

List of publications using the NightShade LB 985, classified by plant species



Updated May 2021

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.

Arabidopsis-related

1. The Arabidopsis Mediator Complex Subunit 8 Regulates Oxidative Stress Responses. *The Plant Cell*, <https://doi.org/10.1093/plcell/koab079>.
2. Bacterial effector targeting of a plant iron sensor facilitates iron acquisition and pathogen colonization. *The Plant Cell*, <https://doi.org/10.1093/plcell/koab075>.
3. The SNL-HDA19 histone deacetylase complex antagonizes HY5 activity to repress photomorphogenesis in Arabidopsis. *New phytologist*, Volume229, Issue 6, March 2021, Pages 3221-3236.
4. MicroRNA775 regulates intrinsic leaf size and reduces cell wall pectin levels by targeting a galactosyltransferase gene in Arabidopsis. *The Plant Cell*, <https://doi.org/10.1093/plcell/koaa049>
5. BIC 1 acts as a transcriptional coactivator to promote brassinosteroid signaling and plant growth. *EMBO J*, (2021)40:e104615. <https://doi.org/10.15252/emj.2020104615>
6. An auxin-regulable oscillatory circuit drives the root clock in Arabidopsis. *Science Advances*, 01 Jan 2021: Vol. 7, no. 1, eabd4722.
7. Magnesium maintains the length of the circadian period in Arabidopsis. *Plant Physiology*, 2021: 0: 1-14.
8. 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:10.1111/pce.13996>
9. Enhancer-mediated reporter gene expression in Arabidopsis thaliana: a forward genetic screen. *The Plant Journal*, Pub Date : 2021-02-06 , <https://doi:10.1111/tpj.15189>.
10. Mitochondrial Transcription Termination Factor 27 Is Required for Salt Tolerance in Arabidopsis thaliana. *Int. J. Mol. Sci.*, 2021 Feb 2;22(3):1466.
11. IAA3-Mediated repression of PIF proteins coordinates light and auxin signaling in Arabidopsis. *PLOS Genetics*, 2021 Feb 18;17(2):e1009384.
12. FHY3 interacts with phytochrome B and regulates seed dormancy and germination. *Plant Physiology*, 2021, kiab147, <https://doi.org/10.1093/plphys/kiab147>.
13. The TAZ domain-containing proteins play important role in the heavy metals stress biology in plants. *Environmental Research*, Available online 24 March 2021, 111030.
14. 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.
15. A Defense Pathway Linking Plasma Membrane and Chloroplasts and Co-opted by Pathogens. *Cell*, 2020 Sep 3;182(5):1109-1124.e25.
16. GOLVEN peptide signalling through RGI receptors and MPK6 restricts asymmetric cell division during lateral root initiation. *Nature Plants*, volume 6, pages533–543(2020).
17. The Tetracentron genome provides insight into the early evolution of eudicots and the formation of vessel elements. *Genome Biology*, 2020, volume 21, Article number: 291 (2020).

18. 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).
19. Phosphorylation of ATG18a by BAK1 suppresses autophagy and attenuates plant resistance against necrotrophic pathogens. *Autophagy*, 2020 Aug 26;1-18.
20. 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.
21. Transcription Factors FHY3 and FAR1 Regulate Light-induced CIRCADIAN CLOCK ASSOCIATED1 Gene Expression in *Arabidopsis*. *The Plant Cell*, 2020 May;32(5):1464-1478.
22. FATTY ACID DESATURASE5 is required to induce autoimmune responses in gigantic chloroplast mutants of *Arabidopsis*. *The Plant Cell*, 2020 Oct;32(10):3240-3255.
23. 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.
24. PIF1 and RVE1 form a transcriptional feedback loop to control light-mediated seed germination in *Arabidopsis*. *Journal of Integrative Plant Biology*, 2020, Volume 62, Issue 9, 1372–1384.
25. 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.
26. Photoexcited phytochrome B interacts with brassinazole resistant 1 to repress brassinosteroid signaling in *Arabidopsis*. *Journal of Integrative Plant Biology*, 2020; 62 (5) ; 652.
27. SIZ1-Mediated SUMO Modification of SEUSS Regulates Photomorphogenesis in *Arabidopsis*. *Plant Communications*, 2020 Jun 2;1(5):100080.
28. 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>. Online ahead of print.
29. Combining GAL4 GFP enhancer trap with split luciferase to measure spatiotemporal promoter activity in *Arabidopsis*. *The Plant Journal*, Volume102, Issue 1, April 2020, Pages 187-198.
30. A bacterial effector protein prevents MAPK-mediated phosphorylation of SGT1 to suppress plant immunity. *Plos Pathogens*, 2020 Sep 25;16(9):e1008933.
31. ABCB-mediated auxin transport in outer root tissues regulates lateral root spacing in *Arabidopsis*. *bioRxiv*, <https://doi.org/10.1101/2020.07.22.206300>.
32. MicroRNA775 Promotes Intrinsic Leaf Size and Reduces Cell Wall Pectin Level via a Target Galactosyltransferase in *Arabidopsis*. *bioRxiv*, doi: <https://doi.org/10.1101/2020.09.17.301705>.
33. *Populus euphratica* WRKY1 binds the promoter of H⁺-ATPase gene to enhance gene expression and salt tolerance. *Journal of Experimental Botany*, 01 Feb 2020, 71(4):1527-1539.
34. Magnesium regulates circadian period in *Arabidopsis*. *bioRxiv*, 2020/5/14.
35. 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.
36. The Evening Complex and the Chromatin-Remodeling Factor PICKLE Coordinately Control Seed Dormancy by Directly Repressing DOG1 in *Arabidopsis*. *Plant Communications*, 2019 Dec 3;1(2):100011.
37. ESCRT-I component VPS23A sustains salt tolerance by strengthening the SOS module in *Arabidopsis*. *Molecular Plant*, 18 May 2020, 13(8):1134-1148.
38. Proteasome subunit RPT2a promotes PTGS through repressing RNA quality control in *Arabidopsis*. *Nature Plants*, volume 5, pages1273–1282(2019).
39. *Arabidopsis* FHY3 and FAR1 Regulate the Balance between Growth and Defense Responses under Shade Conditions. *The Plant Cell*, 2019 Sep; 31(9): 2089–2106.
40. 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.

