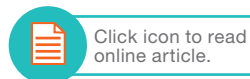


Publication List: ZOE™ Fluorescent Cell Imager



Cell Imaging

Bulletin 6865



Click icon to read
online article.

2016

Bodas M et al. (2016).

Nicotine exposure induces bronchial epithelial cell apoptosis and senescence via ROS mediated autophagy-impairment.
Free Radic Biol Med 97, 441–453.

Category: Fluorescent signal colocalization



Hayrabyan S et al. (2016).

Sertoli cells have a functional NALP3 inflammasome that can modulate autophagy and cytokine production.
Sci Rep 6, 18896.

Category: Fluorescent signal colocalisation



Kudryashova TV et al. (2016).

HIPPO-integrin linked kinase crosstalk controls self-sustaining proliferation and survival in pulmonary hypertension.
Am J of Respir Crit Care Med 194, 866–877.

Category: H&E staining of tissues, immunohistochemistry



Ramesan S et al. (2016).

Acoustically-driven thread-based tuneable gradient generators.
Lab Chip 16, 2,820–2,828.

Category: Fluorescence imaging of organelle staining, brightfield imaging



Sekihara S et al. (2016).

RNA interference directed against the Transglutaminase gene triggers dysbiosis of gut microbiota in *Drosophila*.

J Biol Chem [published online ahead of print October 19, 2016]. Accessed November 1, 2016.

Category: Immunohistochemistry



Vij N et al. (2016).

Cigarette smoke induced autophagy-impairment accelerates lung aging, COPD-emphysema exacerbations and pathogenesis.

Am J Physiol Cell Physiol Jul 13:ajpcell.00110.2016.

Category: Immunohistochemistry; fluorescence imaging of organelle staining



BIO-RAD



Walters HE et al. (2016).

Reversal of phenotypes of cellular senescence by pan-mTOR inhibition.

Aging 8, 231–243.

Category: Fluorescence imaging of organelle staining; brightfield imaging



Walworth K et al. (2016).

Dendrimer-based selective proteostasis-inhibition strategy to control NSCLC growth and progression.

PLoS ONE 11, e0158507.

Category: Immunocytochemistry



Zhang Q et al. (2016).

The effect of increasing concentrations of dl-methionine and 2-hydroxy-4 (methylthio) butanoic acid on hepatic genes controlling methionine regeneration and gluconeogenesis.

J Dairy Sci 99, 8,451–8,460.

Category: Brightfield imaging



Zhou Y et al. (2016).

microRNAs with AAGUGC seed motif constitute an integral part of an oncogenic signaling network.

Oncogene [published online ahead of print August 1, 2016]. Accessed November 1, 2016.

Category: Cell line validation



2015

Bennet JA et al. (2015).

Deficiency in aryl hydrocarbon receptor (AHR) expression throughout aging alters gene expression profiles in murine long-term hematopoietic stem cells.

PLoS One 10, e0133791.

Category: Estimation of transduction efficiency



Elçin S et al. (2015).

Synthesis and evaluation of fluorescence properties of Cu²⁺ selective azocalix[4] arenes and their application in living cell imaging.

Sensors and Actuators B: Chemical 234, 345–352.

Categories: Fluorescence probe imaging



Shivalingappa PC et al. (2015).

Airway exposure to e-cigarette vapors impairs autophagy and induces aggresome formation.

Antioxid Redox Signal 24, 186–204.

Category: Immunocytochemistry; brightfield imaging



Visit www.bio-rad.com/ZOEpublications for more information.



**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666
New Zealand 64 9 415 2280 **Norway** 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

