone, G. L. (2008). [DEVD-NucView488: a novel class of enzyme substrates for real-time detection of caspase-3 activity in live cells](#). *FASEB J* 22, 2243-2252. DOI:10.1096/fj.07-099234 (Cell lines: Jurkat (human T lymphocyte), HeLa (human cervical cancer))
- Cheli, V. T., Santiago Gonzalez, D. A., Spreuer, V., and Paez, P. M. (2015). [Voltage-gated Ca²⁺ entry promotes oligodendrocyte progenitor cell maturation and myelination in vitro](#). *Exp Neurol* 265, 69-83. 10.1016/j.expneurol.2014.12.012 (Cell type: mouse primary oligodendrocyte progenitor cells)
- Chen, C. H., Pan, J., Di, Y. Q., Liu, W., Hou, L., Wang, J. X., and Zhao, X. F. (2017). [Protein kinase C delta phosphorylates ecdysone receptor B1 to promote gene expression and apoptosis under 20-hydroxyecdysone regulation](#). *Proc Natl Acad Sci U S A* 114, E7121-E7130. 10.1073/pnas.1704999114 (Cell line: HaEpi (insect cells))
- Cheng, G., Tse, J., Jain, R. K., and Munn, L. L. (2009). [Micro-environmental mechanical stress controls tumor spheroid size and morphology by suppressing proliferation and inducing apoptosis in cancer cells](#). *PLoS One* 4, e4632. DOI:10.1371/journal.pone.0004632 (Cell line: 67NR (mouse mammary carcinoma))
- Chiu, S. H., Gedda, G., Girma, W. M., Chen, J. K., Ling, Y. C., Ghule, A. V., Ou, K. L., and Chang, J. Y. (2016). [Rapid fabrication of carbon quantum dots as multifunctional nanovehicles for dual-modal targeted imaging and chemotherapy](#). *Acta Biomater* 46, 151-164. 10.1016/j.actbio.2016.09.027 (Cell lines: HeLa, HepG2)
- Choo, B. H.A., and Fong, W.J. (2013) [Use of markers of undifferentiated pluripotent stem cells](#). (2013) United States Patent Application US 8906678 B2 (Cell type: human stem cells)
- Cribbes, S., Kessel, S., McMenemy, S., Qiu, J., and Chan, L. L. (2017). [A Novel Multiparametric Drug-Scoring Method for High-Throughput Screening of 3D Multicellular Tumor Spheroids Using the Celigo Image Cytometer](#). *SLAS Discov* 22, 547-557. 10.1177/2472555217689884 (Cell line: U-87 MG spheroids)

- Daya, S., Roberts, M., Isherwood, B., Ingleston-Orme, A., Caie, P., Teobald, I., Eagle, R., and Carragher, N. (2010). [Integrating an automated in vitro combination screening platform with live-cell and endpoint phenotypic assays to support the testing of drug combinations](#). Paper presented at: SBS 16th Annual Conference & Exhibition (Phoenix, Arizona.)
- De, P., Carlson, J. H., Leyland-Jones, B., Williams, C., and Dey, N. (2018). [Triple Fluorescence staining to Evaluate Mechanism-based Apoptosis following Chemotherapeutic and Targeted Anti-cancer Drugs in Live Tumor Cells](#). *Sci Rep* 8, 13192. 10.1038/s41598-018-31575-3 (Cell line: OVK18)
- Dereli-Korkut, Z., Akaydin, H. D., Ahmed, A. H., Jiang, X., and Wang, S. (2014). [Three Dimensional Microfluidic Cell Arrays for ex Vivo Drug Screening with Mimicked Vascular Flow](#). *Anal Chem*. 10.1021/ac403899j (Cell line: PC9 (human non-small cell lung cancer))
- Dewerth, A., Wonner, T., Lieber, J., Ellerkamp, V., Warmann, S. W., Fuchs, J., and Armeanu-Ebinger, S. (2012). [In vitro evaluation of the Aurora kinase inhibitor VX-680 for Hepatoblastoma](#). *Pediatr Surg Int*. DOI:10.1007/s00383-012-3086-6 (Cell line: HepT1 and HUH6 (human hepatoblastoma))
- Dufer, M., Neye, Y., Horth, K., Krippeit-Drews, P., Hennige, A., Widmer, H., McClafferty, H., Shipston, M. J., Haring, H. U., Ruth, P., et al. (2011). [BK channels affect glucose homeostasis and cell viability of murine pancreatic beta cells](#). *Diabetologia* 54, 423-432. DOI:10.1007/s00125-010-1936-0 (Primary rat pancreatic beta cells)
- Eisenhoffer, G. T., Loftus, P. D., Yoshigi, M., Otsuna, H., Chien, C. B., Morcos, P. A., and Rosenblatt, J. (2012). [Crowding induces live cell extrusion to maintain homeostatic cell numbers in epithelia](#). *Nature*. 10.1038/nature10999 (MDCK)
- Eriksson, A., Hermanson, M., Wickstrom, M., Lindhagen, E., Ekholm, C., Jenmalm Jensen, A., Lothgren, A., Lehmann, F., Larsson, R., Parrow, V., et al. (2012). [The novel tyrosine kinase inhibitor AKN-028 has significant antileukemic activity in cell lines and primary cultures of acute myeloid leukemia](#). *Blood Cancer J* 2, e81. 10.1038/bcj.2012.28 (Cell line: MV4-11 (primary human leukemia))
- Gao, F., Zheng, W., Gao, L., Cai, P., Liu, R., Wang, Y., Yuan, Q., Zhao, Y., and Gao, X. (2017). [Au Nanoclusters and Photosensitizer Dual Loaded Spatiotemporal Controllable Liposomal Nanocomposites Enhance Tumor Photodynamic Therapy Effect by Inhibiting Thioredoxin Reductase](#). *Adv Healthc Mater* 6. 10.1002/adhm.201601453 (Cell line: MDA-MB-361)
- Gao, L., Liu, R., Gao, F., Wang, Y., Jiang, X., and Gao, X. (2014). [Plasmon-mediated generation of reactive oxygen species from near-infrared light excited gold nanocages for photodynamic therapy in vitro](#). *ACS Nano* 8, 7260-7271. 10.1021/nn502325j (Cell line: HeLa)
- Geng, L., Rachakonda, G., Morre, D. J., Morre, D. M., Crooks, P. A., Sonar, V. N., Roti, J. L., Rogers, B. E., Greco, S., Ye, F., et al. (2009). [Indolyl-quinuclidinols inhibit ENOX activity and endothelial cell morphogenesis while enhancing radiation-mediated control of tumor vasculature](#). *FASEB J* 23, 2986-2995. 10.1096/fj.09-130005 (Cell line: HUVEC (Human Umbilical Vascular Epithelial))
- Gier, B., Krippeit-Drews, P., Sheiko, T., Aguilar-Bryan, L., Bryan, J., Dufer, M., and Drews, G. (2009). [Suppression of KATP channel activity protects murine pancreatic beta cells against oxidative stress](#). *J Clin Invest* 119, 3246-3256. 10.1172/JCI38817 (Cell types: primary mouse pancreatic islet cells)
- Granot, Z., Henke, E., Comen, E. A., King, T. A., Norton, L., and Benezra, R. (2011). [Tumor entrained neutrophils inhibit seeding in the premetastatic lung](#). *Cancer Cell* 20, 300-314. 10.1016/j.ccr.2011.08.012 (Cell line: 4T1 (mammary tumor cells))
- Gualda, E. J., Simao, D., Pinto, C., Alves, P. M., and Brito, C. (2014). [Imaging of human differentiated 3D neural aggregates using light sheet fluorescence microscopy](#). *Front Cell Neurosci* 8, 221. 10.3389/fncel.2014.00221 (Cell type: neural aggregates)
- Guardia Clausi, M., Paez, P. M., Campagnoni, A. T., Pasquini, L. A., and Pasquini, J. M. (2012). [Intranasal administration of aTf protects and repairs the neonatal white matter after a cerebral hypoxic-ischemic event](#). *Glia*. 10.1002/glia.22374 (Cell type: primary mouse oligodendrocyte precursor cells)