41. Identification of Novel Inhibitors of Auxin-Induced Ca²⁺ Signaling via a Plant-Based Chemical Screen. *Plant Physiology*, May 2019, Vol. 180, pp. 480–496.
42. The B3-Domain Transcription Factor VAL1 Regulates the Floral Transition by Repressing FLOWERING LOCUS T. *Plant Physiology*, September 2019, Vol. 181, pp. 236–248.
43. The central circadian clock proteins CCA1 and LHY regulate iron homeostasis in Arabidopsis. *Journal of Integrative Plant Biology*, 2019 Feb;61(2):168-181.
44. Chromatin-remodeling factor PICKLE antagonizes polycomb repression of FT to promote flowering. *Plant Physiology*, 2019 Oct;181(2):656-668.
45. Lateral Inhibition by a Peptide Hormone-Receptor Cascade during Arabidopsis Lateral Root Founder Cell Formation. *Developmental Cell*, 2019 Jan 7;48(1):64-75.e5.
46. B-BOX DOMAIN PROTEIN28 Negatively Regulates Photomorphogenesis by Repressing the Activity of Transcription Factor HY5 and Undergoes COP1-Mediated Degradation. *The Plant Cell*, Vol. 30: 2006–2019, September 2018.
47. Benefits of using genomic insulators flanking transgenes to increase expression and avoid positional effects. *Scientific Reports*, 9, Article number: 8474 (2019).
48. A Regulatory Module Controlling GA-Mediated Endosperm Cell Expansion Is Critical for Seed Germination in Arabidopsis. *Molecular Plant*, P1–15, 2018
49. A novel pathway linking plasma membrane and chloroplasts is co-opted by pathogens to suppress salicylic acid-dependent defences. *bioRxiv*, doi: <https://doi.org/10.1101/837955>.
50. Integrated Regulation of Apical Hook Development by Transcriptional Coupling of EIN3/EIL1 and PIFs in Arabidopsis. *The Plant Cell*, 2018 Sep;30(9):1971-1988.
51. Circadian Entrainment in Arabidopsis by the Sugar-Responsive Transcription Factor bZIP63. *Current Biology*. 28, 2597–2606, August 20, 2018.
52. BIG Regulates Dynamic Adjustment of Circadian Period in Arabidopsis thaliana. *Plant Physiology*, sept 2018, vol 178, 358-371.
53. Gateway-compatible inducible vector set for the functional analysis of transcription factors in plants. *Planta*, May 2018, Volume 247, Issue 5, pp 1261–1266.
54. Synthetic hormone-responsive transcription factors can monitor and reprogram plant development. *eLife*, 2018;7:e34702.
55. The Arabidopsis thaliana Mediator subunit MED8 regulates plant immunity to Botrytis Cinerea through interacting with the basic helix-loop-helix (bHLH) transcription factor FAMA. *PLoS ONE*, 2018,13(3),e0193458.
56. 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).
57. BRASSINOSTEROID-SIGNALING KINASE1 Phosphorylates MAPKKK5 to Regulate Immunity in Arabidopsis. *Plant Physiol.*, Published April 2018. DOI: <https://doi.org/10.1104/pp.17.01757>.
58. Reporter-based forward genetic screen to identify bundle sheath anatomy mutants in A. thaliana. *The Plant Journal*, First published: 17 November 2018.
59. STERILE APETALA modulates the stability of a repressor protein complex to control organ size in Arabidopsis thaliana. *PLoS Genet*, 2018, 14(2): e1007218
60. Elongator is required for root stem cell maintenance by regulating SHORT ROOT transcription. *Plant Physiology*, Published on November 6, 2018, as DOI:10.1104/pp.18.00534.
61. 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.
62. Dynamic regulation of PIF5 by COP1–SPA complex to optimize photomorphogenesis in Arabidopsis. *The Plant Journal*, 2018,96,260-273.

