

**Viability/Cytotoxicity Assay Kit for Animal Live & Dead Cells References (with Cell Type)**  
**[Click on the blue link to see journal article](#)**

1. In Vitro Cell Dev Biol Animal (2007) 43, 361–370. <https://doi.org/10.1007/s11626-007-9051-3> (661W photoreceptor cells)
2. Molecular Vision (2007) 13, 1397-405. <http://www.molvis.org/molvis/v13/a154/> (whole rat lens)
3. Biomaterials (2010) 31, 4278–4285. [doi:10.1016/j.biomaterials.2010.01.059](https://doi.org/10.1016/j.biomaterials.2010.01.059) (human mesenchymal stem cells, scaffold culture)
4. J Biomed Biotechnol (2010) 2010, 858094. [doi: 10.1155/2010/858094](https://doi.org/10.1155/2010/858094) (microencapsulated cardiomyoblasts)
5. Cell Transplantation (2011) 20, 1881–1899. DOI: <http://dx.doi.org/10.3727/096368911X566181> (bone marrow mesenchymal stem cells, gelatin scaffold culture)
6. J Biomed Mat Res (2011) 98A(2), 245-256. <https://doi.org/10.1002/jbm.a.33089> (C2C12 cells, cellular microarray)
7. Biomed Aging Pathol (2012) 2(4), 143-150. <https://doi.org/10.1016/j.biomag.2012.10.003> (661W retinal cells)
8. Neurosci Lett (2012) 510(1) 14-19. <https://doi.org/10.1016/j.neulet.2011.12.061> (mouse cerebellar granule neurons)
9. Br J Nut (2013) 110, 20–28. [doi:10.1017/S0007114512004679](https://doi.org/10.1017/S0007114512004679) (primary salmon liver cells)
10. Cell Prolif (2013) 46, 685–694. [doi: 10.1111/cpr.12072](https://doi.org/10.1111/cpr.12072) (RAW 264.7 macrophages, TiO2 nanotube cultures)
11. Histol Histopathol (2013) 28, 1577-1583. DOI: [10.14670/HH-28.1577](https://doi.org/10.14670/HH-28.1577) (ARPE19 retinal cells cocultured with adipose-derived mesenchymal stem cells; Transwell inserts)
12. J Int Med Res (2013) 4(3), 688-696. <https://doi.org/10.1177/0300060513476591> (rabbit meniscal fragments)
13. J Mat Sci Mat Med (2013) 24, 189–198. <https://doi.org/10.1007/s10856-012-4785-1> (mesenchymal stem cell microarray)
14. J Mat Sci Mat Med (2013) 24, 1079–1091. <https://doi.org/10.1007/s10856-013-4869-6> (MC3T3-E1 cells, TiO2 nanotube cultures)
15. J Orthop Res (2013) 31, 1514–1519. DOI [10.1002/jor.22405](https://doi.org/10.1002/jor.22405) (porcine meniscal organ culture)
16. J Pharm Pharmacol (2013) 65(7), 1005-1013. <https://doi.org/10.1111/jphp.12054> (rat bone marrow stromal cells)
17. Transducers & Eurosensors XXVII (2013) 641-644. DOI: [10.1109/Transducers.2013.6626848](https://doi.org/10.1109/Transducers.2013.6626848) (HUVEC)
18. BBA Mol Basis Disease (2014) 1842(9), 1794-1805. <https://doi.org/10.1016/j.bbadis.2014.06.028> (rat lens, human lens epithelial cells (LECs))
19. BMC Plant Bio (2014) 14, 140. [doi:10.1186/1471-2229-14-140](https://doi.org/10.1186/1471-2229-14-140) (Caterpillar ventral eversible glands)
20. Cell Tissue Res (2014) 358, 705–716. (porcine RPE Transwell inserts, human AT-MSC co-culture)
21. Free Radic Biol Med (2014) 72, 134–148. [doi:10.1016/j.freeradbiomed.2014.04.010](https://doi.org/10.1016/j.freeradbiomed.2014.04.010) (mouse, human lens epithelial cells (LECS))
22. J Invest Dermatol (2014) 134(4), 1083-1090. <https://doi.org/10.1038/jid.2013.479> (primary human keratinocytes, melanocytes, dermal fibroblasts)
23. Oral Dis (2014) 20(5), 430-439. <https://doi.org/10.1111/odi.12148> (rabbit adipose-derived stem cells)