- Hafner, M., Niepel, M., Chung, M., and Sorger, P. K. (2016). [Growth rate inhibition metrics correct for confounders in measuring sensitivity to cancer drugs](#). *Nat Methods* 13, 521-527. 10.1038/nmeth.3853 (Cell line: MCF10a)
- Hamzeloo-Moghadam, M., Khalaj, A., and Malekmohammadi, M. (2015). [Cytotoxic Activity and Apoptosis Induction of Hypericum scabrum L.](#) *Iran Red Crescent Med J* 17, e19453. 10.5812/ircmj.19453 (Cell line: MCF-7)
- Handa, P., Tupper, J. C., Jordan, K. C., and Harlan, J. M. (2011). [FLIP \(Flice-like inhibitory protein\) suppresses cytoplasmic double-stranded-RNA-induced apoptosis and NF-kappaB and IRF3-mediated signaling](#). *Cell Commun Signal* 9, 16. 10.1186/1478-811X-9-16 (mouse embryonic fibroblasts)
- Hanley, P. J., Kronlage, M., Kirschning, C., Del Rey, A., Di Virgilio, F., Leipziger, J., Chessell, I. P., Sargin, S., Filippov, M. A., Lindemann, O., et al. (2012). [Transient P2x7 Receptor Activation Triggers Macrophage Death Independent of Tlr2/4, Casp1 and Panx1](#). *J Biol Chem* 287, 10650-10663. 10.1074/jbc.M111.332676 (Primary mouse peritoneal macrophages)
- He, J., Ma, C., Liu, W., and Wang, J. (2014). [On-chip monitoring of skeletal myoblast transplantation for the treatment of hypoxia-induced myocardial injury](#). *Analyst* 139, 4482-4490. 10.1039/c4an00697f (Cell line: H9C2 (myocardial cells))
- Hofmeister-Brix, A., Lenzen, S., and Baltrusch, S. (2013). [The ubiquitin-proteasome system regulates the stability and activity of the glucose sensor glucokinase in pancreatic beta-cells](#). *Biochem J* 456, 173-184. 10.1042/BJ20130262 (Cell line: MIN6)
- Isherwood, B. (2011). [Live Cell in Vitro and in Vivo Imaging Applications: Accelerating Drug Discovery](#). *Pharmaceutics* 3
- Ishii, K., Adachi, T., Imamura, K., Takano, S., Usui, K., Suzuki, K., Hamamoto, H., Watanabe, T., and Sekimizu, K. (2012). [Serratia marcescens induces apoptotic cell death in host immune cells via a lipopolysaccharide- and flagella-dependent mechanism](#). *J Biol Chem*. 10.1074/jbc.M112.399667 (Cell type: silkworm hemocytes)
- Jangamreddy, J. R., Ghavami, S., Grabarek, J., Kratz, G., Wiechec, E., Fredriksson, B. A., Rao Pariti, R. K., Cieslar-Pobuda, A., Panigrahi, S., and Los, M. J. (2013). [Salinomycin induces activation of autophagy, mitophagy and affects mitochondrial polarity: differences between primary and cancer cells](#). *Biochim Biophys Acta* 1833, 2057-2069. 10.1016/j.bbamcr.2013.04.011
- Jechlinger, M., Podsypanina, K., and Varmus, H. (2009). [Regulation of transgenes in three-dimensional cultures of primary mouse mammary cells demonstrates oncogene dependence and identifies cells that survive deinduction](#). *Genes Dev* 23, 1677-1688. 10.1101/gad.1801809 (Cell types: primary mouse mammary cells)
- Johnson, A., Bhattacharya, N., Hanna, M., Pennington, J. G., Schuh, A. L., Wang, L., Otegui, M. S., Stagg, S. M., and Audhya, A. (2015). [TFG clusters COPII-coated transport carriers and promotes early secretory pathway organization](#). *EMBO J* 34, 811-827. 10.15252/embj.201489032 (Cell line: RPE-1)
- Johnston, H. J., Boys, S. K., Makda, A., Carragher, N. O., and Hulme, A. N. (2016). [Naturally Inspired Peptide Leads: Alanine Scanning Reveals an Actin-Targeting Thiazole Analogue of Bisebromoamide](#). *Chembiochem* 17, 1621-1627. 10.1002/cbic.201600257 (Cell line: HCT116)
- Jones, S. (2011). [Cell-Penetrating Peptides: Methods and Protocols](#). *Methods in Molecular Biology* 683, 291-303
- Kamynina, A. V., Holmstrom, K. M., Koroev, D. O., Volpina, O. M., and Abramov, A. Y. (2013). [Acetylcholine and antibodies against the acetylcholine receptor protect neurons and astrocytes against beta-amyloid toxicity](#). *Int J Biochem Cell Biol* 45, 899-907. 10.1016/j.biocel.2013.01.011 (Cell types: primary rat hippocampal & cortical neurons, glia)

