

RedDot™2 References
Click a blue link to open article

1. BMC Evolutionary Biol (2013) 13, 230. <https://doi.org/10.1186/1471-2148-13-230> (whole tardigrade staining)
2. Environ Mol Mutagen (2013) 54, 737–746. <https://doi.org/10.1002/em.21817> (In Cell Western®)
3. Human Gene Ther (2013) 24(3), 259-269. <http://doi.org/10.1089/hum.2012.247>
4. PLoS ONE (2013) 8(10), e77157. <https://doi.org/10.1371/journal.pone.0077157> (In Cell Western®)
5. PLoS Pathog (2013) 9(7), e1003437. <https://doi.org/10.1371/journal.ppat.1003437>
6. Cellular Microbiol. (2015) 17(3), 303-317. <https://doi.org/10.1111/cmi.12375>
7. Cytometry Part A (2014) 85A, 179-187. <https://doi.org/10.1002/cyto.a.22410> (dead cell staining)
8. Nat Neurosci (2014) 17, 822–831. <https://doi.org/10.1038/nn.3721> (real-time imaging dead cells)
9. PLoS Pathog (2014) 10(5), e1004162. [doi: 10.1371/journal.ppat.1004162](https://doi.org/10.1371/journal.ppat.1004162)
10. Biochem Biophys Acta (2015) 1853, 1749-1758. <https://doi.org/10.1016/j.bbamcr.2015.04.009>
11. Cell Signal (2015) 27(5), 989-96. <https://doi.org/10.1016/j.cellsig.2015.02.009>
12. Dev Biol (2015) 407, 57-67. <https://doi.org/10.1016/j.ydbio.2015.08.008>
13. Infect Immun (2015) 83(9), 3704-11. [doi: 10.1128/IAI.00716-15](https://doi.org/10.1128/IAI.00716-15) (In Cell Western®)
14. J Biomed Mater Res A (2015) 103(6), 2193-7. <https://doi.org/10.1002/jbm.a.35352>
15. J Comp Neurol (2015) 523(13), 1865-1885. <https://doi.org/10.1002/cne.23767>
16. Mol Neurodegen (2015) 10, 57. <https://doi.org/10.1186/s13024-015-0053-4>
17. Nat Commun (2015) 6, 7222. [doi: 10.1038/ncomms8222](https://doi.org/10.1038/ncomms8222)
18. Arch Toxicol (2016) 90, 1983–1995. <https://doi.org/10.1007/s00204-015-1599-1>
19. Histopathology (2016) 69(6), 985-992. <https://doi.org/10.1111/his.13022> (500-1000 um tissue sections, solvent clearing)
20. J Neurotrauma (2016) 33(2), 183-93. <https://doi.org/10.1089/neu.2015.3899>
21. Modern Pathology (2016) 29, 557–569. <https://doi.org/10.1038/modpathol.2016.52>
22. Mutagenesis (2016) 31(1) 83–96. <https://doi.org/10.1093/mutage/gev058> (In Cell Western®)
23. Free Radical Biol Med (2016) 93, 155-164. <https://doi.org/10.1016/j.freeradbiomed.2016.01.028> (In Cell Western®)
24. Acta Neuropathol Commun (2017) 5, 81. <https://doi.org/10.1186/s40478-017-0481-1>
25. Cell Physiol Biochem (2017) 42,2281–2294. <https://doi.org/10.1159/000480001>
26. Cell Rep (2017) 20(1), 236-250. <http://dx.doi.org/10.1016/j.celrep.2017.06.010> (whole organ/whole body CUBIC-cancer analysis)

27. mBio (2017) 8, e00007-17. <https://doi.org/10.1128/mBio.00007-17> (In Cell Western®)
28. Mol Carcinog (2017) 56(3), 1117-1126. <https://doi.org/10.1002/mc.22578>
29. Oncotarget (2017) Jun 8(24), 38581-38591. <https://doi.org/10.18632/oncotarget.16934>
30. Pathogens Disease (2017) 75, ftx066. <https://doi.org/10.1093/femspd/ftx066>
31. Bio-protocol (2018) 8(6), e2771. [DOI: 10.21769/BioProtoc.2771](https://doi.org/10.21769/BioProtoc.2771)
32. Cell Rep (2018) 24, 2196–2210. <https://doi.org/10.1016/j.celrep.2018.07.056> (CUBIC)
33. Environ Mol Mutagen (2018) 59(6), 516-528. <https://doi.org/10.1002/em.22197>
34. Mol Cell Neurosci (2018) 92, 149-163. [doi: 10.1016/j.mcn.2018.08.004](https://doi.org/10.1016/j.mcn.2018.08.004)
35. Stem Cell Research (2018), <https://doi.org/10.1016/j.scr.2018.08.006>
36. Toxicol Letters (2018) 287, 100–107. <https://doi.org/10.1016/j.toxlet.2018.02.007> (In Cell Western®)
37. J Med Food (2019) 22(12), 1280-1293. [http://doi.org/10.1089/jmf.2019.0032](https://doi.org/10.1089/jmf.2019.0032) (In Cell Western®)
38. Nat Protoc (2019) 14, 3506–3537. <https://doi.org/10.1038/s41596-019-0240-9> (whole organ, CUBIC)
39. Particle Fibre Toxicol (2020) 17, 11. <https://doi.org/10.1186/s12989-020-00342-6>
40. Cell Metabolism (2020). <https://doi.org/10.1016/j.cmet.2020.09.006> (whole mouse lung, CUBIC)

In Cell Western is a registered trademark of LI-COR Inc.