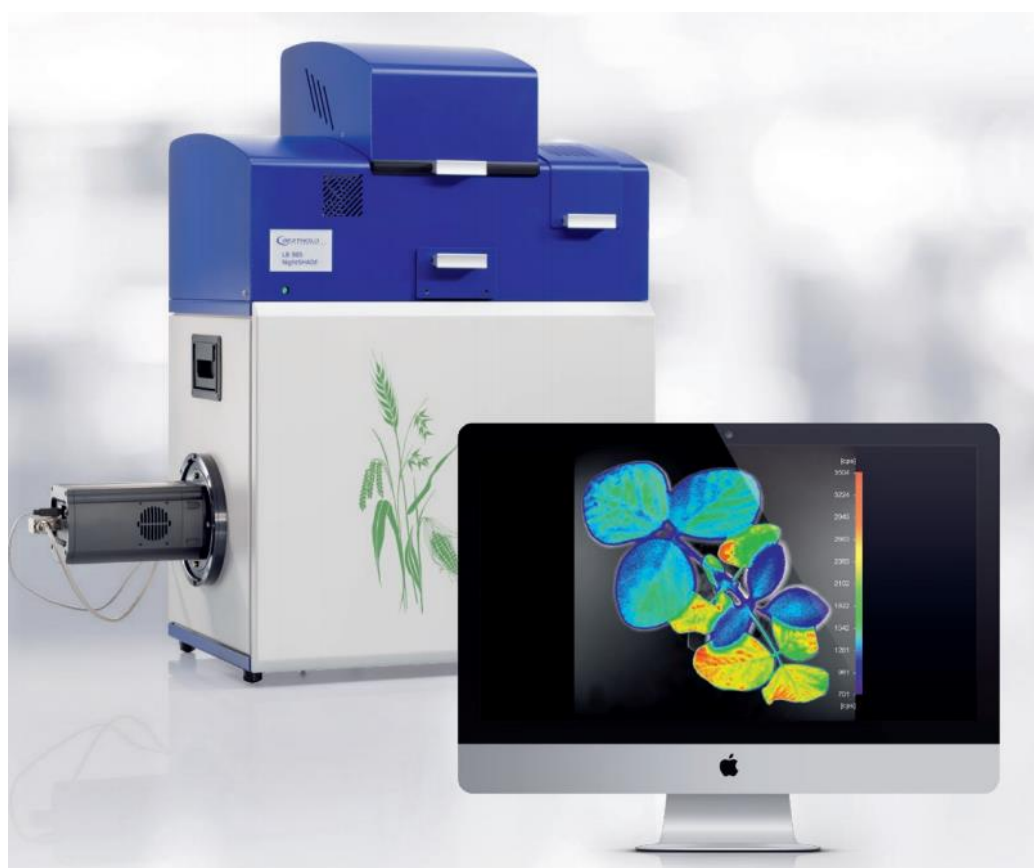


**List of publications using the NightSHADE LB 985,
classified by application**



Updated August 2023

Notes:

1. The **green** font for the journal name indicates its impact factor is over 10.
2. The **red** font for the journal name indicates its impact factor is between 5 and 10.
3. Some older publications use the NightOWL instead of the NightShade.

Signal transduction regulation

1. The D14-SDEL1-SPX4 cascade integrates the strigolactone and phosphate signalling networks in rice. *New Phytologist*, 2023, <https://doi.org/10.1111/nph.18963>
2. Efficient proteome-wide identification of transcription factors targeting Glu-1: A case study for functional validation of TaB3-2A1 in wheat. *Plant Biotechnology Journal*, 2023, <https://doi.org/10.1111/pbi.14103>
3. Transcriptomic and metabolomic analysis reveals a protein module involved in preharvest apple peel browning. *Plant Physiology*, 2023, kiad064, <https://doi.org/10.1093/plphys/kiad0649>
4. The origin and evolution of salicylic acid signaling and biosynthesis in plants. *Molecular Plant*, Volume 16, Issue 1, 2 January 2023, Pages 245-259, <https://doi.org/10.1016/j.molp.2022.12.00>
5. High-Nitrate-Supply-Induced Transcriptional Upregulation of Ascorbic Acid Biosynthetic and Recycling Pathways in Cucumber. *Plants*, 2023, 12(6), 1292; <https://doi.org/10.3390/plants12061292>
6. Kinase MxMPK4-1 and calmodulin binding protein MxIQM3 enhance apple root acidification during Fe deficiency. *Plant Physiology*, 2022, <https://doi.org/10.1093/plphys/kiac587>
7. Transcription Factor CmNAC34 Regulated CmLCYB-Mediated β -Carotene Accumulation during Oriental Melon Fruit Ripening. *International journal of molecular sciences*, 2022, 23(17), 9805; <https://doi.org/10.3390/ijms23179805>
8. Guard cell anion channel PbrSLAC1 regulates stomatal closure through PbrSnRK2.3 protein kinases. *Plant Science*, 2022, Volume 325, 111487, <https://doi.org/10.1016/j.plantsci.2022.111487>
9. The RNA helicase UAP56 and the E3 ubiquitin ligase COP1 coordinately regulate alternative splicing to repress photomorphogenesis in Arabidopsis. *The Plant Cell*, 2022, <https://doi.org/10.1093/plcell/koac235>
10. The Carboxy Terminal Transmembrane Domain of SPL7 Mediates Interaction with RAN1 at the Endoplasmic Reticulum to Regulate Ethylene Signaling in Arabidopsis. *New Phytologist*, 2022, <https://doi.org/10.1111/nph.18376>
11. Identification and Characterization of the Core Region of ZmDi19-5 Promoter Activity and Its Upstream Regulatory Proteins. *International Journal of Molecular Science*, 2022, 23(13), 7390; <https://doi.org/10.3390/ijms23137390>
12. HSFA1 proteins mediate heat-induced accumulation of CPT7-derived polyprenols affecting thylakoid organization. *bioRxiv*, 2021, <https://doi.org/10.1101/2021.12.22.473876>
13. The blue light receptor CRY1 interacts with GID1 and DELLA proteins to repress gibberellin signaling and plant growth. *Plant Communications*, Volume 2, Issue 6, 8 November 2021, 100245, <https://doi.org/10.1016/j.xplc.2021.100245>
14. A reciprocal inhibitory module for Pi and iron signaling. *Molecular Plant*, Volume 15, Issue 1, 3 January 2022, Pages 138-150, <https://doi.org/10.1016/j.molp.2021.09.011>
15. The role of light-harvesting complex I in excitation energy transfer from LHClI to photosystem I in Arabidopsis. *Plant Physiology*, Volume 188, Issue 4, April 2022, Pages 2241–2252, <https://doi.org/10.1093/plphys/kiab579>
16. Phytochrome B Conveys Low Ambient Temperature Cues to the Ethylene-Mediated Leaf Senescence in Arabidopsis. *Plant Cell Physiology*, 2022, 63(3): 326–339, <https://doi.org/10.1093/pcp/pcab178>
17. Plants utilise ancient conserved peptide upstream open reading frames in stress-responsive translational