- Kasim, N. R., Kuzelova, K., Holoubek, A., and Model, M. A. (2013). [Live fluorescence and transmission-through-dye microscopic study of actinomycin D-induced apoptosis and apoptotic volume decrease](#). *Apoptosis* 18, 521-532. 10.1007/s10495-013-0804-z (Cell line: HeLa)
- Kathuria, M., Bhattacharjee, A., Sashidhara, K. V., Singh, S. P., and Mitra, K. (2014). [Induction of Mitochondrial Dysfunction and Oxidative Stress in Leishmania donovani by Orally Active Clerodane Diterpene](#). *Antimicrob Agents Chemother* 58, 5916-5928. 10.1128/AAC.02459-14 (Cell type: leishmania)
- Katsoulis, E., Mabley, J. G., Samai, M., Sharpe, M. A., Green, I. C., and Chatterjee, P. K. (2010). [Lipotoxicity in renal proximal tubular cells: relationship between endoplasmic reticulum stress and oxidative stress pathways](#). *Free Radic Biol Med* 48, 1654-1662. 10.1016/j.freeradbiomed.2010.03.021 (Cell line: NRK-52E)
- Kawasaki, K., Ushioda, R., Ito, S., Ikeda, K., Masago, Y., and Nagata, K. (2015). [Deletion of the collagen-specific molecular chaperone Hsp47 causes endoplasmic reticulum stress-mediated apoptosis of hepatic stellate cells](#). *J Biol Chem* 290, 3639-3646. 10.1074/jbc.M114.592139 (Cell type: mouse hepatic stellate cells (HSC))
- King, M. A., Ganley, I. G., and Flemington, V. (2016). [Inhibition of cholesterol metabolism underlies synergy between mTOR pathway inhibition and chloroquine in bladder cancer cells](#). *Oncogene*. 10.1038/onc.2015.511 (Cell line: RT112)
- Klotz, D. M., Nelson, S. A., Kroboth, K., Newton, I. P., Radulescu, S., Ridgway, R. A., Sansom, O. J., Appleton, P. L., and Nathke, I. S. (2012). [The microtubule poison vinorelbine kills cells independently of mitotic arrest and targets cells lacking the APC tumour suppressor more effectively](#). *J Cell Sci*. 10.1242/jcs.091843 (U2OS (osteosarcoma))
- Kobos, R., Nagai, M., Tsuda, M., Merl, M. Y., Saito, T., Lae, M., Mo, Q., Olshen, A., Lianoglou, S., Leslie, C., et al. (2013). [Combining integrated genomics and functional genomics to dissect the biology of a cancer-associated, aberrant transcription factor, the ASPSCR1-TFE3 fusion oncoprotein](#). *J Pathol* 229, 743-754. 10.1002/path.4158 (Cell line: FU-UR-1 (renal cell carcinoma))
- Koerner, C., Keklikoglou, I., Bender, C., Woerner, A., Muenstermann, E., and Wiemann, S. (2013). [microRNA-31 sensitizes human breast cells to apoptosis by direct targeting of protein kinase C epsilon \(PKC{varepsilon}\)](#). *J Biol Chem*. 10.1074/jbc.M112.414128 (Cell lines: MCF10A, MDA-MB-231 (human breast cancer))
- Kovac, S., Domijan, A. M., Walker, M. C., and Abramov, A. Y. (2014). [Seizure activity results in calcium- and mitochondria-independent ROS production via NADPH and xanthine oxidase activation](#). *Cell Death Dis* 5, e1442. 10.1038/cddis.2014.390 (Cell type: cortical neurons)
- Kujjo, L. L., Ronningen, R., Ross, P., Pereira, R. J., Rodriguez, R., Beyhan, Z., Goissis, M. D., Baumann, T., Kagawa, W., Camsari, C., et al. (2012). [RAD51 plays a crucial role in halting cell death program induced by ionizing radiation in bovine oocytes](#). *Biol Reprod* 86, 76. 10.1095/biolreprod.111.092064 (Cell types: bovine and mouse oocytes)
- Lanz, H. L., Zimmerman, R. M., Brouwer, J., Noteborn, M. H., and Backendorf, C. (2013). [Mitotic catastrophe triggered in human cancer cells by the viral protein apoptin](#). *Cell Death Dis* 4, e487. 10.1038/cddis.2013.2 (Cell lines: U2OS, Saos-2 (human osteosarcoma))
- Lee, Y. C., Jin, J. K., Cheng, C. J., Huang, C. F., Song, J. H., Huang, M., Brown, W. S., Zhang, S., Yu-Lee, L. Y., Yeh, E. T., et al. (2013). [Targeting Constitutively Activated beta1 Integrin inhibits Prostate Cancer Metastasis](#). *Mol Cancer Res*. 10.1158/1541-7786.MCR-12-0551
- Lefort, R., Pozueta, J., and Shelanski, M. (2012). [Cross-Linking of Cell Surface Amyloid Precursor Protein Leads to Increased beta-Amyloid Peptide Production in Hippocampal Neurons: Implications for Alzheimer's Disease](#). *J Neurosci* 32, 10674-10685. 10.1523/JNEUROSCI.6473-11.2012 (Cell type: primary rat hippocampal neurons)

