

# GEN

## Genetic Engineering & Biotechnology News

March 15, 2015  
www.GENengnews.com

TOOLS TECHNOLOGIES TECHNIQUES

## PEOPLE GEN

### Mike McMullen PROFILE

is the new CEO of Agilent. Most recently, he was president and COO at the company. Before that, he was president of Agilent's chemical analysis group. McMullen has 30 years of industry and senior management experience at Agilent. Before joining Agilent's predecessor HP, he worked at DuPont.



### Robin L. Carmichael PROFILE

was promoted to president and COO at Helix BioMedix. Most recently, she was vp of marketing and business development at the company, COO of DERMAdoctor, and vp of marketing at ProCyt. Carmichael has over 20 years of industry experience.



### Kevin Hrusovsky PROFILE

is now CEO of Quantarix. Most recently, he was executive chairman of Quantarix. Before that, Hrusovsky was president of PerkinElmer Life Sciences &



Technology, CEO of Caliper Life Sciences, and CEO of Zymark. He is on the board of directors at Quantarix, 908 Devices, SynapDx, and Solest Energy.

Catalyst Pharmaceuticals named Richard J. Daly and Donald A. Denkhau to its board of directors...3-V Biosciences welcomed Rekha Hemrajani as CFO, svp of business and financial operations...Steven J. Swanson, Ph.D., joined ImmunoCellular Therapeutics as svp of research...Pathway Genomics

hired Anja Kammesheid, Ph.D., as vp of research and development...CannaPharmaRx added Christopher P. Schnitker as CFO...Regen BioPharma hired Todd Caven as CFO...Bob Baltera is now CEO of Laguna Pharmaceuticals...John F. Glenn is the new CFO of Armethen.

Vigilant Biosciences hired Greg Stanley as its vp of global sales and marketing and elected Timothy C. Barabe to its board of directors...GTx appointed Robert J. Wills, Ph.D., to its board of directors...Mitra Biotech named Eric Rowinsky, M.D., to its scientific advisory board...Cosmederm Bioscience added Steve Massah as COO...Perry Calias, Ph.D., joined Pharming as CSO...InVivo added Lorianne Masuoka, M.D., as CMO.

Vaccinogen added Peter Morsing, Ph.D., as head of business development...VistaGen Therapeutics elected Gerard Sanacora Ph.D., M.D., to its clinical and scientific advisory board...Dynavax appointed Laura Brege to its board of directors and chair of the audit

committee...Dance Biopharm welcomed J. Tyler Martin, M.D., as president and member of its board of directors.

DelMar Pharmaceuticals named Lynda Cranston to its board of directors...Aduro hired Sylvia R. Wheeler as svp, corporate affairs and investor relations...Tom Waddell, Ph.D., is the new vp of product development at Xagenic...Lightlake Therapeutics added Arvind Agrawal as evp of medical affairs...Matthew Dallas joined CoLucid Pharmaceuticals as CFO and treasurer...Optivia Biotechnology elected Jean-Michel Scherrmann, to its scientific advisory board.

Exact Sciences appointed Michael S. Wygza to its board of directors...Cancer Genetics hired Randy Goodman, Ph.D., as director of reimbursement and clinical revenue cycle management...Circasia named Lota S. Zoth as nonexecutive director, a member of its remuneration committee, and chair of its audit committee...Peter Radovich is now vp of program leadership and business strategy at Global Blood Therapeutics.

PerkinElmer elected Sylvie Grégoire to its board of directors...Cerenis Therapeutics welcomed Renee Benghozi, M.D., as CMO and Christian Chavy, Michael Davidson, M.D., and Marc Riviere, M.D., to its board of directors...Richard VanBibbe, Ph.D., joined Aortica as vp of clinical affairs...Cidara Therapeutics added Paul Daruwala as CCO and James Balkovec, Ph.D., as svp, research...Clin-

ical Genomics hired Warren Bingham as vp, Asia Pacific...Labcyte welcomed Chris Grimley as vp of marketing...Saad Usmani, M.D., is now acting CMO at SkylineDx.

Sera Prognostics elected Arnold L. Oronsky, Ph.D., to its board of directors...Biofortuna named Ian Johnson chairman to its board of directors...Glyconics Limited has appointed Jeremy Williams to its board of directors...TapImmune hired Patrick D. Yermanian, M.D., as consulting medical director...Ronnie Meninger is now CEO, Hagai Stadler is director on the board of directors, and Sammy Boussiba Ph.D., is chief scientist to the board of directors at Algatech...Symic Biomedical named Barbara Yanni and Bill Newell to its board of directors...Enumeral elected Paul J. Sekhri to its board of directors...Felix Karim, Ph.D., joined Compugen as vp of corporate and business development.

Send your People Announcements to  
people@GENengnews.com

Vaccinogen appointed Hakan Edstrom and Ronald W. Kaiser to its board of directors. Edstrom is CEO of MannKind Corporation, before MannKind, he was president and senior corporate vp at Bausch & Lomb. Edstrom has more than 40 years



of experience. Kaiser is a board member of the NACD. In the past, he was director at Tangoe. He has more than 35 years experience.