

Selected Publications Using YourSeq RNA-Seq Kits and Services under Amaryllis Nucleics

- Irulappan V, Kandpal M, Saini K, Rai A, Ranjan A, Sinharoy S, Senthil-Kumar M. **Drought stress exacerbates fungal colonization and endodermal invasion and dampens defense responses to increase dry root rot in chickpea.** Mol Plant Microbe Interact. 2022 Mar 6. [doi: 10.1094/MPMI-07-21-0195-FI](https://doi.org/10.1094/MPMI-07-21-0195-FI)
 - YourSeq RNAseq Library Kit (FT)
 - Chickpea root
- Jathar V, Saini K, Chauhan A, Rani R, Ichihashi Y, Ranjan A. **Spatial control of cell division by GA-OsGRF7/8 module in a leaf explaining the leaf length variation between cultivated and wild rice.** New Phytol. 2022 May;234(3):867-883. [doi: 10.1111/nph.18029](https://doi.org/10.1111/nph.18029).
 - YourSeq RNAseq Library Kit (FT)
 - Rice leaf
- Boehme, J.T., Morris, C.J., Chiacchia, S.R. *et al.* **HIF-1 α promotes cellular growth in lymphatic endothelial cells exposed to chronically elevated pulmonary lymph flow.** *Sci Rep* **11**, 1468 (2021). <https://doi.org/10.1038/s41598-020-80882-1>
 - YourSeq RNAseq Service
 - Primary lymphatic endothelial cells (LECs)
- Hsiao-Pei Yang, Marius Wenzel, Duncan A Hauser, Jessica M Nelson, Xia Xu, Marek Eliáš, Fay-Wei Li, ***Monodopsis* and *Vischeria* Genomes Shed New Light on the Biology of Eustigmatophyte Algae, *Genome Biology and Evolution***, Volume 13, Issue 11, November 2021, evab233, <https://doi.org/10.1093/gbe/evab233>
 - YourSeq RNAseq Library Kit (FT)
 - Eustigmatophyte algae
- Sota T, Nomura S. **Evolutionary changes in gene expression profiles associated with the coevolution of male and female genital parts among closely related ground beetle species.** Research Square; 2022. DOI: [10.21203/rs.3.rs-1531925/v1](https://doi.org/10.21203/rs.3.rs-1531925/v1)
 - YourSeq RNAseq Library Kit (FT)
 - Carabus (Ohomopterus) ground beetle
- Drobniitch ST, Comas LH, Flynn N, et al. **Drought-Induced Root Pressure in *Sorghum bicolor*.** *Front Plant Sci.* 2021;12:571072. Published 2021 Feb 3. [doi:10.3389/fpls.2021.571072](https://doi.org/10.3389/fpls.2021.571072)
 - YourSeq RNAseq Service
 - Sorghum root

- Chodasiewicz M, Sokolowska EM, Nelson-Dittrich AC, Masiuk A, Beltran JCM, Nelson ADL, Skirycz A. **Identification and Characterization of the Heat-Induced Plastidial Stress Granules Reveal New Insight Into *Arabidopsis* Stress Response**. Front Plant Sci. 2020 Oct 30;11:595792. [doi: 10.3389/fpls.2020.595792](https://doi.org/10.3389/fpls.2020.595792).
 - YourSeq RNAseq library Kit (FT)
 - *Arabidopsis thaliana* chloroplast stress granules (cpSGs)
- Dew-Budd K, Cheung J, Palos K, Forsythe ES, Beilstein MA. **Evolutionary and biochemical analyses reveal conservation of the Brassicaceae telomerase ribonucleoprotein complex**. PLoS One. 2020 Apr 9;15(4):e0222687. [doi: 10.1371/journal.pone.0222687](https://doi.org/10.1371/journal.pone.0222687).
 - YourSeq RNAseq library Kit (FT)
 - *Arabidopsis thaliana* telomerase complex
- Chunmei Li, Nozue K, Maloof JN. **MYCs and PIFs Act Independently in Arabidopsis Growth Regulation**. *G3 Genes/Genomes/Genetics*, Volume 10, Issue 5, 1 May 2020, Pages 1797–1807. <https://doi.org/10.1534/g3.120.401188>
 - YourSeq RNAseq library Kit
 - *Arabidopsis thaliana* seeds
- Shi Y, Hung ST, Rocha G, Lin S, Linares GR, Staats KA, Seah C, Wang Y, Chickering M, Lai J, Sugawara T, Sagare AP, Zlokovic BV, Ichida JK. **Identification and therapeutic rescue of autophagosome and glutamate receptor defects in C9ORF72 and sporadic ALS neurons**. JCI Insight. 2019 Jul 16;5(15):e127736. [doi: 10.1172/jci.insight.127736](https://doi.org/10.1172/jci.insight.127736).
 - YourSeq RNAseq Service (3'DGE)
 - Human C9ORF72 ALS patient (C9-ALS) iMNs
- Babos KN, Galloway KE, Kisler K, Zitting M, Li Y, Shi Y, Quintino B, Chow RH, Zlokovic BV, Ichida JK. **Mitigating Antagonism between Transcription and Proliferation Allows Near-Deterministic Cellular Reprogramming**. Cell Stem Cell. 2019 Oct 3;25(4):486-500.e9. [doi: 10.1016/j.stem.2019.08.005](https://doi.org/10.1016/j.stem.2019.08.005).
 - YourSeq RNAseq Service
 - Mouse *Hb9::GFP+* cells
- Warren JG, Mercado J, Grace D. **Occurrence of Hop Latent Viroid Causing Disease in *Cannabis sativa* in California**. Plant Disease 2019 Aug 21; 103(10): 2699. <https://doi.org/10.1094/PDIS-03-19-0530-PDN>
 - YourSeq RNAseq library Kit (FT)
 - *Cannabis sativa*
- Bellmann J, Monette A, Tripathy V, Sójka A, Abo-Rady M, Janosh A, Bhatnagar R, Bickle M, Mouland AJ, Sternecker J. **Viral Infections Exacerbate FUS-ALS Phenotypes in iPSC-**

Derived Spinal Neurons in a Virus Species-Specific Manner. Front Cell Neurosci. 2019 Oct 22;13:480. [doi: 10.3389/fncel.2019.00480](https://doi.org/10.3389/fncel.2019.00480)

- YourSeq RNAseq Service
- induced pluripotent stem cell (iPSC)-derived spinal neurons (SNs)
- Ichida JK, Staats KA, Davis-Dusenbery BN, Clement K, Galloway KE, Babos KN, Shi Y, Son EY, Kiskinis E, Atwater N, Gu H, Gnirke A, Meissner A, Eggan K. **Comparative genomic analysis of embryonic, lineage-converted and stem cell-derived motor neurons.** Development. 2018 Nov 21;145(22):dev168617. [doi: 10.1242/dev.168617](https://doi.org/10.1242/dev.168617).
 - YourSeq RNAseq library Kit (3'DGE)
 - Embryonic motor neurons