- Leuenroth, S. J., Okuhara, D., Shotwell, J. D., Markowitz, G. S., Yu, Z., Somlo, S., and Crews, C. M. (2007). [Triptolide is a traditional Chinese medicine-derived inhibitor of polycystic kidney disease](#). *Proc Natl Acad Sci U S A* 104, 4389-4394. 10.1073/pnas.0700499104 (Cell types: primary mouse kidney epithelial cells)
- Levy, D., Ruiz, J. L., Celestino, A. T., Silva, S. F., Ferreira, A. K., Isaac, C., and Bydlowski, S. P. (2014). [Short-term effects of 7-ketocholesterol on human adipose tissue mesenchymal stem cells in vitro](#). *Biochem Biophys Res Commun*. 10.1016/j.bbrc.2014.01.132 (Cell type: human adipose mesenchymal stem cells)
- Lim, S. O., Li, C. W., Xia, W., Cha, J. H., Chan, L. C., Wu, Y., Chang, S. S., Lin, W. C., Hsu, J. M., Hsu, Y. H., et al. (2016). [Deubiquitination and Stabilization of PD-L1 by CSN5](#). *Cancer Cell*. 10.1016/j.ccell.2016.10.010 (Cell line: BT549)
- Liu, W., Tian, C., Yan, M., Zhao, L., Ma, C., Li, T., Xu, J., and Wang, J. (2016). [Heterotypic 3D tumor culture in a reusable platform using pneumatic microfluidics](#). *Lab Chip* 16, 4106-4120. 10.1039/c6lc00996d (Cell lines: HepG2, U-251 MG, MKN-45, NIH 3T3)
- Liu, W., Xu, J., Li, T., Zhao, L., Ma, C., Shen, S., and Wang, J. (2015). [Monitoring tumor response to anticancer drugs using stable three-dimensional culture in a recyclable microfluidic platform](#). *Anal Chem* 87, 9752-9760. 10.1021/acs.analchem.5b01915 (Cell line: U251 (glioblastoma))
- Lu, B., Al-Ramahi, I., Valencia, A., Wang, Q., Berenshteyn, F., Yang, H., Gallego-Flores, T., Ichcho, S., Lacoste, A., Hild, M., et al. (2013). [Identification of NUB1 as a suppressor of mutant Huntingtin toxicity via enhanced protein clearance](#). *Nat Neurosci*. 10.1038/nn.3367 (Cell type: mouse striatal cells)
- Majewska, E., Wittek, N., Rysz, J., and Baj, Z. (2012). [The influence of uremic high cystatin C concentration on neutrophil apoptosis and selected neutrophil functions isolated from healthy subjects](#). *Med Sci Monit* 18, CR667-673 (Cell type: Primary human neutrophils)
- Mankus, C., Rich, C., Minns, M., and Trinkaus-Randall, V. (2011). [Corneal Epithelium Expresses a Variant of P2X\(7\) Receptor in Health and Disease](#). *PLoS One* 6, e28541. 10.1371/journal.pone.0028541
- Mau, M., Kalbe, C., Viergutz, T., Nurnberg, G., and Rehfeldt, C. (2008). [Effects of dietary isoflavones on proliferation and DNA integrity of myoblasts derived from newborn piglets](#). *Pediatr Res* 63, 39-45. 10.1203/PDR.0b013e31815b8e60 (Cell type: primary porcine myoblasts)
- Merlet, E., Atassi, F., Motiani, R. K., Mougnot, N., Jacquet, A., Nadaud, S., Capiod, T., Trebak, M., Lompre, A. M., and Marchand, A. (2013). [miR-424/322 regulates vascular smooth muscle cell phenotype and neointimal formation in the rat](#). *Cardiovasc Res*. 10.1093/cvr/cvt045 (Cell type: primary rat vascular smooth muscle)
- Moore, M. J., Wang, Q., Kennedy, C. J., and Silver, P. A. (2010). [An alternative splicing network links cell-cycle control to apoptosis](#). *Cell* 142, 625-636. 10.1016/j.cell.2010.07.019 (Cell line: HeLa)
- Naghibi, F., Khalaj, A., Mosaddegh, M., Malekmohamadi, M., and Hamzeloo-Moghadam, M. (2014). [Cytotoxic activity evaluation of some medicinal plants, selected from Iranian traditional medicine Pharmacopoeia to treat cancer and related disorders](#). *J Ethnopharmacol*. 10.1016/j.jep.2014.05.025 (Cell line: MCF7)
- Nakayama Wong, L. S., Lame, M. W., Jones, A. D., and Wilson, D. W. (2010). [Differential cellular responses to protein adducts of naphthoquinone and monocrotaline pyrrole](#). *Chem Res Toxicol* 23, 1504-1513. 10.1021/tx1002436 (Cell type: bronchial epithelial)
- Olivera-Martinez, I., Harada, H., Halley, P. A., and Storey, K. G. (2012). [Loss of FGF-Dependent Mesoderm Identity and Rise of Endogenous Retinoid Signalling Determine Cessation of Body Axis Elongation](#). *PLoS Biol* 10, e1001415. 10.1371/journal.pbio.1001415 (Chick embryos)