63. Effect of transcription terminator usage on the establishment of transgene transcriptional gene silencing. *BMC Res. Notes*, 2018,11,511
64. 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.
65. Phosphorylation and negative regulation of CONSTITUTIVELY PHOTOMORPHOGENIC 1 by PINOID in Arabidopsis. *PNAS*, June 20, 2017 114 (25) 6617-6622.
66. Light-Dependent Degradation of PIF3 by SCFEBF1/2 Promotes a Photomorphogenic Response in Arabidopsis. *Current Biology*, 27, 2420–2430, August 21, 2017.
67. Sucrose and Ethylene Signaling Interact to Modulate the Circadian Clock. *Plant Physiology*, 2017 Oct;175(2):947-958.
68. MicroRNA319-regulated TCPs interact with FBHs and PFT1 to activate CO transcription and control flowering time in Arabidopsis. *PLoS Genetics*, 2017, 13(5): e1006833.
69. The GYF domain protein PSIG1 dampens the induction of cell death during plant-pathogen interactions. *PLoS Genetics*, 2017 Oct 26;13(10):e1007037.
70. Salicylic Acid-Dependent Plant Stress Signaling via Mitochondrial Succinate Dehydrogenase. *Plant Physiology*, April 2017, Vol. 173, pp. 2029–2040.
71. Adjustment of the Arabidopsis circadian oscillator by sugar signaling dictates the regulation of starch metabolism. *Scientific REPORTS*, 2017 Aug 16;7(1):8305.
72. Phosphorylation and negative regulation of CONSTITUTIVELY PHOTOMORPHOGENIC 1 by PINOID in Arabidopsis. *PNAS*, June 20, 2017 114 (25) 6617-6622.
73. 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.
74. Spatial Auxin Signaling Controls Leaf Flattening in Arabidopsis. *Current Biology*, 27, 2940–2950, October 9, 2017
75. Modeling the photoperiodic entrainment of the plant circadian clock. *Journal of Theoretical Biology*, Volume 420, 7 May 2017, Pages 220-231.
76. Kinetics of the LOV domain of ZEITLUPE determine its circadian function in Arabidopsis. *eLife*, <https://doi.org/10.7554/eLife.21646.001>.
77. The Arabidopsis NADPH oxidases RbohD and RbohF display differential expression patterns and contributions during plant immunity. *Journal of Experimental Botany*, Volume 67, Issue 6, 1 March 2016, Pages 1663–1676.
78. Control of seed dormancy in Arabidopsis by a cis-acting noncoding antisense transcript. *PNAS*. November 29, 2016 113 (48) E7846-E7855.
79. Mediator subunit MED25 links the jasmonate receptor to transcriptionally active chromatin. *PNAS*, published ahead of print October 2, 2017.
80. Antisense transcription represses Arabidopsis seed dormancy QTL DOG1 to regulate drought tolerance. *EMBO reports*, (2017) 18, 2186-2196.
81. AKIN10, a representative Arabidopsis SNF1-related protein kinase 1 (SnRK1), phosphorylates and downregulates plant HMG-CoA reductase. *FEBS Letter*, Volume591, Issue8, April 2017, Pages 1159-1166.
82. The Transcription Factor MYB29 Is a Regulator of ALTERNATIVE OXIDASE1a. *Plant Physiology*, March 2017, Vol. 173, pp. 1824–1843.
83. 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.
84. Identification of novel transcriptional regulators of Zat12 using comprehensive yeast one-hybrid screens. *Physiologia Plantarum*. 2016; volume 157;422-441
85. Phosphorylation of CBP20 Links MicroRNA to Root Growth in the Ethylene Response. *PLoS Genet.*, 2016.12(11): e1006437.

86. The Arabidopsis NADPH oxidases RbohD and RbohF display differential expression patterns and contributions during plant immunity. *Journal of Experimental Botany*, Vol. 67, No. 6 pp. 1663–1676, 2016.
87. Jasmonate-Activated MYC2 Represses ETHYLENE INSENSITIVE3 Activity to Antagonize Ethylene Promoted Apical Hook Formation in Arabidopsis. *The Plant Cell*, Vol. 26: 1105–1117, March 2014.
88. A SUMO Ligase AtMMS21 Regulates the Stability of the Chromatin Remodeler BRAHMA in Root Development. *Plant Physiology*, March 2017, Vol. 173, pp. 1574–1582.
89. A Protein-Based Genetic Screening Uncovers Mutants Involved in Phytochrome Signaling in Arabidopsis. *Front. Plant Sci.*, 22 July 2016 | <https://doi.org/10.3389/fpls.2016.01086>
90. FHY3 and FAR1 Act Downstream of Light Stable Phytochromes. *Front Plant Sci.* 22 February 2016 | <https://doi.org/10.3389/fpls.2016.00175>.
91. Chloroplast retrograde signal regulates flowering. *PNAS*, September 20, 2016 113 (38) 10708-10713.
92. Transcriptional Mechanism of Jasmonate Receptor COI1-Mediated Delay of Flowering Time in Arabidopsis. *The Plant Cell*, Vol. 27: 2814–2828, October 2015.
93. ELF3 Controls Thermoresponsive Growth in Arabidopsis. *Current Biology*, 2015, 25, 194–199.
94. 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.
95. Arabidopsis DELLA and Two HD-ZIP Transcription Factors Regulate GA Signaling in the Epidermis through the L1 Box cis-Element. *The Plant Cell*, Vol. 27: 2437–2454, September 2015.
96. Red Light-Mediated Degradation of CONSTANS by the E3 Ubiquitin Ligase HOS1 Regulates Photoperiodic Flowering in Arabidopsis. *The Plant Cell*, Vol. 27: 2437–2454, September 2015.
97. Redox rhythm reinforces the circadian clock to gate immune response. *Nature*, Vol. 523, 472-476, June 2015. <https://doi.org/10.1038/nature14449>
98. A reversible decrease in ribulose 1,5-bisphosphate carboxylase/oxygenase carboxylation activity caused by the aggregation of the enzyme's large subunit is triggered in response to the exposure of moderate irradiance-grown plants to low irradiance. *Physiologia Plantarum*, 154: 591–608. 2015.
99. Arabidopsis DAYU/ABERRANT PEROXISOME MORPHOLOGY9 Is a Key Regulator of Peroxisome Biogenesis and Plays Critical Roles during Pollen Maturation and Germination in Planta. *The Plant Cell*, Published February 2014. DOI: <https://doi.org/10.1105/tpc.113.121087>.
100. Polycomb-Group Proteins and FLOWERING LOCUS T Maintain Commitment to Flowering in Arabidopsis thaliana. *The Plant Cell*, Vol. 26: 2457–2471, June 2014.
101. Assessing the Impact of Photosynthetic Sugars on the Arabidopsis Circadian Clock. *Environmental Responses in Plants*. pp 133-140.
102. Overexpression of a Novel Component Induces HAK5 and Enhances Growth in Arabidopsis. *IRSN Botany*, volume 2014, Article ID 490252.
103. 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.
104. Epigenetic reprogramming that prevents transgenerational inheritance of the vernalized state. *Nature*. volume 515, pages 587–590(2014).
105. Photosynthetic entrainment of the Arabidopsis thaliana circadian clock. *Nature*. volume 502, pages689–692(2013)
106. 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.
107. Long-Term In Vivo Imaging of Luciferase-Based Reporter Gene Expression in Arabidopsis Roots. *Root Development* pp 177-190. (Protocol)
108. Cesium Inhibits Plant Growth through Jasmonate Signaling in Arabidopsis thaliana. *Int. J. Mol. Sci.* 2013, 14, 4545-4559.