24. PLoS ONE (2014) 9(8), e104771. [doi:10.1371/journal.pone.0104771](https://doi.org/10.1371/journal.pone.0104771) (H9c2 cardiomyocytes)
25. J Am Heart Assoc (2015) 4(6), e002043. [doi: 10.1161/JAHA.115.002043](https://doi.org/10.1161/JAHA.115.002043) (primary human cardiac pericytes CorMatrix scaffold culture)
26. J Biomed Optics (2015) 20(9), 095005. [DOI: 10.1117/1.JBO.20.9.095005](https://doi.org/10.1117/1.JBO.20.9.095005) (human, rabbit, porcine ex vivo whole eye)
27. J Mat Chem B (2015) 3(45), 8749–8912. <https://doi.org/10.1039/C5TB01440A> (mouse bone marrow stromal cells, Fe-calcium phosphate discs)
28. Tissue Engineering Part C Meth (2015) 21(8), 852-861. <http://doi.org/10.1089/ten.tec.2014.0376> (human iPSC-derived cardiomyocyte spheroids)
29. Tissue Eng Part C Meth (2015) 21(11),1148-61. [doi: 10.1089/ten.TEC.2014.0709](https://doi.org/10.1089/ten.TEC.2014.0709) (MS-1 endothelial cells)
30. Toxicol in Vitro (2015) 30(1B), 492-505. <https://doi.org/10.1016/j.tiv.2015.09.012> (salmon hepatocytes)
31. ASAIO J (2016) 62(3), 340-348. [doi: 10.1097/MAT.0000000000000373](https://doi.org/10.1097/MAT.0000000000000373) (MS-1 endothelial cells)
32. J Neural Eng (2016) 13(1), 016006. [doi:10.1088/1741-2560/13/1/016006](https://doi.org/10.1088/1741-2560/13/1/016006) (rabbit lacrimal glands)
33. PeerJ (2016) 4, e2148. [DOI 10.7717/peerj.2148](https://doi.org/10.7717/peerj.2148) (human gingival fibroblasts, co-culture with *C. albicans* (calcofluor white co-stain))
34. PLoS One (2016) 11(6), e0157188. [doi:10.1371/journal.pone.0157188](https://doi.org/10.1371/journal.pone.0157188) (human gingival fibroblasts, co-culture with *C. albicans* (calcofluor white co-stain))
35. Proc SPIE (2016) 9693, 969319. [doi: 10.1117/12.2213171](https://doi.org/10.1117/12.2213171) (rabbit retina)
36. Sci Rep (2016) 6, 22020. [DOI: 10.1038/srep22020](https://doi.org/10.1038/srep22020) (human progenitor skin cells, NP/JATH fibroblasts)
37. Biophys J (2017) 113, 472–480. <http://dx.doi.org/10.1016/j.bpj.2017.06.014> (U-87, U251, DBTRG, NHA, hydrogels)
38. Xenotransplantation (2017) 24(1). [doi:10.1111/xen.12275](https://doi.org/10.1111/xen.12275) (porcine pancreatic islets)
39. Adv Healthcare Mater (2018) 7(9), 1701249. [DOI: 10.1002/adhm.201701249](https://doi.org/10.1002/adhm.201701249) (MCF7 spheroids, hydrogels)
40. Front Cell Infect Microbiol (2018) 8, 225. <https://doi.org/10.3389/fcimb.2018.00225> (porcine monocytes, flow cytometry)
41. IJP Drugs Drug Resist (2018) 8, 394–402. <https://doi.org/10.1016/j.ijpddr.2018.08.001> (PMJ2R cells, flow cytometry)
42. J Mat Sci Mat Med (2018) 29, 141. <https://doi.org/10.1007/s10856-018-6138-1> (MC3T3-E1 cells, TiO2 nanotubes)
43. Mol Neurobiol (2018) 55, 3611–3626. [DOI 10.1007/s12035-017-0478-6](https://doi.org/10.1007/s12035-017-0478-6) (HBMVP, bEnd3 cells, EVOS FL imager)
44. Polymer Chem (2018) 9(27), 3798-3807. [doi: 10.1039/C8PY00676H](https://doi.org/10.1039/C8PY00676H) (C3H10T1/2 cells)
45. BBRC (2019) 508, 654-659. <https://doi.org/10.1016/j.bbrc.2018.11.160> (hMSCs, HUVECs, HEK spheroids)
46. Biophys J (2019) 117, 975–986. <https://doi.org/10.1016/j.bpj.2019.07.041> (NIH-3T3 microtissues in collagen)
47. Colloids Surfaces B Biointerfaces (2019) 177, 121–129. <https://doi.org/10.1016/j.colsurfb.2019.01.054> (human corneal epithelial cells)
48. Life Sci (2019) 221, 249-258. <https://doi.org/10.1016/j.lfs.2019.02.035> (H9c2 cardiomyocytes)

49. Mat Sci Eng C (2019) 99, 1199-1212. <https://doi.org/10.1016/j.msec.2019.02.042> (mBMSC and HUVEC cells)
50. ACS Appl Mat Interfaces (2020) 12(4), 4308-4322. (MDA-MB-231 spheroids in Matrigel®)
51. Ceramics International (2020) 46, 353–364. <https://doi.org/10.1016/j.ceramint.2019.08.269> (hBMSCs)
52. Int J Pharmaceut (2020) 577, 118995. <https://doi.org/10.1016/j.ijpharm.2019.118995> (A549 spheroids)
53. Front Cell Infect Microbiol (2020) 9, 426. [doi: 10.3389/fcimb.2019.00426](https://doi.org/10.3389/fcimb.2019.00426) (*P. salmonis* LF-89 cells)