- Ousingsawat, J., Wanitchakool, P., Kmit, A., Romao, A. M., Jantarajit, W., Schreiber, R., and Kunzelmann, K. (2015). [Anoctamin 6 mediates effects essential for innate immunity downstream of P2X7 receptors in macrophages](#). *Nat Commun* 6, 6245. 10.1038/ncomms7245 (Cell type: primary mouse macrophages)
- Overmeyer, J. H., Kaul, A., Johnson, E. E., and Maltese, W. A. (2008). [Active ras triggers death in glioblastoma cells through hyperstimulation of macropinocytosis](#). *Mol Cancer Res* 6, 965-977. 10.1158/1541-7786.MCR-07-2036 (Cell types: U251)
- Paez, P. M., Fulton, D. J., Spreuer, V., Handley, V., Campagnoni, C. W., and Campagnoni, A. T. (2009). [Regulation of store-operated and voltage-operated Ca²⁺ channels in the proliferation and death of oligodendrocyte precursor cells by golli proteins](#). *ASN Neuro* 1. 10.1042/AN20090003 (Cell types: Primary mouse oligodendrocytes)
- Paez, P. M., Spreuer, V., Handley, V., Feng, J. M., Campagnoni, C., and Campagnoni, A. T. (2007). [Increased expression of golli myelin basic proteins enhances calcium influx into oligodendroglial cells](#). *J Neurosci* 27, 12690-12699. 10.1523/JNEUROSCI.2381-07.2007 (Cell types: N19 conditionally immortalized mouse oligodendrocyte cell line)
- Palorini, R., Votta, G., Pirola, Y., De Vitto, H., De Palma, S., Airoidi, C., Vasso, M., Ricciardiello, F., Lombardi, P. P., Cirulli, C., et al. (2016). [Protein Kinase A Activation Promotes Cancer Cell Resistance to Glucose Starvation and Anoikis](#). *PLoS Genet* 12, e1005931. 10.1371/journal.pgen.1005931 (Cell line: MDA-MB-231)
- Paromov, V., Brannon, M., Kumari, S., Samala, M., Qui, M., Smith, M., and Stone, W. L. (2011). [Sodium pyruvate modulates cell death pathways in HaCaT keratinocytes exposed to half-mustard gas](#). *Int J Toxicol* 30, 197-206. 10.1177/1091581810390824 (Cell line: HaCaT (keratinocytes))
- Pathak, S., Dorfmüller, H. C., Borodkin, V. S., and van Aalten, D. M. (2008). [Chemical dissection of the link between streptozotocin, O-GlcNAc, and pancreatic cell death](#). *Chem Biol* 15, 799-807. 10.1016/j.chembiol.2008.06.010 (Cell types: Min6)
- Puigvert, J. C., Huvneers, S., Fredriksson, L., op het Veld, M., van de Water, B., and Danen, E. H. (2009). [Cross-talk between integrins and oncogenes modulates chemosensitivity](#). *Mol Pharmacol* 75, 947-955. 10.1124/mol.108.051649 (Cell types: GE11)
- Ren, L., Liu, W., Wang, Y., Wang, J. C., Tu, Q., Juan, X., Liu, R., Shen, S., and Wang, J. (2012). [Investigation of Hypoxia-induced Myocardial Injury Dynamics in a Tissue Interface Mimicking Microfluidic Device](#). *Anal Chem*. 10.1021/ac3025812 (Cell line: H9C2)
- Renjini, A. P., Titus, S., Narayan, N., Murali, M., Jha, R. K., and Laloraya, M. (2014). [STAT3 and Mcl-1 unite to cause mesenchymal epithelial transition](#). *J Cell Sci*. 10.1242/jcs.138214
- Riches, K., Angelini, T. G., Mudhar, G. S., Kaye, J., Clark, E., Bailey, M. A., Sohrabi, S., Korossis, S., Walker, P. G., Scott, D. J., et al. (2013). [Exploring smooth muscle phenotype and function in a bioreactor model of abdominal aortic aneurysm](#). *J Transl Med* 11, 208. 10.1186/1479-5876-11-208 (human porcine smooth muscle cells; Incucyte)
- Salem, O., Erdem, N., Jung, J., Munstermann, E., Worner, A., Wilhelm, H., Wiemann, S., and Korner, C. (2016). [The highly expressed 5'isomiR of hsa-miR-140-3p contributes to the tumor-suppressive effects of miR-140 by reducing breast cancer proliferation and migration](#). *BMC Genomics* 17, 566. 10.1186/s12864-016-2869-x (Cell lines: MDA-MB-231, MDA-MB-468)
- Salpeter, S. J., Pozniak, Y., Merquiol, E., Ben-Nun, Y., Geiger, T., and Blum, G. (2015). [A novel cysteine cathepsin inhibitor yields macrophage cell death and mammary tumor regression](#). *Oncogene*. 10.1038/onc.2015.51 (Cell type: primary mouse macrophages)
- Santo, V. E., Estrada, M. F., Rebelo, S. P., Abreu, S., Silva, I., Pinto, C., Veloso, S. C., Serra, A. T., Boghaert, E., Alves, P. M., et al. (2016). [Adaptable stirred-tank culture strategies for large scale production of multicellular spheroid-based tumor cell models](#). *J Biotechnol* 221, 118-129. 10.1016/j.jbiotec.2016.01.031 (Cell lines: multiple human cancer cell lines; spheroids)