- regulation. *Plant Cell Environment*, 2022 Feb 6, <https://doi.org/10.1111/pce.14277>.
18. A transcriptional complex of FtMYB102 and FtbHLH4 coordinately regulates the accumulation of rutin in *Fagopyrum tataricum*. *bioRxiv*, 2022, <https://doi.org/10.1101/2022.05.02.490379>
 19. Phosphorylation of the plasma membrane H⁺-ATPase AHA2 by BAK1 is required for ABA-induced stomatal closure in Arabidopsis. *The Plant Cell*, 2022, <https://doi.org/10.1093/plcell/koac106>
 20. The CDC48 complex mediates ubiquitin-dependent degradation of intra-chloroplast proteins in plants. *Cell reports*, VOLUME 39, ISSUE 2, 110664, APRIL 12, 2022, <https://doi.org/10.1016/j.celrep.2022.110664>
 21. Inducible epigenome editing probes for the role of histone H3K4 methylation in Arabidopsis heat stress memory. *Plant Physiology*, 2022 Mar 14; <https://doi.org/10.1093/plphys/kiac113>
 22. Integrated transcriptome, small RNA and degradome analysis provide insights into the transcriptional regulatory networks underlying cold acclimation in jojoba. *Scientia Horticulturae*, Volume 299, 1 June 2022, 111050, <https://doi.org/10.1016/j.scienta.2022.111050>
 23. Mobility of FLOWERING LOCUS T protein as a systemic signal in trifoliolate orange and its low accumulation in grafted juvenile scions. *Horticulture Research*, 2022, <https://doi.org/10.1093/hr/uhac056>
 24. Kn1 participates in spindle assembly checkpoint signaling in maize. *PNAS*, May 18, 2021 118 (20) e2022357118; <https://doi.org/10.1073/pnas.2022357118>
 25. A transcriptional regulatory module controls lipid accumulation in soybean. *New phytologist*, 231:661–678, <http://doi.org/10.1111/nph.17401>
 26. IAA3-Mediated repression of PIF proteins coordinates light and auxin signaling in Arabidopsis. *PLOS Genetics*, 2021 Feb 18;17(2):e1009384, <https://doi.org/10.1371/journal.pgen.1009384>
 27. A Defense Pathway Linking Plasma Membrane and Chloroplasts and Co-opted by Pathogens. *Cell*, 2020 Sep 3;182(5):1109–1124.e25, <https://doi.org/10.1016/j.cell.2020.07.020>
 28. GOLVEN peptide signalling through RGI receptors and MPK6 restricts asymmetric cell division during lateral root initiation. *Nature Plants*, volume 6, pages533–543(2020), <https://doi.org/10.1038/s41477-020-0645-z>
 29. Genome-wide MNase hypersensitivity assay unveils distinct classes of open chromatin associated with H3K27me3 and DNA methylation in Arabidopsis thaliana. *Genome Biology*, volume 21, Article number: 24 (2020), <https://doi.org/10.1186/s13059-020-1927-5>
 30. Salicylic Acid Suppresses Apical Hook Formation via NPR1-Mediated Repression of EIN3 and EIL1 in Arabidopsis. *The Plant Cell*, 2020 Mar;32(3):612–629, <https://doi.org/10.1105/tpc.19.00658>
 31. The plasma-membrane polyamine transporter PUT3 is regulated by the Na⁺/H⁺ antiporter SOS1 and protein kinase SOS2. *New phytologist*, Volume226, Issue3, May 2020, 785–797, <https://doi.org/10.1111/nph.16407>
 32. Photoexcited phytochrome B interacts with brassinazole resistant 1 to repress brassinosteroid signaling in Arabidopsis. *Journal of Integrative Plant Biology*, 2020; 62 (5) ; 652, <https://doi.org/10.1111/jipb.12822>
 33. 植物转录因子与 DNA 互作研究技术. *植物学报 Chinese Bulletin of Botany*, 2020, 55 (4): 468–474.
 34. Uptake of graphene enhanced the photophosphorylation performed by chloroplasts in rice plants. *Nano Research*, volume 13, pages 3198–3205 (2020), <https://doi.org/10.1007/s12274-020-2862-1>
 35. Two RING-Finger Ubiquitin E3 Ligases Regulate the Degradation of SPX4, An Internal Phosphate Sensor, for Phosphate Homeostasis and Signaling in Rice. *Molecular Plant*, Volume 12, Issue 8, 5 August 2019, Pages 1060–1074.
 36. Proteasome subunit RPT2a promotes PTGS through repressing RNA quality control in Arabidopsis. *Nature Plants*, volume 5, pages1273–1282(2019).
 37. The Blue-Light Receptor CRY1 Interacts with BZR1 and BIN2 to Modulate the Phosphorylation and Nuclear Function of BZR1 in Repressing BR Signaling in Arabidopsis. *Molecular Plant*, Volume 12, Issue 5, 6 May 2019, Pages 689–703.