109. Phosphorylation-Coupled Proteolysis of the Transcription Factor MYC2 Is Important for Jasmonate Signaled Plant Immunity. *PLOS Genetics*, April 2013 | Volume 9 | Issue 4 | e1003422.
110. 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.
111. Cyclin-dependent Kinase E1 (CDKE1) Provides a Cellular Switch in Plants between Growth and Stress Responses. *Journal of Biological Chemistry*, 2013, VOL. 288, NO. 5, pp. 3449–3459.
112. Temporal-Spatial Interaction between Reactive Oxygen Species and Abscisic Acid Regulates Rapid Systemic Acclimation in Plants. *The Plant Cell*, Vol. 25: 3553–3569.
113. 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.
114. Feedback inhibition by thiols outranks glutathione depletion: a luciferase-based screen reveals glutathione-deficient γ -ECS and glutathione synthetase mutants impaired in cadmium-induced sulfate assimilation. *The Plant Journal*, (2012) 70, 783–795.
115. Expression of the central growth regulator BIG BROTHER is regulated by multiple cis-elements. *BMC Plant Biology*, 12, Article number: 41 (2012).
116. 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
117. The Basic Helix-Loop-Helix Transcription Factor MYC2 Directly Represses PLETHORA Expression during Jasmonate-Mediated Modulation of the Root Stem Cell Niche in Arabidopsis. *The Plant Cell*, Vol. 23: 3335–3352, September 2011.
118. 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.
119. Extracellular Nucleotides and Apyrases Regulate Stomatal Aperture in Arabidopsis. *Plant Physiology*, August 2011, Vol. 156, pp. 1740–1753.
120. Ambient temperature response establishes ELF3 as a required component of the core Arabidopsis circadian clock. *PNAS*, February 16, 2010 107 (7) 3257-3262.
121. 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.
122. The CURLY LEAF Interacting Protein BLISTER Controls Expression of Polycomb-Group Target Genes and Cellular Differentiation of Arabidopsis thaliana. *The Plant Cell*, Vol. 22: 2291–2305, July 2010.
123. Altered interactions within FY/AtCPSF complexes required for Arabidopsis FCA-mediated chromatin silencing. *PNAS*, May 26, 2009 106 (21) 8772-8777.
124. The Screening for Novel Inhibitors of Auxin-Induced Ca²⁺ Signaling. *Plant Chemical Genomics*. pp 89-98 (book)
125. ARABIDOPSIS TRITHORAX1 Dynamically Regulates FLOWERING LOCUS C Activation via Histone 3 Lysine 4 Trimethylaion. *The Plant Cell*, Vol. 20: 580–588, March 2008.
126. Arabidopsis FHY3 Specifically Gates Phytochrome Signaling to the Circadian Clock. *The Plant Cell*, Vol. 18, 2506–2516, October 2006.
127. 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.
128. Independent Roles for EARLY FLOWERING 3 and ZEITLUPE in the Control of Circadian Timing,

- Hypocotyl Length, and Flowering Tim. *Plant Physiology*, November 2005, Vol. 139, pp. 1557–1569.
129. 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.
130. Signs of the time: environmental input to the circadian clock. *Journal of Experimental Botany*, Volume 53, Issue 374, 1 July 2002, Pages 1535–1550.
131. A Role for LKP2 in the Circadian Clock of Arabidopsis. *The Plant Cell*, Vol. 13, 2659–2670, December 2001.

