

JC-1 Mitochondrial Membrane Potential Detection Kit References

[Click on the blue link to see journal article](#)

1. Med Mol Morphol (2008) 41(4), 211-20. [doi: 10.1007/s00795-008-0414-y](https://doi.org/10.1007/s00795-008-0414-y)
2. Cancer Res (2009) 69, 9. [DOI: 10.1158/0008-5472.CAN-08-2747](https://doi.org/10.1158/0008-5472.CAN-08-2747)
3. Mitochondrion 10 (2010) 137–142. [doi:10.1016/j.mito.2009.12.144](https://doi.org/10.1016/j.mito.2009.12.144)
4. Breast Cancer Res Treat (2011) 128, 337–346. [DOI 10.1007/s10549-010-1085-7](https://doi.org/10.1007/s10549-010-1085-7)
5. Polish J Microbiol (2011) 60(3), 203–207. [PMID: 22184926](https://pubmed.ncbi.nlm.nih.gov/22184926/)
6. Synapse (2011) 65(7), 562-571. <https://doi.org/10.1002/syn.20875>
7. Ann Neurol (2012) 71(5), 687–698. [doi:10.1002/ana.23536](https://doi.org/10.1002/ana.23536)
8. BBRC (2012) 423(2), 436–440. [doi:10.1016/j.bbrc.2012.05.154](https://doi.org/10.1016/j.bbrc.2012.05.154)
9. BMC Neurosci (2012) 13, 11. <https://doi.org/10.1186/1471-2202-13-11>
10. Evid Based Complement Alt Med (2012) 517965. [doi: 10.1155/2012/517965](https://doi.org/10.1155/2012/517965)
11. J Biosci (2012) 37(6) 1005–1016. [DOI 10.1007/s12038-012-9265-8](https://doi.org/10.1007/s12038-012-9265-8)
12. Oncol Rep (2012) 28, 127-132. [DOI: 10.3892/or.2012.1775](https://doi.org/10.3892/or.2012.1775)
13. Resp Res (2012) 13, 73. [doi:10.1186/1465-9921-13-73](https://doi.org/10.1186/1465-9921-13-73)
14. Antimicrob Agents Chemother (2013) 57(1), 326 –332. [doi:10.1128/AAC.01366-12](https://doi.org/10.1128/AAC.01366-12)
15. BBRC (2013) 442, 33–37. <http://dx.doi.org/10.1016/j.bbrc.2013.10.164>
16. Cell Death Dis (2013) 4, e783. [doi:10.1038/cddis.2013.314](https://doi.org/10.1038/cddis.2013.314)
17. Cell Mol Neurobiol (2013) 33, 625–635. [DOI 10.1007/s10571-013-9921-z](https://doi.org/10.1007/s10571-013-9921-z)
18. Mol Neurobiol (2013) 48, 257–269. [DOI 10.1007/s12035-013-8458-y](https://doi.org/10.1007/s12035-013-8458-y)
19. Mol Vision (2013) 19, 25-38. [PMCID: PMC3541045](https://pubmed.ncbi.nlm.nih.gov/PMC3541045/)
20. Retina (2013) 33(10), 2155-2161. [doi: 10.1097/IAE.0b013e31828b91e6](https://doi.org/10.1097/IAE.0b013e31828b91e6)
21. BMC Comp Alt Med (2014) 14, 370. [doi:10.1186/1472-6882-14-370](https://doi.org/10.1186/1472-6882-14-370)
22. Br J Ophthalmol (2014) 98, i11–i16. [doi:10.1136/bjophthalmol-2014-305302](https://doi.org/10.1136/bjophthalmol-2014-305302)
23. Cell Death Dis (2014) 5, e1150. [doi:10.1038/cddis.2014.53](https://doi.org/10.1038/cddis.2014.53)
24. Indian J Ophthal (2014) 62(1), 16–22. [doi:10.4103/0301-4738.126168](https://doi.org/10.4103/0301-4738.126168)
25. PLoS ONE (2014) 9(4), e95500. [doi:10.1371/journal.pone.0095500](https://doi.org/10.1371/journal.pone.0095500)
26. Trans Am Ophthalmol Soc (2014) 112, 116-141. [PMCID: PMC4311675](https://pubmed.ncbi.nlm.nih.gov/PMC4311675/)
27. ACS Chem Neurosci (2015) 6(4), 588–598. [doi:10.1021/cn500323q](https://doi.org/10.1021/cn500323q)
28. Clin Exp Ophthalmol (2015) 43, 820–829. [doi: 10.1111/ceo.12568](https://doi.org/10.1111/ceo.12568)
29. Sci Rep (2015) 5, 15967. [DOI: 10.1038/srep15967](https://doi.org/10.1038/srep15967)