- Saw, T. B., Doostmohammadi, A., Nier, V., Kocgozlu, L., Thampi, S., Toyama, Y., Marcq, P., Lim, C. T., Yeomans, J. M., and Ladoux, B. (2017). [Topological defects in epithelia govern cell death and extrusion](#). *Nature* 544, 212-216. 10.1038/nature21718 (Cell line: MDCK)
- Schmid, U., Stopper, H., Schweda, F., Queisser, N., and Schupp, N. (2008). [Angiotensin II induces DNA damage in the kidney](#). *Cancer Res* 68, 9239-9246. 10.1158/0008-5472.CAN-08-1310 (Cell types: primary mouse kidney epithelial cells)
- Schmitt, H., Lenzen, S., and Baltrusch, S. (2011). [Glucokinase mediates coupling of glycolysis to mitochondrial metabolism but not to beta cell damage at high glucose exposure levels](#). *Diabetologia*. 10.1007/s00125-011-2133-5 (RINm5F and primary rat beta cells)
- Shaw, T. N., Stewart-Hutchinson, P. J., Strangward, P., Dandamudi, D. B., Coles, J. A., Villegas-Mendez, A., Gallego-Delgado, J., van Rooijen, N., Zindy, E., Rodriguez, A., et al. (2015). [Perivascular Arrest of CD8+ T Cells Is a Signature of Experimental Cerebral Malaria](#). *PLoS Pathog* 11, e1005210. 10.1371/journal.ppat.1005210 (Ex vivo mouse brains)
- Sirianant, L., Wanitchakool, P., Ousingsawat, J., Benedetto, R., Zormpa, A., Cabrita, I., Schreiber, R., and Kunzelmann, K. (2016). [Non-essential contribution of LRRC8A to volume regulation](#). *Pflugers Arch*. 10.1007/s00424-016-1789-6 (Cell line: HeLa)
- Smith, G. S., Voyer-Grant, J. A., and Harauz, G. (2012). [Monitoring cleaved caspase-3 activity and apoptosis of immortalized oligodendroglial cells using live-cell imaging and cleaveable fluorogenic-dye substrates following potassium-induced membrane depolarization](#). *J Vis Exp*. 10.3791/3422 (Cell line: N19 (oligodendrocyte))
- Son, M. Y., Choi, H., Han, Y. M., and Cho, Y. S. (2013). [Unveiling the critical role of REX1 in the regulation of human stem cell pluripotency](#). *Stem Cells* 31, 2374-2387. 10.1002/stem.1509 (Cell types: embryonic stem cells, HFFs)
- Soriano-Romani, L., Garcia-Posadas, L., Lopez-Garcia, A., Paraoan, L., and Diebold, Y. (2015). [Thrombospondin-1 induces differential response in human corneal and conjunctival epithelial cells lines under in vitro inflammatory and apoptotic conditions](#). *Exp Eye Res* 134, 1-14. 10.1016/j.exer.2015.03.004 (Cell lines: HCE(human corneal epithelial), IOBA-NHC(conjunctival epithelial))
- Standiford, L. R., Standiford, T. J., Newstead, M. J., Zeng, X., Ballinger, M. N., Kovach, M. A., Reka, A. K., and Bhan, U. (2012). [TLR4-dependent GM-CSF protects against lung injury in Gram-negative bacterial pneumonia](#). *Am J Physiol Lung Cell Mol Physiol* 302, L447-454. 10.1152/ajplung.00415.2010 (mouse alveolar epithelial cells)
- Stopper, H., Schupp, N., Fazeli, G., Dietel, B., Queisser, N., Walitza, S., and Gerlach, M. (2009). [Genotoxicity of the neurotransmitter dopamine in vitro](#). *Toxicol In Vitro* 23, 640-646. 10.1016/j.tiv.2009.03.001 (Cell types: TK6 (human lymphoblastoid), NRK (rat kidney), PC12 (rat pheochromocytoma))
- Tadokoro, T., Gao, X., Hong, C. C., Hotten, D., and Hogan, B. L. (2016). [BMP signaling and cellular dynamics during regeneration of airway epithelium from basal progenitors](#). *Development*. 10.1242/dev.126656 (Ex vivo mouse trachea)
- Tang, Y., Xie, T., Florian, S., Moerke, N., Shamu, C., Benes, C., and Mitchison, T. J. (2013). [Differential determinants of cancer cell insensitivity to antimetabolic drugs discriminated by a one-step cell imaging assay](#). *J Biomol Screen* 18, 1062-1071. 10.1177/1087057113493804 (5637 cells; paclitaxel treatment)
- Teng, H., Zhang, Z. G., Wang, L., Zhang, R. L., Zhang, L., Morris, D., Gregg, S. R., Wu, Z., Jiang, A., Lu, M., et al. (2008). [Coupling of angiogenesis and neurogenesis in cultured endothelial cells and neural progenitor cells after stroke](#). *J Cereb Blood Flow Metab* 28, 764-771. 10.1038/sj.jcbfm.9600573 (Cell types: Primary rat subventricular zone neural progenitors)
- Tyciakova, S., Matuskova, M., Bohovic, R., Polakova, K., Toro, L., Skolekova, S., and Kucerova, L. (2015). [Genetically engineered mesenchymal stromal cells producing TNFalpha have tumour suppressing effect on human melanoma xenograft](#). *J Gene Med* 17, 54-67. 10.1002/jgm.2823 (Cell lines: A375, SKBR3, MDA-MB-231, HT29, SK-OV-3, U-87 MG)