38. Identification of Novel Inhibitors of Auxin-Induced Ca²⁺ Signaling via a Plant-Based Chemical Screen. *Plant Physiology*, May 2019, Vol. 180, pp. 480–496.
39. Gateway-compatible inducible vector set for the functional analysis of transcription factors in plants. *Planta*, May 2018, Volume 247, Issue 5, pp 1261–1266.
40. Phosphorylation and negative regulation of CONSTITUTIVELY PHOTOMORPHOGENIC 1 by PINOID in Arabidopsis. *PNAS*, June 20, 2017 114 (25) 6617-6622.
41. Salicylic Acid-Dependent Plant Stress Signaling via Mitochondrial Succinate Dehydrogenase. *Plant Physiology*, April 2017, Vol. 173, pp. 2029–2040.
42. Phosphorylation and negative regulation of CONSTITUTIVELY PHOTOMORPHOGENIC 1 by PINOID in Arabidopsis. *PNAS*, June 20, 2017 114 (25) 6617-6622.
43. AKIN10, a representative Arabidopsis SNF1-related protein kinase 1 (SnRK1), phosphorylates and downregulates plant HMG-CoA reductase. *FEBS Letter*, Volume 591, Issue 8, April 2017, Pages 1159-1166.
44. The Transcription Factor MYB29 Is a Regulator of ALTERNATIVE OXIDASE1a. *Plant Physiology*, March 2017, Vol. 173, pp. 1824–1843.
45. Optimization of the pollen-tube pathway method of plant transformation using the Yellow Cameleon 3.6 calcium sensor in *Solanum lycopersicum*. *Biologia*, 72/10: 1147–1155, 2017.
46. ESCRT-I Component VPS23A Affects ABA Signaling by Recognizing ABA Receptors for Endosomal Degradation. *Molecular Plant*, Volume 9, Issue 12, 5 December 2016, Pages 1570-1582.
47. A Small G Protein as a Novel Component of the Rice Brassinosteroid Signal Transduction. *Molecular Plant*, Volume 9, Issue 9, 6 September 2016, Pages 1260-1271.
48. A Functional Antagonistic Relationship between Auxin and Mitochondrial Retrograde Signaling Regulates Alternative Oxidase1a Expression in Arabidopsis. *Plant Physiology*, July 2014, Vol. 165, pp. 1233–1254.
49. Epigenetic reprogramming that prevents transgenerational inheritance of the vernalized state. *Nature*. volume 515, pages 587–590(2014).
50. Arabidopsis DELLA and Two HD-ZIP Transcription Factors Regulate GA Signaling in the Epidermis through the L1 Box cis-Element. *The Plant Cell*, Vol. 26: 2905–2919, July 2014.
51. Long-Term In Vivo Imaging of Luciferase-Based Reporter Gene Expression in Arabidopsis Roots. *Root Development* pp 177-190.
52. Temporal-Spatial Interaction between Reactive Oxygen Species and Abscisic Acid Regulates Rapid Systemic Acclimation in Plants. *The Plant Cell*, Vol. 25: 3553–3569.
53. The Screening for Novel Inhibitors of Auxin-Induced Ca²⁺ Signaling. *Plant Chemical Genomics*. pp 89-98
54. Mitochondrial localization and putative signaling function of sucrose synthase in maize. *Journal of Biological Chemistry*, VOL. 281, NO. 23, pp. 15625–15635, June 9, 2006.

Studies on circadian rhythms

1. ABCB-mediated shootward auxin transport feeds into the root clock. *EMBO Reports*, (2023)24:e56271 <https://doi.org/10.15252/embr.202256271>
2. The Na⁺/H⁺ antiporter SALT OVERLY SENSITIVE 1 regulates salt compensation of circadian rhythms by stabilizing GIGANTEA in Arabidopsis. *PNAS*, 2022 Aug 16; 119(33):e2207275119. <https://doi.org/10.1073/pnas.2207275119>
3. Universal Stress Protein (USP) regulates the circadian rhythm of central oscillator genes in Arabidopsis. *FEBS Letters*, 2022, <https://doi.org/10.1002/1873-3468.14410>.

