

Forget-Me-Not™ EvaGreen® Master Mix References

1. Int J Antimicrob Agents (2017) 50, 572–580. <https://doi.org/10.1016/j.ijantimicag.2017.04.007>
2. J Biol Chem (2017) 292(50), 20494–20508. DOI 10.1074/jbc.M117.807677
3. Immunity (2018) 49, 1090–1102. <https://doi.org/10.1016/j.immuni.2018.10.009>
4. Oncotarget (2018) 9(17), 13244–13253. doi: 10.18632/oncotarget.24284
5. Sci Rep (2018) 8, 13152. DOI:10.1038/s41598-018-31409-2
6. Biophys J (2019) 116(10), 1803-1814. <https://doi.org/10.1016/j.bpj.2019.04.011>
7. bioRxiv (2019) doi: <https://doi.org/10.1101/702076>
8. Environ Sci Technol (2019) 53, 3599–3609. DOI: 10.1021/acs.est.9b00798
9. Gut Microbes (2019) 10(4), 439-446. DOI: 10.1080/19490976.2018.1554962
10. Int J Mol Sci (2019) 20(8), 1955. doi:10.3390/ijms20081955
11. J Apicultural Res (2019) DOI: 10.1080/00218839.2019.1676999
12. J Virol (2019) 93(20) e00364-19. DOI: 10.1128/JVI.00364-19
13. microPublication Biology (2019). <https://doi.org/10.17912/micropub.biology.000136>
14. Oxidative Med Cellular Longev (2019), 4019619. <https://doi.org/10.1155/2019/4019619>
15. Oxidative Med Cellular Longev (2019), 7189854. <https://doi.org/10.1155/2019/7189854>
16. Parasites Vectors (2019) 12, 523. <https://doi.org/10.1186/s13071-019-3781-4> (HRM)
17. Physiol Mol Biol Plants (2019) 25, 1497–1506. <https://doi.org/10.1007/s12298-019-00707-y>
18. Physiologia Plantarum (2019) 165(2), 356-368. <https://doi.org/10.1111/ppl.12866>
19. Phytomedicine (2019) 57, 282-291. <https://doi.org/10.1016/j.phymed.2018.12.044>
20. Trees (2019) 33, 587–597. <https://doi.org/10.1007/s00468-018-1803-z>
21. Food Chem Toxicol (2020) 138, 111250. <https://doi.org/10.1016/j.fct.2020.111250>
22. Front Cell Infect Microbiol (2020) 9,426. <https://doi.org/10.3389/fcimb.2019.00426>
23. Mol Biol Rep (2020). <https://doi.org/10.1007/s11033-020-05323-8>
24. Mol Genet Genomic Med (2020) 00, e1221. DOI: 10.1002/mgg3.1221 (HRM)
25. Plant PhysiolBiochem (2020). 151, 264-270. <https://doi.org/10.1016/j.plaphy.2020.03.031>
26. Sci Rep (2020) 10, 1551. <https://doi.org/10.1038/s41598-020-58411-x>
27. Ticks Tick-borne Dis (2020) 11, 101312. <https://doi.org/10.1016/j.ttbdis.2019.101312> (HRM)
28. Exp Therapeut Med (2021) 21, 403. DOI: 10.3892/etm.2021.9834 (HRM)