



Join Us for this Live Event

Genomics, transcriptomics, proteomics, and metabolomics are widely used and have greatly contributed to current biological understanding in many areas. However, each omic has challenges and presents only a partial picture of the organism under study. Agilent is a leader in analytical tools, methods, reagents and software for each omic and is at the forefront of developing solutions to integrate multi-omics data. This monthly webinar series will cover topics of interest to researchers using genomics, transcriptomics, proteomics, and metabolomics. Join us to hear how other researchers are applying Agilent's solutions to their research.

ALS genetics: of yeasts and men

Thursday, June 16, 2016

11:00 a.m. EST (New York)
8:00 a.m. PST (San Francisco)
5:00 p.m. CET (Berlin)

Speaker



Julien Couthouis, PhD
Research Associate
Stanford University

[Register now](#)

For Research Use Only. Not for use in diagnostic procedures.

CONNECT WITH AGILENT



[Agilent Home](#) | [About Agilent](#) | [Legal](#) | [Subscriptions](#) | 5301 Stevens Creek Blvd, Santa Clara CA 95051, USA

Agilent is committed to respecting and protecting your privacy. To view the promises that we make to you regarding your personal information, please view [Agilent's privacy statement](#). For questions related to privacy, please send your inquiry to privacy_advocate@agilent.com. If you prefer not to receive promotional e-mails from Agilent in the future, you can [unsubscribe](#) at any time. You can also [update](#) your e-mail address, contact information, or interest profile to ensure you receive the most relevant information.

© Agilent Technologies, Inc. 2016