4. An auxin-regulable oscillatory circuit drives the root clock in Arabidopsis. *Science Advances*, 01 Jan 2021: Vol. 7, no. 1, <https://doi.org/10.1126/sciadv.abd4722>
5. Magnesium maintains the length of the circadian period in Arabidopsis. *Plant Physiology*, 2021: 0: 1-14, <https://doi.org/10.1101/2020.05.14.096537>
6. Arabidopsis sirtuins and poly ADP ribose polymerases regulate gene expression in the day but do not affect circadian rhythms. *Plant, Cell & Environment*, 2021 Jan 19. <https://doi.org/10.1111/pce.13996>
7. Transcription Factors FHY3 and FAR1 Regulate Light-induced CIRCADIAN CLOCK ASSOCIATED1 Gene Expression in Arabidopsis. *The Plant Cell*, 2020 May;32(5):1464-1478, <https://doi.org/10.1105/tpc.19.00981>
8. Magnesium regulates circadian period in Arabidopsis. *bioRxiv*, 2020/5/14, <https://doi.org/10.1101/2020.05.14.096537>
9. The Rice Circadian Clock Regulates Tiller Growth and Panicle Development Through Strigolactone Signaling and Sugar Sensing. *The Plant Cell*, August 13, 2020, <https://doi.org/10.1105/tpc.20.00289>
10. The central circadian clock proteins CCA1 and LHY regulate iron homeostasis in Arabidopsis. *Journal of Integrative Plant Biology*, 2019 Feb;61(2):168-181. <https://doi.org/10.1111/jipb.12696>
11. LHY2 Integrates Night-Length Information to Determine Timing of Poplar Photoperiodic Growth. *Current Biology*, 2019 Jul 22;29(14):2402-2406.
12. Circadian Entrainment in Arabidopsis by the Sugar-Responsive Transcription Factor bZIP63. *Current Biology*. 28, 2597–2606, August 20, 2018.
13. BIG Regulates Dynamic Adjustment of Circadian Period in Arabidopsis thaliana. *Plant Physiology*, sept 2018, vol 178, 358-371.
14. Sucrose and Ethylene Signaling Interact to Modulate the Circadian Clock. *Plant Physiology*, 2017 Oct;175(2):947-958.
15. Adjustment of the Arabidopsis circadian oscillator by sugar signalling dictates the regulation of starch metabolism. *Scientific REPORTS*, 2017 Aug 16;7(1):8305.
16. Modeling the photoperiodic entrainment of the plant circadian clock. *Journal of Theoretical Biology*, volume 420, 7 May 2017, Pages 220-231.
17. Kinetics of the LOV domain of ZEITLUPE determine its circadian function in Arabidopsis. *eLife*, <https://doi.org/10.7554/eLife.21646.001>
18. Assessing the Impact of Photosynthetic Sugars on the Arabidopsis Circadian Clock. *Environmental Responses in Plants*, pp 133-140.
19. Photosynthetic entrainment of the Arabidopsis thaliana circadian clock. *Nature*. volume 502, pages 689–692(2013)
20. Arabidopsis FHY3 Specifically Gates Phytochrome Signaling to the Circadian Clock. *The Plant Cell*, Vol. 18, 2506–2516, October 2006.
21. Ambient temperature response establishes ELF3 as a required component of the core Arabidopsis circadian clock. *PNAS*, February 16, 2010 107 (7) 3257-3262.
22. Real-Time Reporting of Circadian-Regulated Gene Expression by Luciferase Imaging in Plants and Mammalian Cells. *Methods in Enzymology*, Volume 393, 2005, Pages 269-288.
23. Independent Roles for EARLY FLOWERING 3 and ZEITLUPE in the Control of Circadian Timing, Hypocotyl Length, and Flowering Time. *Plant Physiology*, November 2005, Vol. 139, pp. 1557–1569.
24. The F-Box Protein ZEITLUPE Confers Dosage-Dependent Control on the Circadian Clock, Photomorphogenesis, and Flowering Time. *The Plant Cell*, Vol. 16, 769–782, March 2004.
25. Signs of the time: environmental input to the circadian clock. *Journal of Experimental Botany*, Volume 53, Issue 374, 1 July 2002, Pages 1535–1550.

26. A Role for LKP2 in the Circadian Clock of Arabidopsis. *The Plant Cell*, Vol. 13, 2659-2670, December 2001.

Gene Expression Studies

1. Characterization and Functional Explorations of O-glycosylation Enzymes SECRET AGENT and SPINDLY in *Pyrus bretschneideri*. *Journal of Plant Biology*, 2022-1-4, [10.1007/s12374-021-09345-y](https://doi.org/10.1007/s12374-021-09345-y)
2. Rapid Validation of Transcriptional Enhancers Using a Transient Reporter Assay. *Modeling Transcriptional Regulation* 2021, pp 253-25, https://doi.org/10.1007/978-1-0716-1534-8_16
3. Downregulation of Polyamine and Diamine Oxidases in Silicon-Treated Cucumber. *Plants*, 2021, 10(6), 1248, <https://doi.org/10.3390/plants10061248>
4. Enhancer-mediated reporter gene expression in *Arabidopsis thaliana*: a forward genetic screen. *The Plant Journal*, Pub Date: 2021-02-06, <https://doi.org/10.1111/tbj.15189>
5. Combining GAL4 GFP enhancer trap with split luciferase to measure spatiotemporal promoter activity in *Arabidopsis*. *The Plant Journal*, Volume102, Issue1, April 2020, Pages 187-198, <https://doi.org/10.1111/tbj.14603>
6. The VvSUPERMAN-like Gene is Differentially Expressed Between Bicarpellate and Tricarpellate Florets of *Vitis vinifera* L. cv. 'Xiangfei' and its Heterologous Expression Reduces Carpel Number in Tomato. *Plant and Cell Physiology*, Volume 61, Issue 10, October 2020, Pages 1760–1774, <https://doi.org/10.1093/pcp/pcaa103>
7. Differential Expression of Fungal Genes Determines the Lifestyle of *Plectosphaerella* Strains During *Arabidopsis thaliana* Colonization. *Molecular Plant*, 2020, vol. 33, No. 11, pp. 1299-1314, <https://doi.org/10.1094/mpmi-03-20-0057-r>
8. Introns mediate post-transcriptional enhancement of nuclear gene expression in the green microalga *Chlamydomonas reinhardtii*. *PLOS Genetics*, 2020 Jul; 16(7): e1008944, <https://doi.org/10.1371/journal.pgen.1008944>
9. Differential Expression of Fungal Genes Determines the Lifestyle of *Plectosphaerella* Strains During *Arabidopsis thaliana* Colonization. *Molecular Plant-microbe Interactions*, 1 Oct 2020, <https://doi.org/10.1094/MPMI-03-20-0057-R>
10. The light-induced transcription factor FtMYB116 promotes accumulation of rutin in *Fagopyrum tataricum*. *Plant, Cell & Environment*, 2019,42(4), 1340-1351.
11. Engineering transkingdom signalling in plants to control gene expression in rhizosphere bacteria. *Nature Communications*, volume 10, Article number: 3430 (2019).
12. Benefits of using genomic insulators flanking transgenes to increase expression and avoid positional effects. *Scientific Reports*, 9, Article number: 8474 (2019).
13. Overexpression and VIGS system for functional gene validation in oriental melon (*Cucumis melo* var. *makuwa* Makino). *Plant Cell, Tissue and Organ Culture (PCTOC)*, volume 137, pages 275–284(2019).
14. Effect of transcription terminator usage on the establishment of transgene transcriptional gene silencing. *BMC Res. Notes*, 2018,11,511
15. Engineered Fusion Proteins for Efficient Protein Secretion and Purification of a Human Growth Factor from the Green Microalga *Chlamydomonas reinhardtii*. *ACS Synth. Biol.*, 2018, 7, 11, 2547–2557.
16. Expression of BpIAA10 from *Betula platyphylla* (birch) is differentially regulated by different hormones and light intensities. *Plant Cell, Tissue and Organ Culture (PCTOC)*. 2018, Volume 132, Issue 2, pp 371–381.
17. Analysis of the promoter features of BpCUC2 in *Betula platyphylla* × *Betula pendula*. *Plant Cell, Tissue and Organ Culture (PCTOC)*. 2018, Volume 132, Issue 1, pp 191–199.

