

Select Publications

Amaryllis Nucleics Technology

- [Drobnitch ST, Comas LH, Flynn N, Ibarra Caballero J, Barton RW, et al. \(2021\). Drought-Induced Root Pressure in Sorghum bicolor. *Front Plant Sci* 12: 571072.](#)
- [Boehme JT, Morris CJ, Chiacchia SR, Gong W, Wu KY, et al. \(2021\). HIF-1 \$\alpha\$ promotes cellular growth in lymphatic endothelial cells exposed to chronically elevated pulmonary lymph flow. *Sci Rep* 11\(1\): 1468.](#)
- [Chodasiewicz M, Sokolowska EM, Nelson-Dittrich AC, Masiuk A, Beltran JCM, et al. \(2020\). Identification and Characterization of the Heat-Induced Plastidial Stress Granules Reveal New Insight Into Arabidopsis Stress Response. *Front Plant Sci* 11: 595792.](#)
- [Dew-Budd K, Cheung J, Palos K, Forsythe ES, Beilstein MA \(2020\). Evolutionary and biochemical analyses reveal conservation of the Brassicaceae telomerase ribonucleoprotein complex. *PLoS One* 15\(4\): e0222687.](#)
- [Li C, Nozue K, Maloof JN \(2020\). MYCs and PIFs Act Independently in Arabidopsis Growth Regulation. *G3 \(Bethesda\)*.](#)
- [Traubenik S, Blanco F, Zanetti ME, Reynoso MA \(2020\). TRAP-SEQ of Eukaryotic Translatomes Applied to the Detection of Polysome-Associated Long Noncoding RNAs. *Methods Mol Biol* 2166: 451-472.](#)
- [Shi Y, Hung ST, Rocha G, Lin S, Linares GR, et al. \(2019\). Identification and therapeutic rescue of autophagosome and glutamate receptor defects in C9ORF72 and sporadic ALS neurons. *JCI Insight* 5.](#)
- [Babos KN, Galloway KE, Kisler K, Zitting M, Li Y, et al. \(2019\). Mitigating Antagonism between Transcription and Proliferation Allows Near-Deterministic Cellular Reprogramming. *Cell Stem Cell* 25\(4\): 486-500.e9.](#)
- [Jiang J, Zeng L, Ke H, De La Cruz B, Dehesh K \(2019\). Orthogonal regulation of phytochrome B abundance by stress-specific plastidial retrograde signaling metabolite. *Nat Commun* 10\(1\): 2904.](#)

- [Ito T, Nishio H, Tarutani Y, Emura N, Honjo MN, et al. \(2019\). Seasonal Stability and Dynamics of DNA Methylation in Plants in a Natural Environment. *Genes \(Basel\)* 10\(7\).](#)
- [Warren JG, Mercado J, Grace D \(2019\). Occurrence of Hop Latent Viroid Causing Disease in Cannabis sativa in California. *Plant Disease* 103\(10\): 2699.](#)
- [Wu J, Ichihashi Y, Suzuki T, Shibata A, Shirasu K, et al. \(2019\). Abscisic acid-dependent histone demethylation during postgermination growth arrest in Arabidopsis. *Plant Cell Environ* 42\(7\): 2198-2214.](#)
- [Amini H, Naghavi MR, Shen T, Wang Y, Nasiri J, et al. \(2019\). Tissue-Specific Transcriptome Analysis Reveals Candidate Genes for Terpenoid and Phenylpropanoid Metabolism in the Medicinal Plant Ferula assafoetida. *G3 \(Bethesda\)* 9\(3\): 807-816.](#)
- [Bellmann J, Monette A, Tripathy V, Sójka A, Abo-Rady M, et al. \(2019\). Viral Infections Exacerbate FUS-ALS Phenotypes in iPSC-Derived Spinal Neurons in a Virus Species-Specific Manner. *Front Cell Neurosci* 13: 480.](#)
- [Motomura K, Arae T, Uramoto HA, Suzuki Y, Takeuchi H, et al. \(2019\). AtNOT1 is a novel regulator of gene expression during pollen development. *Plant Cell Physiol*.](#)
- [Hinckley WE, Keymanesh K, Cordova JA, Brusslan JA \(2019\). The HAC1 histone acetyltransferase promotes leaf senescence and regulates the expression of ERF022. *Plant Direct* 3\(8\): e00159.](#)
- [Gaudinier A, Rodriguez-Medina J, Zhang L, Olson A, Liseron-Monfils C, et al. \(2018\). Transcriptional regulation of nitrogen-associated metabolism and growth. *Nature* 563\(7730\): 259-264.](#)
- [Ichihashi Y, Kusano M, Kobayashi M, Suetsugu K, Yoshida S, et al. \(2018\). Transcriptomic and Metabolomic Reprogramming from Roots to Haustoria in the Parasitic Plant, Thesium chinense. *Plant Cell Physiol* 59\(4\): 724-733.](#)
- [Kim JS, Yamaguchi-Shinozaki K, Shinozaki K \(2018\). ER-Anchored Transcription Factors bZIP17 and bZIP28 Regulate Root Elongation. *Plant Physiol* 176\(3\): 2221-2230.](#)
- [Uemura A, Yamaguchi N, Xu Y, Wee W, Ichihashi Y, et al. \(2018\). Regulation of floral meristem activity through the interaction of AGAMOUS, SUPERMAN, and CLAVATA3 in Arabidopsis. *Plant Reprod* 31\(1\): 89-105.](#)

- [Li R, Jeong K, Davis JT, Kim S, Lee S, et al. \(2018\). Integrated QTL and eQTL Mapping Provides Insights and Candidate Genes for Fatty Acid Composition, Flowering Time, and Growth Traits in a F2 Population of a Novel Synthetic Allopolyploid Brassica napus. Front Plant Sci 9: 1632.](#)
- [Ichida JK, Staats KA, Davis-Dusenbery BN, Clement K, Galloway KE, et al. \(2018\). Comparative genomic analysis of embryonic, lineage-converted and stem cell-derived motor neurons. Development 145\(22\).](#)
- [Song YH, Kubota A, Kwon MS, Covington MF, Lee N, et al. \(2018\). Molecular basis of flowering under natural long-day conditions in Arabidopsis. Nat Plants 4\(10\): 824-835.](#)
- [Townsley BT, Covington MF, Ichihashi Y, Zumstein K, Sinha NR \(2015\). BrAD-seq: Breath Adapter Directional sequencing: a streamlined, ultra-simple and fast library preparation protocol for strand specific mRNA library construction. Front Plant Sci 6: 366.](#)
- [Chitwood DH, Kumar R, Ranjan A, Pelletier JM, Townsley BT, et al. \(2015\). Light-Induced Indeterminacy Alters Shade-Avoiding Tomato Leaf Morphology. Plant Physiol 169\(3\): 2030-47.](#)