Rice-related

1. Post-Golgi Trafficking of Rice Storage Proteins Requires the Small GTPase Rab7 Activation Complex MON1-CCZ1. *Plant Physiology*, <https://doi.org/10.1093/plphys/kiab175>.
2. DHD4, a CONSTANS-like family transcription factor, delays heading date by affecting the formation of the FAC complex in rice. *Molecular Plant*, Volume 14, Issue 2, 1 February 2021, Pages 330–343.
3. Transcriptional activation and phosphorylation of OsCNGC9 confer enhanced chilling tolerance in rice. *Molecular Plant*, Volume 14, Issue 2, 1 February 2021, Pages 315–329.
4. The URL1–ROC5–TPL2 transcriptional repressor complex represses the ACL1 gene to modulate leaf rolling in rice. *Plant Physiology*, kiaa121, <https://doi.org/10.1093/plphys/kiaa121>.
5. Uptake of graphene enhanced the photophosphorylation performed by chloroplasts in rice plants. *Nano Research*, volume 13, pages3198–3205(2020).
6. OsmiR530 acts downstream of OsPIL15 to regulate grain yield in rice. *New phytologist*, (2020) 226: 823–837.
7. Rice stripe virus suppresses jasmonic acid-mediated resistance by hijacking brassinosteroid signaling pathway in rice. *Plos Pathogens*, 2020 Aug 31;16(8):e1008801.
8. The Rice Circadian Clock Regulates Tiller Growth and Panicle Development Through Strigolactone Signaling and Sugar Sensing. *The Plant Cell*, August 13, 2020, doi:10.1105/tpc.20.00289
9. OsCYCP4s coordinate phosphate starvation signaling with cell cycle progression in rice. *Journal of Integrative Plant Biology*, Volume62, Issue7, July 2020, Pages 1017–1033.
10. A cyclic nucleotide-gated channel mediates cytoplasmic calcium elevation and disease resistance in rice. *Cell Research*, volume 29, pages820–831(2019).
11. 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.
12. A stress-responsive bZIP transcription factor OsbZIP62 improves drought and oxidative tolerance in rice. *BMC Plant Biology*, 19, Article number: 260 (2019).
13. OsBBX14 promotes photomorphogenesis in rice by activating OsHY5L1 expression under blue light conditions. *Plant Science*, Volume 284, July 2019, Pages 192–202.
14. 基于瞬时表达系统的水稻 miRNA 靶基因快速验证系统的建立. *Biotechnology Bulletin 生物技术通报*, 2019, 35(10): 57–63.
15. INDETERMINATE SPIKELET1 Recruits Histone Deacetylase and a Transcriptional Repression Complex to Regulate Rice Salt Tolerance. *Plant Physiol.* Vol. 178, 2018.
16. Brassinosteroids Regulate OFP1, a DLT Interacting Protein, to Modulate Plant Architecture and Grain Morphology in Rice. *Front. Plant Sci.*, 27 September 2017.
17. 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.

18. Rice Dwarf Virus P2 Protein Hijacks Auxin Signaling by Directly Targeting the Rice OsIAA10 Protein, Enhancing Viral Infection and Disease Development. *Plos Pathogens*, Published: September 8, 2016 <https://doi.org/10.1371/journal.ppat.1005847>.
19. The Starch Granule-Associated Protein EARLY STARVATION1 Is Required for the Control of Starch Degradation in Arabidopsis thaliana Leaves. *The Plant Cell*, Vol. 28: 1472–1489, June 2016.
20. Retention of OsNMD3 in the cytoplasm disturbs protein synthesis efficiency and affects plant development in rice. *Journal of Experimental Botany*, Volume 65, Issue 12, 1 July 2014, Pages 3055–3069.
21. Leaf Rolling Controlled by the Homeodomain Leucine Zipper Class IV Gene Roc5 in Rice. *Plant Physiology*, July 2011, Vol. 156, pp. 1589–1602.
22. Quantitative Measurements of Xanthomonas Oryzae pv. Oryzae Distribution in Rice Using Fluorescent-Labeling. *Journal of Plant Biology*, August 2011, 54:269.
23. Functional analysis of the structural domain of ARF proteins in rice (Oryza sativa L.). *Journal of Experimental Botany*, Vol. 61, No. 14, pp. 3971–3981, 2010.

Wheat- and Barley-related

1. 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.
2. Thinopyrum intermedium TiAP1 interacts with a chitin deacetylase from Blumeria graminis f. sp. tritici and increases the resistance to Bgt in wheat. bioRxiv, doi: <https://doi.org/10.1101/2021.02.08.430348>.
3. 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>.
4. 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): e0240470.
5. TaDA1, a conserved negative regulator of kernel size, has an additive effect with TaGW2 in common wheat (Triticum aestivum L.). *Plant Biotechnology Journal*, 2020 May;18(5):1330-1342.
6. 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>
7. JAZ proteins modulate seed germination through interaction with ABI5 in bread wheat and Arabidopsis. *New phytologist*, (2019) 223: 246–260.
8. Comparative transcriptome and histological analyses of wheat in response to phytotoxic aphid Schizaphis graminum and non-phytotoxic aphid Sitobion avenae feeding. *BMC Plant Biology*, volume 19, Article number: 547 (2019).
9. Heterodera avenae GLAND5 Effector Interacts with Pyruvate Dehydrogenase Subunit of Plant to Promote Nematode Parasitism. *Frontiers in Microbiology*. 10:1241. <https://doi.org/10.3389/fmicb.2019.01241>.
10. TaCOLD1 defines a new regulator of plant height in bread wheat. *Plant Biotechnology Journal*, 2019 Mar; 17(3): 687–699.
11. TaZIM-A1 negatively regulates flowering time in common wheat (Triticum aestivum L.). *Journal of Integrative Plant Biology*, March 2019 | Volume 61 | Issue 3 | 359–376.
12. 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.

13. 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>
14. A rapid and efficient method for uniform gene expression using the barley stripe mosaic virus. *Plant Methods*, (2017) 13:24.
15. A rapid and efficient method for uniform gene expression using the barley stripe mosaic virus. *Plant Methods*, 13, Article number: 24 (2017)
16. Metabolism of ¹⁴C-labeled polychlorinated biphenyl congeners by wheat cell suspension cultures. *Journal of Environmental Science and Health, Part B. Pesticides, Food Contaminants, and Agricultural Wastes*. Volume 50, 2015 - Issue 2, Pages 69-80.