18. Arabidopsis SWI/SNF chromatin remodeling complex binds both promoters and terminators to regulate gene expression. *Nucleic Acids Research*, 2017, Vol. 45, No. 6, 3116–3129.
19. Mediator subunit MED25 links the jasmonate receptor to transcriptionally active chromatin. *PNAS*, published ahead of print October 2, 2017.
20. miR156-Targeted SBP-Box Transcription Factors Interact with DWARF53 to Regulate TEOSINTE BRANCHED1 and BARREN STALK1 Expression in Bread Wheat. *Plant Physiology*, July 2017, Vol. 174, pp. 1931–1948.
21. miR156-regulated TaSPLs interact with TaD53 to regulate TaTB1 and TaBA1 expression in bread wheat. *Plant Physiology*, Published May 2017. DOI: <https://doi.org/10.1104/pp.17.00445>
22. Identification of novel transcriptional regulators of Zat12 using comprehensive yeast one-hybrid screens. *Physiologia Plantarum*. 2016; volume 157;422-441
23. Functional conservation and divergence of GmCHLI genes in polyploid soybean. *The Plant Journal*, (2016) 88, 584–596.
24. Comparison of Intact Arabidopsis thaliana Leaf Transcript Profiles during Treatment with Inhibitors of Mitochondrial Electron Transport and TCA Cycle. *PLOS ONE*, September 2012 | Volume 7 | Issue 9 | e44339
25. BSCTV C2 Attenuates the Degradation of SAMDC1 to Suppress DNA Methylation-Mediated Gene Silencing in Arabidopsis. *The Plant Cell*, Vol. 23: 273–288, January 2011.
26. The gene encoding the catalytically inactive α -amylase BAM4 involved in starch breakdown in Arabidopsis leaves is expressed preferentially in vascular tissues in source and sink organs. *Journal of Plant Physiology*, 167 (2010) 890–895.
27. Altered interactions within FY/AtCPSF complexes required for Arabidopsis FCA-mediated chromatin silencing. *PNAS*, May 26, 2009, 106 (21) 8772-8777.

Stress resistance research

1. The TaWRKY13–TaNHX2 pathway positively regulates cadmium tolerance by inhibiting the expression of TaHMA2. *Environmental and Experimental Botany*, Volume 211, July 2023, 105378, <https://doi.org/10.1016/j.envexpbot.2023.105378>
2. Allelic variation in transcription factor PtoWRKY68 contributes to drought tolerance in Populus. *Plant Physiology*, 2023, <https://doi.org/10.1093/plphys/kiad315>
3. An ethylene-induced NAC transcription factor acts as a multiple abiotic stress responder in conifer. *Horticulture Research*, 2023, <https://doi.org/10.1093/hr/uhad130>
4. The TabHLH094–TaMYC8 complex mediates the cadmium response in wheat. *Molecular Breeding*, volume 43, Article number: 57 (2023), <https://doi.org/10.1007/s11032-023-01404-1>
5. TaTIP41 and TaTAP46 positively regulate drought tolerance in wheat by inhibiting PP2A activity. *Journal of Integrative Plant Biology*, 2023, <https://doi.org/10.1111/jipb.13542>
6. Thermosensitive SUMOylation of TaHsfA1 defines a dynamic ON/OFF molecular switch for the heat stress response in wheat. *The Plant Cell*, 2023, <https://doi.org/10.1093/plcell/koad192>
7. Wall-associated kinase BrWAK1 confers resistance to downy mildew in Brassica rapa. *Plant Biotechnology Journal*, 2023, <https://doi.org/10.1111/pbi.14118>
8. Light and jasmonic acid coordinately regulate the phosphateresponses under shade and phosphate starvation conditions in Arabidopsis. *Plant Direct*, 2023, <https://doi.org/10.1002/pld3.504>
9. Populus euphratica GLABRA3 Binds PLD δ Promoters to Enhance Salt Tolerance. *International journal of*