- Urbinati, G., Ali, H. M., Rousseau, Q., Chapuis, H., Desmaele, D., Couvreur, P., and Massaad-Massade, L. (2015). [Antineoplastic Effects of siRNA against TMPRSS2-ERG Junction Oncogene in Prostate Cancer](#). *PLoS One* 10, e0125277. 10.1371/journal.pone.0125277 (Cell line: VCaP)
- Virtanen, S. S., Kukkonen-Macchi, A., Vainio, M., Elima, K., Harkonen, P. L., Jalkanen, S., and Yegutkin, G. G. (2014). [Adenosine Inhibits Tumor Cell Invasion via Receptor-independent Mechanisms](#). *Mol Cancer Res*. 10.1158/1541-7786.MCR-14-0302-T (Cell line: PC3)
- Volosin, M., Trotter, C., Cragolini, A., Kenchappa, R. S., Light, M., Hempstead, B. L., Carter, B. D., and Friedman, W. J. (2008). [Induction of proneurotrophins and activation of p75NTR-mediated apoptosis via neurotrophin receptor-interacting factor in hippocampal neurons after seizures](#). *J Neurosci* 28, 9870-9879. 10.1523/JNEUROSCI.2841-08.2008 (Cell type: rat hippocampal neurons)
- Vuorinen, K., Gao, F., Oury, T. D., Kinnula, V. L., and Myllarniemi, M. (2007). [Imatinib mesylate inhibits fibrogenesis in asbestos-induced interstitial pneumonia](#). *Exp Lung Res* 33, 357-373. 10.1080/01902140701634827 (Cell types: CCL-190, primary human IPF fibroblasts)
- Wang, S., Ren, W., Liu, J., Lahat, G., Torres, K., Lopez, G., Lazar, A. J., Hayes-Jordan, A., Liu, K., Bankson, J., et al. (2010). [TRAIL and doxorubicin combination induces proapoptotic and antiangiogenic effects in soft tissue sarcoma in vivo](#). *Clin Cancer Res* 16, 2591-2604. 10.1158/1078-0432.CCR-09-2443 (Human soft tissue sarcoma cell lines)
- Wetzel-Smith, M. K., Hunkapiller, J., Bhangale, T. R., Srinivasan, K., Maloney, J. A., Atwal, J. K., Sa, S. M., Yaylaoglu, M. B., Foreman, O., Ortman, W., et al. (2014). [A rare mutation in UNC5C predisposes to late-onset Alzheimer's disease and increases neuronal cell death](#). *Nat Med* 20, 1452-1457. 10.1038/nm.3736 (Cell line: HEK293, time-lapse)
- Wu, D., Qiao, K., Feng, M., Fu, Y., Cai, J., Deng, Y., Tachibana, H., and Cheng, X. (2018). [Apoptosis of Acanthamoeba castellanii Trophozoites Induced by Oleic Acid](#). *J Eukaryot Microbiol* 65, 191-199. 10.1111/jeu.12454 (Cell type: Acanthamoeba)
- Wu, Y., and MacRae, T. H. (2010). [Truncation attenuates molecular chaperoning and apoptosis inhibition by p26, a small heat shock protein from Artemia franciscana](#). *Biochem Cell Biol* 88, 937-946. 10.1139/O10-143 (Cell line: 293H)
- Yamada, K., Asai, K., Nagayasu, F., Sato, K., Ijiri, N., Yoshii, N., Imahashi, Y., Watanabe, T., Tochino, Y., Kanazawa, H., et al. (2016). [Impaired nuclear factor erythroid 2-related factor 2 expression increases apoptosis of airway epithelial cells in patients with chronic obstructive pulmonary disease due to cigarette smoking](#). *BMC Pulm Med* 16, 27. 10.1186/s12890-016-0189-1 (Cell line: A549, in Incucyte)
- Yancey, P. G., Blakemore, J., Ding, L., Fan, D., Overton, C. D., Zhang, Y., Linton, M. F., and Fazio, S. (2010). [Macrophage LRP-1 controls plaque cellularity by regulating efferocytosis and Akt activation](#). *Arterioscler Thromb Vasc Biol* 30, 787-795. 10.1161/ATVBAHA.109.202051 (Primary mouse peritoneal macrophages)
- Yang, D., Elner, S. G., Chen, X., Field, M. G., Petty, H. R., and Elner, V. M. (2011). [MCP-1-activated monocytes induce apoptosis in human retinal pigment epithelium](#). *Invest Ophthalmol Vis Sci* 52, 6026-6034. 10.1167/iovs.10-7023 (Cell type: human retinal pigmented epithelial)
- Yang, D., Elner, S. G., Lin, L. R., Reddy, V. N., Petty, H. R., and Elner, V. M. (2009). [Association of superoxide anions with retinal pigment epithelial cell apoptosis induced by mononuclear phagocytes](#). *Invest Ophthalmol Vis Sci* 50, 4998-5005. 10.1167/iovs.09-3620 (Cell types: mouse RPE)
- Yao, Y., Cui, X., Al-Ramahi, I., Sun, X., Li, B., Hou, J., Difiglia, M., Palacino, J., Wu, Z. Y., Ma, L., et al. (2015). [A striatal-enriched intronic GPCR modulates huntingtin levels and toxicity](#). *Elife* 4. 10.7554/eLife.05449 (Cell type: human iPS-derived neurons)
- Yoshioka, Y., Fujibayashi, H., Kameda, K., Kan, D., Tone, S., and Tamura, M. (2015). [Induction of apoptosis in Caco-2 cells by exogenously added O\(2\)\(-\) produced by a nanodevice](#). *Exp Cell Res* 331, 408-415. 10.1016/j.yexcr.2014.12.007 (Cell line: Caco-2)