Maize-related

1. Characterization of novel regulatory modules controlling leaf angle in maize. *bioRxiv*, doi: <https://doi.org/10.1101/2021.02.02.429433>.
2. Genome-wide selection and genetic improvement during modern maize breeding. *Nature Genetics*, volume 52, pages565–571(2020).
3. Genome Editing Enables Next-Generation Hybrid Seed Production Technology. *Molecular Plant*. 2020,13,1-8.
4. CRISPR/Cas9-mediated knockout and overexpression studies reveal a role of maize phytochrome C in regulating flowering time and plant height. *Plant Biotechnology Journal*, 2020 Dec;18(12):2520-2532.
5. Nuclear-Encoded Maturase Protein 3 is Required for the Splicing of Various Group II Introns in Mitochondria during Maize (*zea mays* L.) Seed Development. *Plant and Cell Physiology*, 2020 Dec 30;pcaa161.
6. Creation two hyperactive variants of phytochrome B1 for attenuating shade avoidance syndrome in maize. *Journal of Integrative Agriculture*, 2020-19-01.
7. Characterization of Maize Phytochrome-Interacting Factors in Light Signaling and Photomorphogenesis. *Plant Physiology*, October 2019, Vol. 181, pp. 789–803.
8. Oxalotrophy, a widespread trait of plant-associated Burkholderia species, is involved in successful root colonization of lupin and maize by Burkholderia phytofirmans. *Frontiers in Microbiology*, 09 January 2014 | <https://doi.org/10.3389/fmicb.2013.00421>.
9. 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.
10. Effect of Corn Plant on Survival and Phenanthrene Degradation Capacity of Pseudomonas Sp. UG14Lr in Two Soils. *International Journal of Phytoremediation*, 14:585–595, 2012.
11. 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.

Soybean related

1. Spatial Divergence of PHR-PHT1 Modules Maintains Phosphorus Homeostasis in Soybean Nodules. *Plant Physiology*, 2020 Sep;184(1):236-250.
2. GmGPA3 is involved in post-Golgi trafficking of storage proteins and cell growth in soybean cotyledons. *Plant Science*, Volume 294, May 2020, 110423.
3. GmNFYA13 improves salt and drought tolerance in transgenic soybean plants. *Frontiers in Plant Science*, 2020 Oct 23;11:587244.



4. 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.
5. FT/FD-GRF5 repression loop directs growth to increase soybean yield. *bioRxiv*, October 2018, DOI: 10.1101/437053.
6. Functional conservation and divergence of GmCHLI genes in polyploid soybean. *The Plant Journal*, (2016) 88, 584–596.

Cotton related

1. The Cotton Apoplastic Protein CRR1 Stabilizes Chitinase 28 to Facilitate Defense against the Fungal Pathogen *Verticillium dahliae*. *The Plant Cell*, Vol. 31: 520–536, February 2019.
2. Cotton plant defence against a fungal pathogen is enhanced by expanding BLADE-ON-PETIOLE1 expression beyond lateral-organ boundaries. *Communications Biology*, volume 2, Article number: 238 (2019).
3. Mutation of SELF-PRUNING homologs in cotton promotes short-branching plant architecture. *Journal of Experimental Botany*, Volume 69, Issue 10, 27 April 2018, Pages 2543–2553.

Nicotiana Benthamiana related

1. Carbon-based nanomaterials suppress Tobacco Mosaic Virus (TMV) infection and induce resistance in *Nicotiana benthamiana*. *Journal of Hazardous Materials*, Volume 404, Part A, 15 February 2021, 124167.
2. ASCORBATE PEROXIDASE6 delays the onset of age-dependent leaf senescence. *Plant Physiology*, 2020: 0: 1–16.
3. Rapid validation of transcriptional enhancers using agrobacterium-mediated transient assay. *Plant Methods*, 15, Article number: 21 (2019).
4. Luciferase Complementation Assay for Protein-Protein Interactions in Plants. *Current Protocols in Plant Biology*, 3:42–50, March 2018.
5. A rapid assay to quantify the cleavage efficiency of custom-designed nucleases in plants. *Plant Molecular Biology*, 2013, Volume 82, Issue 3, pp 207–221.

Fruit related

1. Ethylene increases the cold tolerance of apple via the MdERF1B–MdCibHLH1 regulatory module. *The plant Journal*, First published: 26 January 2021 <https://doi.org/10.1111/tpj.15170>
2. A basic/helix–loop–helix transcription factor controls leaf shape by regulating auxin signaling in apple. *New phytologist*. First published: 26 July 2020 <https://doi.org/10.1111/nph.16828>.
3. 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.
4. Acidovorax citrulli Type III Effector AopP Suppresses Plant Immunity by Targeting the Watermelon Transcription Factor WRKY6. *Front Plant Sci.*, 2020 Nov 20;11:579218.
5. ETHYLENE RESPONSE FACTOR39–MYB8 complex regulates low-temperature-induced lignification of loquat fruit. *Journal of Experimental Botany*, Volume 71, Issue 10, 30 May 2020, Pages 3172–3184,
6. High CO₂/hypoxia-induced softening of persimmon fruit is modulated by DkERF8/16 and DkNAC9 complexes. *Journal of Experimental Botany*, Volume 71, Issue 9, 9 May 2020, Pages 2690–2700.