- molecular sciences*, 2023, 24(9), 8208; <https://doi.org/10.3390/ijms24098208>
10. Kinase MxMPK4-1 and calmodulin-binding protein MxIQM3 enhance apple root acidification during Fe deficiency. *Plant Physiology*, Volume 191, Issue 3, March 2023, Pages 1968–1984, <https://doi.org/10.1093/plphys/kiac587>
 11. Growth-regulating factor 15-mediated gene regulatory network enhances salt tolerance in poplar. *Plant Physiology*, Volume 191, Issue 4, April 2023, Pages 2367–2384, <https://doi.org/10.1093/plphys/kiac600>
 12. Promoter-pervasive transcription causes RNA polymerase II pausing to boost DOG1 expression in response to salt. *The EMBO Journal*, (2023)42:e112443, <https://doi.org/10.15252/embj.2022112443>
 13. Tetratricopeptide-containing SMALL KERNEL 11 is essential for the assembly of cytochrome c oxidase in maize mitochondria. *Plant Physiology*, Volume 192, Issue 1, May 2023, Pages 170–187, <https://doi.org/10.1093/plphys/kiad062>
 14. Pepper clade A PP2C, CaSIP1, negatively modulates drought resistance by suppressing CaSnRK2.6 kinase activity. *Environmental and Experimental Botany*, Volume 209, May 2023, 105275, <http://doi.org/10.1016/j.envexpbot.2023.105275>
 15. The basic helix-loop-helix transcription factor gene, OsHHLH38, plays a key role in controlling rice salt tolerance. *Journal of Integrative Plant Biology*, 2023, <https://doi.org/10.1111/jipb.13489>
 16. Combined lncRNA and mRNA Expression Profiles Identified the lncRNA–miRNA–mRNA Modules Regulating the Cold Stress Response in *Ammopiptanthus nanus*. *International journal of molecular sciences*, 2023, 24(7), 6502; <https://doi.org/10.3390/ijms24076502>
 17. Calcium-dependent activation of CPK12 facilitates its cytoplasm-to-nucleus translocation to potentiate plant hypoxia sensing by phosphorylating ERF-VII transcription factors. *Molecular Plant*, 2023, <https://doi.org/10.1016/j.molp.2023.04.002>
 18. Pepper homeobox abscisic acid signalling-related transcription factor 1, CaHAT1, plays a positive role in drought response. *Plant Cell Environment*, 2023, <https://doi.org/10.1111/pce.14597>
 19. Global crotonylatome and GWAS revealed a TaSRT1-TaPGK model regulating wheat cold tolerance through mediating pyruvate. *Science Advances*, 2023, Vol 9, Issue 19, <https://doi.org/10.1126/sciadv.adg1012>
 20. Functional characterization of dehydrins CpRAB and CpERD and their roles in regulating cold resistance of zucchini fruit under high relative humidity storage. *Postharvest Biology and Technology*, 2023, Volume 202, August 2023, 112387, <https://doi.org/10.1016/j.postharvbio.2023.112387>
 21. Natural variation in Beauty Mark is associated with UV-based geographical adaptation in *Gossypium* species. *BMC Biology*, volume 21, Article number: 106 (2023), <https://doi.org/10.1186/s12915-023-01591-5>
 22. Plant-specific histone deacetylases associate with ARGONAUTE4 to promote heterochromatin stabilization and plant heat tolerance. *New Phytologist*, 2023, <https://doi.org/10.1111/nph.18729>
 23. Promoter-pervasive transcription causes RNA polymerase II pausing to boost DOG1 expression in response to salt. *The EMBO Journal*, 2023, <https://doi.org/10.15252/embj.2022112443>
 24. Involvement of SISTOP1 regulated SIFDH expression in aluminum tolerance by reducing NAD⁺ to NADH in the tomato root apex. *The Plant Journal*, 2022, <https://doi.org/10.1111/tpj.16054>
 25. Genome-wide identification of WOX family members in nine Rosaceae species and a functional analysis of MdWOX13-1 in drought resistance. *Plant Science*, Available online 19 December 2022, 111564, <https://doi.org/10.1016/j.plantsci.2022.111564>
 26. Thioredoxin h2 inhibits MPK5-MPK3 cascade to regulate the CBF-COR signaling pathway in *Citrullus lanatus* suffering chilling stress. *Horticulture Research*, 2022, <https://doi.org/10.1093/hr/uhac256>
 27. MdWRKY39 negatively regulates apple phosphorus-deficiency tolerance by inhibiting MdPHT1;7 expression. *Scientia Horticulturae*, Volume 310, 2023, 111715, <https://doi.org/10.1016/j.scienta.2022.111715>
 28. Transcription factors ABF4 and ABR1 synergistically regulate amylase-mediated starch catabolism in drought

- tolerance. *Plant Physiology*, 2022 Sep 14, <https://doi.org/10.1093/plphys/kiac428>
29. HY5-HDA9 orchestrates the transcription of HsfA2 to modulate salt stress response in Arabidopsis. *Journal of Integrative Plant Biology*, 2022, <https://doi.org/10.1111/jipb.13372>
 30. Wheat TaSnRK2.10 phosphorylates TaERD15 and TaENO1 and confers drought tolerance when overexpressed in rice. *Plant Physiology*, 2022, <https://doi.org/10.1093/plphys/kiac523>
 31. MxMMPK4-1 phosphorylates NADPH oxidase to trigger the MxMMPK6-2-MxbHLH104 pathway mediated Fe deficiency responses in apple. *Plant, Cell & Environment*, 2022 June 23, <https://doi.org/10.1111/pce.14384>
 32. Thinopyrum intermedium TiAP1 interacts with a chitin deacetylase from *Blumeria graminis* f. sp. *tritici* and increases the resistance to Bgt in wheat. *Plant Biotechnology Journal*, 2022, <https://doi.org/10.1101/2021.02.08.430348>
 33. Alternative splicing of REGULATOR OF LEAF INCLINATION 1 modulates phosphate starvation signaling and growth in plants. *The Plant Cell*, 2022, <https://doi.org/10.1002/1873-3468.14410>
 34. Ubiquitin Ligase OsRINGzf1 Regulates Drought Resistance by Controlling the Turnover of OsPIP2;1. *Plant Biotechnology Journal*, 2022, May 19, <https://doi.org/10.1111/pbi.13857>
 35. Regulatory interaction of BcWRKY33A and BcHSFA4A promotes salt tolerance in non-heading Chinese cabbage [*Brassica campestris* (syn. *Brassica rapa*) ssp. *chinensis*]. *Horticulture Research*, 2022, uhac113, <https://doi.org/10.1093/hr/uhac113>
 36. The miR164a-NAM3 module confers cold tolerance by inducing ethylene production in tomato. *The Plant Journal*, 2022 May 15. <https://doi.org/10.1111/tpj.15807>
 37. SISnRK2.3 interacts with SISU1 to modulate high temperature tolerance via Abscisic acid (ABA) controlling stomatal movement in tomato. *Plant Science*, 321 (2022) 111305, <https://doi.org/10.2139/ssrn.4041332>
 38. ERF9 of *Poncirus trifoliata* (L.) Raf. undergoes feedback regulation by ethylene and modulates cold tolerance via regulating a glutathione S-transferase U17 gene. *Plant Biotechnology Journal*, 2022 Jan;20(1):183-200. <https://doi.org/10.1111/pbi.13705>
 39. Interaction of wheat methionine sulfoxide reductase TaMSRB5.2 with glutathione S-transferase TaGSTF3-A contributes to seedling osmotic stress resistance. *Environmental and Experimental Botany*, Volume 194, February 2022, 104731, <https://doi.org/10.1016/j.envexpbot.2021.104731>
 40. SIFHY3 and SIHY5 act compliantly to enhance cold tolerance through the integration of myo-inositol and light signaling in tomato. *New phytologist*, 2022, 233: 2127–2143, <https://doi.org/10.1111/nph.17934>
 41. The miR157-SPL-CNR module acts upstream of bHLH101 to negatively regulate iron deficiency responses in tomato. *Journal of Integrative Plant Biology*, 2022, <https://doi.org/10.1111/jipb.13251>
 42. Sorting Nexin1 negatively modulates phosphate uptake by facilitating Phosphate Transporter1;1 degradation in Arabidopsis. *The Plant Journal*, 2022, <https://doi.org/10.1111/tpj.15778>
 43. The soybean PLATZ transcription factor GmPLATZ17 suppresses drought tolerance by interfering with stress-associated gene regulation of GmDREB5. *The Crop Journal*, <https://doi.org/10.1016/j.cj.2022.03.009>
 44. CsbZIP50 binds to the G-box/ABRE motif in CsRD29A promoter to enhance drought tolerance in cucumber. *Environmental and Experimental Botany*, Volume 199, July 2022, 104884, <https://doi.org/10.1016/j.envexpbot.2022.104884>
 45. Modelling biophoton emission kinetics based on the initial intensity value in *Helianthus annuus* plants exposed to different types of stress. *Scientific Reports*, (2022) 12:2317 <https://doi.org/10.1038/s41598-022-06323-3>
 46. PtrMYB3, a R2R3-MYB Transcription Factor from *Poncirus trifoliata*, Negatively Regulates Salt Tolerance and Hydrogen Peroxide Scavenging. *Antioxidants*, 2021, 10(9), 1388; <https://doi.org/10.3390/antiox10091388>
 47. Nuclear factor Y subunit GmNFYA competes with GmHDA13 for interaction with GmFVE to positively regulate salt tolerance in soybean. *Plant Biotechnology Journal*, Volume19, Issue11, November 2021, Pages 2362-2379,