30. Biomaterials (2016) 76, 115-132. <https://doi.org/10.1016/j.biomaterials.2015.10.044>
31. Carbohydrate Polymers (2016) 146, 187-196. <https://doi.org/10.1016/j.carbpol.2016.03.043>
32. Cell Transplantation (2016) 25, 1967–1977. DOI: <http://dx.doi.org/10.3727/096368916X691457>
33. J Gen Virol (2016) 97, 3183–3192. DOI [10.1099/jgv.0.000625](https://doi.org/10.1099/jgv.0.000625)
34. J Ocular Pharmacol Ther (2016) 32(2), 102-108. DOI: [10.1089/jop.2015.0083](https://doi.org/10.1089/jop.2015.0083)
35. Mol Med Rep (2016) 13, 4461-4468. DOI: [10.3892/mmr.2016.5073](https://doi.org/10.3892/mmr.2016.5073)
36. Oncotarget (2016) 7(51), 84989-84998. doi: [10.18632/oncotarget.13356](https://doi.org/10.18632/oncotarget.13356)
37. Sci Rep (2016) 6, 19720. DOI: [10.1038/srep19720](https://doi.org/10.1038/srep19720)
38. Virology J (2016) 13, 26. DOI [10.1186/s12985-016-0480-z](https://doi.org/10.1186/s12985-016-0480-z)
39. British J Anaesth (2017) 119(3), 481–91. doi: [10.1093/bja/aex071](https://doi.org/10.1093/bja/aex071)
40. J Histochem Cytochem (2017) 65(9), 543–558. <https://doi.org/10.1089/jop.2016.0129>
41. J Ocular Pharmacol Ther (2017) 33(8), 610-619. doi: [10.1089/jop.2016.0129](https://doi.org/10.1089/jop.2016.0129)
42. Plasma Process Polym (2017) 14, e1600162. DOI [10.1002/ppap.201600162](https://doi.org/10.1002/ppap.201600162)
43. Tissue Cell (2017) 49(5), 537-544. <https://doi.org/10.1016/j.tice.2017.06.005>
44. World J Gastroenterol (2017) 23(5), 830-841. DOI: [10.3748/wjg.v23.i5.830](https://doi.org/10.3748/wjg.v23.i5.830)
45. Biomed Sci Letters (2018) 24, 15-22. <https://doi.org/10.15616/BSL.2018.24.1.15>
46. Cancer Sci (2018) 109, 699–709. DOI: [10.1111/cas.13482](https://doi.org/10.1111/cas.13482)
47. Int J Biol Sci (2018) 14(13), 1834-1844. doi: [10.7150/ijbs.27640](https://doi.org/10.7150/ijbs.27640)
48. Molecular Cell (2018) 72(3), 413-425. <https://doi.org/10.1016/j.molcel.2018.08.040>
49. Bioscience Reports (2019) 39, BSR20190069. <https://doi.org/10.1042/BSR20190069>
50. Canadian J Animal Sci (2019) 99(3), 456-464. <https://doi.org/10.1139/cjas-2018-0103>
51. Cancers (2019) 11(1), 64. <https://doi.org/10.3390/cancers11010064>
52. Cells (2019) 8(5), 454. <https://doi.org/10.3390/cells8050454>
53. Front Cell Dev Biol (2019) 7, 396. doi: [10.3389/fcell.2019.00396](https://doi.org/10.3389/fcell.2019.00396)
54. Fungal Biology (2019) <https://doi.org/10.1016/j.funbio.2019.09.011>
55. J Cell Physiol (2020) 235(2), 1531-1542. <https://doi.org/10.1002/jcp.29072>