7. High concentrations of CPPU promotes cucurbitacin B accumulation in melon (*Cucumis melo* var. *makuwa* Makino) fruit by inducing transcription factor CmBt. *Plant Physiology and Biochemistry*, 2020 Sep;154:770-781.
8. 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, pages275–284(2019).
9. An effective artificial microRNA vector based on Fv-miR166 precursor from strawberry. *Scientia Horticulturae*, Volume 256, 15 October 2019, 108643.
10. The regulatory module MdPUB29-MdbHLH3 connects ethylene biosynthesis with fruit quality in apple. *New phytologist*, First published: 04 October 2018 <https://doi.org/10.1111/nph.15511>.
11. Ethylene response factors Pp4ERF24 and Pp12ERF96 regulate blue light-induced anthocyanin biosynthesis in 'Red Zaosu' pear fruits by interacting with MYB114. *Plant Molecular Biology*, First Online: 11 December 2018, pp 1–12.
12. The Jasmonate-Activated Transcription Factor MdMYC2 Regulates ETHYLENE RESPONSE FACTOR and Ethylene Biosynthetic Genes to Promote Ethylene Biosynthesis during Apple Fruit Ripening. *The Plant Cell*, Vol. 29: 1316–1334, June 2017.
13. Melon 13-lipoxygenase CmLOX18 may be involved in C6 volatiles biosynthesis in fruit. *Scientific Reports*, volume 7, Article number: 2816 (2017).
14. Apple MdMYC2 reduces aluminum stress tolerance by directly regulating MdERF3 gene. *Plant and Soil*, 2017, Volume 418, Issue 1–2, pp 255–266.
15. High-level expression of a novel chromoplast phosphate transporter CIPHT4;2 is required for flesh color development in watermelon. *New phytologist*, (2017) 213: 1208–1221.
16. Accumulation of Carotenoids and Expression of Carotenoid Biosynthetic Genes during Maturation in Citrus Fruit. *Plant Physiology*, February 2004, Vol. 134, pp. 824–837.

Vegetable related

1. Tomato YABBY2b controls plant height through regulating indole-3-acetic acid-amido synthetase (GH3.8) expression. *Plant Science*. 2020 Aug; 297:110530
2. 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.
3. Tomato Root Transformation Followed by Inoculation with *Ralstonia Solanacearum* for Straightforward Genetic Analysis of Bacterial Wilt Disease. *Journal of Visualized Experiments*, 2020 Mar 11;(157). doi: 10.3791/60302.
4. 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.
5. A novel biosensor to monitor proline in pea root exudates and nodules under osmotic stress and recovery. *Plant and Soil*, (2020) 452:413–422.
6. Identification of a Major QTL (qRRs-10.1) That Confers Resistance to *Ralstonia solanacearum* in Pepper (*Capsicum annuum*) Using SLAF-BSA and QTL Mapping. *Int. J. Mol. Sci.* 2019, 20(23), 5887.
7. Cucumber Mildew Resistance Locus O Interacts with Calmodulin and Regulates Plant Cell Death Associated with Plant Immunity. *Int. J. Mol. Sci.*, 2019, 20(12), 2995.
8. Pan-genome analysis identifies intersecting roles for *Pseudomonas* specialized metabolites in potato pathogen inhibition. *BioRxiv*, 2019, <https://doi.org/10.1101/783258>.

9. The Two Translationally Controlled Tumor Protein Genes, CsTCTP1 and CsTCTP2, Are Negative Modulators in the Cucumis sativus Defense Response to Sphaerotheca fuliginea. *Front Plant Sci.* 2018; 9: 544.
10. Functional Characterization of BoaMYB51s as Central Regulators of Indole Glucosinolate Biosynthesis in Brassica oleracea var. alboglabra Bailey. *Front Plant Sci.*, 2018; 9: 1599.
11. 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.
12. 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.
13. Evaluation of *Ralstonia solanacearum* Infection Dynamics in Resistant and Susceptible Pepper Lines Using Bioluminescence Imaging. *Plant Disease*. 2017. 101(2): 272-278.
14. The detection of hydrogen peroxide involved in plant virus infection by fluorescence spectroscopy. *Luminescence*, Volume31, Issue5, August 2016, Pages 1158-1165.
15. Role of Tomato Lipoxygenase D in Wound-Induced Jasmonate Biosynthesis and Plant Immunity to Insect Herbivores. *PLOS Genetics.*, 2013 | Volume 9 | Issue 12 | e1003964.
16. 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.
17. Evaluation of a Biocontrol Preparation Consisting of *Enterobacter asburiae* JX1 and a Lytic Bacteriophage Cocktail To Suppress the Growth of *Salmonella Javiana* Associated with Tomatoes. *Journal of Food Protection*, Vol. 72, No. 11, 2009, Pages 2284–2292.
18. Development of a novel inducible bioluminescent and antibiotic resistance tagging system and its use to investigate the role of antibiotic production by *Pectobacterium carotovorum* ssp. *carotovorum* during potato tuber infection. *European Journal of Plant Pathology*, volume 125, Article number: 655 (2009)
19. Internalization of *Escherichia coli* O157:H7 following Biological and Mechanical Disruption of Growing Spinach Plants. *Journal of Food Protection*, Vol. 68, No. 12, 2005, Pages 2506–2509.
20. Interaction of *Escherichia coli* with Growing Salad Spinach Plants. *Journal of Food Protection*, Vol. 66, No. 10, 2003, Pages 1790–1797.
21. Internalization of bioluminescent *Escherichia coli* and *Salmonella* Montevideo in growing bean sprouts. *Journal of Applied Microbiology*, 2003, 95, 719–727.