- <https://doi.org/10.1111/pbi.13668>
48. ThHSFA1 Confers Salt Stress Tolerance through Modulation of Reactive Oxygen Species Scavenging by Directly Regulating ThWRKY4. *International journal of molecular sciences*, 2021, 22, 5048, <https://doi.org/10.3390/ijms22095048>.
 49. The NF-Y-PYR module integrates the abscisic acid signal pathway to regulate plant stress tolerance. *Plant Biotechnology Journal*, 2021 Dec; 19(12): 2589–2605, <https://doi.org/10.1111/pbi.13684>
 50. PagWOX11/12a activates PagCYP736A12 gene that facilitates salt tolerance in poplar. *Plant Biotechnology Journal*, 2021 Nov; 19(11): 2249–2260, <https://doi.org/10.1111/pbi.13653>
 51. ERF108 from *Poncirus trifoliata* (L.) Raf. functions in cold tolerance by modulating raffinose synthesis through transcriptional regulation of PtrRafS. *The Plant Journal*, Volume108, Issue3, November 2021, Pages 705-724, <https://doi.org/10.1111/tbj.15465>
 52. Consortia of Plant-Growth-Promoting Rhizobacteria Isolated from Halophytes Improve Response of Eight Crops to Soil Salinization and Climate Change Conditions. *Agronomy*, 2021, 11(8), 1609; <https://doi.org/10.3390/agronomy11081609>
 53. Mitochondrial Transcription Termination Factor 27 Is Required for Salt Tolerance in *Arabidopsis thaliana*. *Int. J. Mol. Sci.*, 2021 Feb 2;22(3):1466, <https://doi.org/10.3390/ijms22031466>
 54. The TAZ domain-containing proteins play important role in the heavy metals stress biology in plants. *Environmental Research*, 24 March 2021, 111030, <https://doi.org/10.1016/j.envres.2021.111030>
 55. ThNAC12 from *Tamarix hispida* directly regulates ThPIP2;5 to enhance salt tolerance by modulating reactive oxygen species. *Plant Physiology and Biochemistry*, Volume 163, June 2021, Pages 27-35, <https://doi.org/10.1016/j.plaphy.2021.03.042>
 56. Transcriptional activation and phosphorylation of OsCNGC9 confer enhanced chilling tolerance in rice. *Molecular Plant*, Volume 14, Issue 2, Pages 315-329, <https://doi.org/10.1016/j.molp.2020.11.022>
 57. TaSnRK2.4 is a vital regulator in control of thousand-kernel weight and response to abiotic stress in wheat. *Journal of Integrative Agriculture*, 2021, 20(1): 46–54, [https://doi.org/10.1016/s2095-3119\(19\)62830-3](https://doi.org/10.1016/s2095-3119(19)62830-3)
 58. Ethylene increases the cold tolerance of apple via the MdERF1B–MdCibHLLH1 regulatory module. *The Plant Journal*, First published: 26 January 2021, <https://doi.org/10.1111/tbj.15170>
 59. Interaction between MdMYB63 and MdERF106 enhances salt tolerance in apple by mediating Na⁺/H⁺ transport. *Plant Physiology and Biochemistry*, Volume 155, October 2020, Pages 464-471, [https://doi.org/10.1016/s2095-3119\(19\)62830-3](https://doi.org/10.1016/s2095-3119(19)62830-3)
 60. OsCYCP4s coordinate phosphate starvation signaling with cell cycle progression in rice. *Journal of Integrative Plant Biology*, Volume62, Issue7, July 2020, Pages 1017-1033, <https://doi.org/10.1111/jipb.12885>
 61. Mutation of DELAYED GREENING1 impairs chloroplast RNA editing at elevated ambient temperature in *Arabidopsis*. *Journal of Genetics and Genomics*, Volume 47, Issue 4, 20 April 2020, Pages 201-212, <https://doi.org/10.1016/j.jgg.2020.03.005>
 62. The *Arabidopsis* phosphatase PP2C49 negatively regulates salt tolerance through inhibition of AtHKT1;1. *Journal of Integrative Plant Biology*, 2020 Sep 2. <https://doi.org/10.1111/jipb.13008>
 63. *Populus euphratica* WRKY1 binds the promoter of H⁺-ATPase gene to enhance gene expression and salt tolerance. *Journal of Experimental Botany*, 2020, 71(4):1527-1539, <https://doi.org/10.1093/jxb/erz493>
 64. ESCRT-I component VPS23A sustains salt tolerance by strengthening the SOS module in *Arabidopsis*. *Molecular Plant*, 18 May 2020, 13(8):1134-1148, <https://doi.org/10.1016/j.molp.2020.05.010>
 65. Efficient expression and function of a receptor-like kinase in wheat powdery mildew defence require an intron-located MYB binding site. *Plant Biotechnology Journal*, 2020 Nov 22. <https://doi.org/10.1111/pbi.13512>
 66. Effect of cadmium stress on certain physiological parameters, antioxidative enzyme activities and biophoton emission of leaves in barley (*Hordeum vulgare* L.) seedlings. *PLoS One*, 2020 Nov 3; 15(11),