Zhang, C., Lee, S., Peng, Y., Bunker, E., Giaime, E., Shen, J., Zhou, Z., and Liu, X. (2014). [PINK1 triggers autocatalytic activation of Parkin to specify cell fate decisions](#). *Curr Biol* 24, 1854-1865. 10.1016/j.cub.2014.07.014 (Cell line: MEFs)

Zhao, W. L., Wang, D., Liu, C. Y., and Zhao, X. F. (2016). [G-protein-coupled receptor kinase 2 terminates G-protein-coupled receptor function in steroid hormone 20-hydroxyecdysone signaling](#). *Sci Rep* 6, 29205. 10.1038/srep29205 (Cell line: HaEpi (insect cells))

Zhou, Y., Sun, K., Ma, Y., Yang, H., Zhang, Y., Kong, X., and Wei, L. (2014). [Autophagy inhibits chemotherapy-induced apoptosis through downregulating Bad and Bim in hepatocellular carcinoma cells](#). *Sci Rep* 4, 5382. 10.1038/srep05382 (Cell line: SMMC-7721 (hepatocellular carcinoma))

Zhu, Q. S., Ren, W., Korchin, B., Lahat, G., Dicker, A., Lu, Y., Mills, G., Pollock, R. E., and Lev, D. (2008). [Soft tissue sarcoma cells are highly sensitive to AKT blockade: a role for p53-independent up-regulation of GADD45 alpha](#). *Cancer Res* 68, 2895-2903. 10.1158/0008-5472.CAN-07-6268 (Cell types: human soft tissue sarcoma cell lines)

NucView™ 488 in other applications

Chen, Y., Lewis, W., Diwan, A., Cheng, E. H., Matkovich, S. J., and Dorn, G. W., 2nd (2010). [Dual autonomous mitochondrial cell death pathways are activated by Nix/BNip3L and induce cardiomyopathy](#). *Proc Natl Acad Sci U S A* 107, 9035-9042. DOI:10.1073/pnas.0914013107 (Primary mouse embryonic fibroblast (MEF) cells)

Choo, B. H.A., and Fong, W.J. (2014) [Use of markers of undifferentiated pluripotent stem cells](#). (2013) United States Patent Application US 8906678 B2

Katsoulieiris, E., Mabley, J. G., Samai, M., Sharpe, M. A., Green, I. C., and Chatterjee, P. K. (2010). [Lipotoxicity in renal proximal tubular cells: relationship between endoplasmic reticulum stress and oxidative stress pathways](#). *Free Radic Biol Med* 48, 1654-1662. 10.1016/j.freeradbiomed.2010.03.021 (NRK-52E)

Monaco, G., Decrock, E., Akl, H., Ponsaerts, R., Vervliet, T., Luyten, T., De Maeyer, M., Missiaen, L., Distelhorst, C. W., De Smedt, H., et al. (2012). [Selective regulation of IP3-receptor-mediated Ca²⁺ signaling and apoptosis by the BH4 domain of Bcl-2 versus Bcl-Xl](#). *Cell Death Differ* 19, 295-309. 10.1038/cdd.2011.97 (Cell line: WEHI 7.2 (murine lymphoid))

Slee, E. A., Adrain, C., and Martin, S. J. (1999). [Serial killers: ordering caspase activation events in apoptosis](#). *Cell Death Differ* 6, 1067-1074. 10.1038/sj.cdd.4400601

Tang, Y., Xie, T., Florian, S., Moerke, N., Shamu, C., Benes, C., and Mitchison, T. J. (2013). [Differential determinants of cancer cell insensitivity to antimetabolic drugs discriminated by a one-step cell imaging assay](#). *J Biomol Screen* 18, 1062-1071. 10.1177/1087057113493804 (5637 cells; paclitaxel treatment)

Tanne, Y., Tanimoto, K., Okuma, S., Kunimatsu, R., Hirose, N., Mitsuyoshi, T., Tanne, K. (2013). [Effects of hyaluronan oligosaccharides on apoptosis of human gingival fibroblasts](#). *Open Journal of Stomatology* 3, 19-24 (Cell type: human gingival fibroblasts)

Uddin, M. I., Evans, S. M., Craft, J. R., Marnett, L. J., Uddin, M. J., and Jayagopal, A. (2015). [Applications of azo-based probes for imaging retinal hypoxia](#). *ACS Med Chem Lett* 6, 445-449. 10.1021/ml5005206 (Cell line: R28)

Verma, M., Shulga, N., and Pastorino, J. G. (2013). [Sirtuin-3 modulates Bak- and Bax-dependent apoptosis](#). *J Cell Sci* 126, 274-288. 10.1242/jcs.115188 (Cell line: HeLa)

Zhang, Y., Dong, L., Yang, X., Shi, H., and Zhang, L. (2011). [alpha-Linolenic acid prevents endoplasmic reticulum stress-mediated apoptosis of stearic acid lipotoxicity on primary rat hepatocytes](#). *Lipids Health Dis* 10, 81. 10.1186/1476-511X-10-81 (primary rat hepatocytes)