Microalga and Bacteria related

1. High cell density cultivation enables efficient and sustainable recombinant polyamine production in the microalga *Chlamydomonas reinhardtii*. *Bioresource Technology*, Volume 323, March 2021, 124542.
2. Rational Promoter Engineering Enables Robust Terpene Production in Microalgae. *ACS Synthetic Biology*, March 25, 2021, <https://doi.org/10.1021/acssynbio.0c00632>.
3. 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.
4. The *Medicago truncatula* Vacuolar iron Transporter-Like proteins VTL4 and VTL8 deliver iron to symbiotic bacteria at different stages of the infection process. *New phytologist*, 2020, First published: 10 June 2020 <https://doi.org/10.1111/nph.16735>
5. Introns mediate post-transcriptional enhancement of nuclear gene expression in the green microalga *Chlamydomonas reinhardtii*. *PLOS Genetics*, 2020 Jul; 16(7): e1008944.

6. A bacterial effector uncovers a metabolic pathway involved in resistance to bacterial wilt disease. bioRxiv, doi: <https://doi.org/10.1101/2020.09.22.307595>.
7. From macro to micro: a combined bioluminescence-fluorescence approach to monitor bacterial localization. *Environmental Microbiology*, First published: 26 October 2020 <https://doi.org/10.1111/1462-2920.15296>
8. 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>.
9. Engineering transkingdom signalling in plants to control gene expression in rhizosphere bacteria. *Nature Communications*, volume 10, Article number: 3430 (2019).
10. 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.
11. Tailored carbon partitioning for phototrophic production of (E)- α -bisabolene from the green microalga Chlamydomonas reinhardtii. *Metabolic Engineering*. Volume 45, January 2018, Pages 211–222.
12. Eukaryotic microalgae as hosts for light-driven heterologous isoprenoid production. *Planta*, 2019 Jan;249(1):155-180.
13. Inducibility of Tn916 conjugative transfer in Enterococcus faecalis by subinhibitory concentrations of ribosome-targeting antibiotics. *J Antimicrob Chemother.* 2017; 72: 2722–2728.
14. Bacterial Biosensors for in Vivo Spatiotemporal Mapping of Root Secretion. *Plant Physiology*, July 2017, Vol. 174, pp. 1289–1306.
15. Rapid construction and screening of artificial microRNA systems in Chlamydomonas reinhardtii. *The Plant Journal*, (2014) 79, 1052–1064.
16. Robust Expression and Secretion of Xylanase1 in Chlamydomonas reinhardtii by Fusion to a Selection Gene and Processing with the FMDV 2A Peptide. *PLOS ONE*, August 24, 2012. <https://doi.org/10.1371/journal.pone.0043349>.
17. Phylogenetics and Taxonomy of the Fungal Vascular Wilt Pathogen Verticillium, with the Descriptions of Five New Species. *PLoS ONE*, 2011,6(12): e28341.
18. Functional imaging of biophoton responses of plants to fungal infection. *Eur J Plant Pathol*, (2011) 130:249–258.
19. Interactions of Escherichia coli O157:H7, Salmonella typhimurium and Listeria monocytogenes plants cultivated in a gnotobiotic system. *International Journal of Food Microbiology*, Volume 99, Issue 1, 1 March 2005, Pages 7-18.
20. Development of a luciferase reporter gene, luxCt, for Chlamydomonas reinhardtii chloroplast. *The Plant Journal*, 2004,37,449-458.

Others

1. TREE1-EIN3-mediated transcriptional repression inhibits shoot growth in response to ethylene. *PNAS*, November 17, 2020 117 (46) 29178-29189;
2. Ginsenosides regulate adventitious root formation in Panax ginseng via a PgCLE45-PgWOX11 regulatory module. *Journal of Experimental Botany*, Volume 71, Issue 20, 22 October 2020, Pages 6396–6407.
3. Polyploidy-mediated divergent light-harvesting and photoprotection strategies under temperature stress in a Mediterranean carnation complex. *Environmental and Experimental Botany*, 171 (2020) 103956.

4. The florigen interactor BdES43 represses flowering in the model temperate grass *Brachypodium distachyon*. *The Plant Journal*, 2020 Apr;102(2):262-275.
5. N-glucosyltransferase GbNGT1 from Ginkgo complement auxin metabolic pathway. *bioRxiv* preprint posted August 31, 2020.
6. The light-induced transcription factor FtMYB116 promotes accumulation of rutin in *Fagopyrum tataricum*. *Plant, Cell & Environment*, 2019,42(4), 1340-1351.
7. LHY2 Integrates Night-Length Information to Determine Timing of Poplar Photoperiodic Growth. *Current Biology*, 2019 Jul 22;29(14):2402-2406.
8. 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.
9. 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.
10. Phenotypic plasticity of polyploid plant species promotes transgressive behaviour in their hybrids. *AoB PLANTS*. Volume 10, Issue 5, October 2018, ply055.
11. Real-time monitoring of PtaHMGB activity in poplar transactivation assays. *Plant Methods*, 13, Article number: 50 (2017).
12. Enhancement of alkaloid production in opium and California poppy by transactivation using heterologous regulatory factors. *Plant Biotechnology Journal*, (2008) 6, pp. 160–175.