<https://doi.org/10.1371/journal.pone.0240470>

67. The barley stripe mosaic virus expression system reveals the wheat C2H2 zinc finger protein TaZFP1B as a key regulator of drought tolerance. *BMC Plant Biology*, (2020) 20:144. <https://doi.org/10.1186/s12870-020-02355-x>
68. Creation two hyperactive variants of phytochrome B1 for attenuating shade avoidance syndrome in maize. *Journal of Integrative Agriculture*, 2020-19-01, [https://doi.org/10.1016/s2095-3119\(20\)63466-9](https://doi.org/10.1016/s2095-3119(20)63466-9)
69. GmNFYA13 improves salt and drought tolerance in transgenic soybean plants. *Frontiers in Plant Science*, 2020 Oct 23;11:587244, <https://doi.org/10.3389/fpls.2020.587244>
70. Oxidative stress level and dehydrin gene expression pattern differentiate two contrasting cucumber F1 hybrids under high fertigation treatment. *International Journal of Biological Macromolecules*, Volume 161, 15 October 2020, Pages 864-874, <https://doi.org/10.1016/j.ijbiomac.2020.06.050>
71. A novel biosensor to monitor proline in pea root exudates and nodules under osmotic stress and recovery. *Plant and Soil*, (2020) 452:413–422, <https://doi.org/10.1007/s11104-020-04577-2>
72. Polyploidy-mediated divergent light-harvesting and photoprotection strategies under temperature stress in a Mediterranean carnation complex. *Environmental and Experimental Botany*, 171 (2020) 103956, <https://doi.org/10.1016/j.envexpbot.2019.103956>
73. GmWRKY54 improves drought tolerance through activating genes in abscisic acid and Ca²⁺ signaling pathways in transgenic soybean. *The Plant Journal*, Volume 100, Issue 2, October 2019, Pages 384-398, <https://doi.org/10.1111/tbj.14449>
74. A stress-responsive bZIP transcription factor OsbZIP62 improves drought and oxidative tolerance in rice. *BMC Plant Biology*, 19, Article number: 260 (2019), <https://doi.org/10.1186/s12870-019-1872-1>
75. INDETERMINATE SPIKELET1 Recruits Histone Deacetylase and a Transcriptional Repression Complex to Regulate Rice Salt Tolerance. *Plant Physiol* 2018, 178 (2), 824-837, <https://doi.org/10.1104/pp.18.00324>
76. The Arabidopsis RNA Polymerase II Carboxyl Terminal Domain (CTD) Phosphatase-Like1 (CPL1) is a biotic stress susceptibility gene. *Scientific Reports*, volume 8, Article number: 13454 (2018).
77. Coordinating the overall stomatal response of plants: Rapid leaf-to-leaf communication during light stress. *Science Signaling*, 20 Feb 2018:Vol. 11, Issue 518, eaam9514.
78. Light and Ethylene Coordinately Regulate the Phosphate Starvation Response through Transcriptional Regulation of PHOSPHATE STARVATION RESPONSE1. *The Plant Cell*, Vol. 29: 2269–2284, September 2017.
79. Antisense transcription represses Arabidopsis seed dormancy QTL DOG1 to regulate drought tolerance. *EMBO reports*, (2017) 18, 2186-2196.
80. Apple MdMYC2 reduces aluminum stress tolerance by directly regulating MdERF3 gene. *Plant and Soil*, 2017, Volume 418, Issue 1–2, pp 255–266.
81. Rapid detoxification via glutathione S-transferase (GST) conjugation confers a high level of atrazine resistance in Palmer amaranth (*Amaranthus palmeri*). *Pest management science*, Volume73, Issue11, November 2017, Pages 2236-2243.
82. ELF3 Controls Thermoresponsive Growth in Arabidopsis. *Current Biology*, 2015, 25, 194–199.
83. Ultra-fast alterations in mRNA levels uncover multiple players in light stress acclimation in plants. *The Plant Journal*, Volume84, Issue4, November 2015, Pages 760-772.
84. Ectopic Overexpression of SsCBF1, a CRT/DRE-Binding Factor from the Nightshade Plant *Solanum lycopersicoides*, Confers Freezing and Salt Tolerance in Transgenic Arabidopsis. *PLOS ONE*, June 2013 | Volume 8 | Issue 6 | e61810.
85. Overexpression of a Maize Transcription Factor ZmPHR1 Improves Shoot Inorganic Phosphate Content and Growth of Arabidopsis under Low-Phosphate Conditions. *Plant Molecular Biology Reporter*, June 2013, Volume 31, Issue 3, pp 665–677.



86. Ectopic Overexpression of SlHsfA3, a Heat Stress Transcription Factor from Tomato, Confers Increased Thermotolerance and Salt Hypersensitivity in Germination in Transgenic Arabidopsis. *PLoS ONE*, January 2013 | Volume 8 | Issue 1 | e54880.
87. Arabidopsis Ubiquitin Conjugase UBC32 Is an ERAD Component That Functions in Brassinosteroid-Mediated Salt Stress Tolerance. *The Plant Cell*, Vol. 24: 233–244, January 2012.