

September 18, 2014
(UVPM: 204,575)

Hiring in the Hub 9/12-9/18: MC10, Nara, RAMP & Unified Inbox Name New Executives By Lauren Landry



Boston's business elite has been busy this past week. Catch up on who's going where below:

Alcohol delivery service [Drizly](#) named **Phuc Truong** its senior vice president of strategic ad solutions. He most recently served as the managing director of [Mobext](#), the mobile marketing agency for [Havas Media](#), where he led mobile and advertising efforts in North America. Truong will report to Drizly Co-founder and CEO Nick Rellas, who shared in a statement:

[Truong's] extensive experience in cultivating strategic partnerships and developing best-in-class ad solutions will help take Drizly to the next level.

[Nara](#) named **Jana Eggers** president of the recommendation startup. Eggers will be responsible for leading Nara's daily operations across all divisions, and comes to the company from nonprofit software supplier [Blackbaud](#), where she served as senior vice president of products and marketing. At Blackbaud, Eggers managed an 850-person team.

Scott Pomerantz was appointed the new chief executive officer and president for flexible electronics maker [MC10](#). He most recently spent seven years working for [Broadcom](#), a producer of semiconductors for wired and wireless communications, where he served as the senior vice president and general manager of wireless connectivity.

[CoachUp](#) announced New England Patriots Wide Receiver **Julian Edelman** [will be joining the private coaching service](#) as a member of its Athlete Advisory Council. Hear more from Edelman below:

Haverhill, Mass.-based [Primary Design](#) promoted **John Schroeder** from his role as long-time director of account services to that of the firm's executive vice president. In his new position, he will be tasked with helping Primary Design expand into new markets, as well as solidifying the company as a one-stop shop for clients "looking for everything from brand design to environment display design and installation to digital application and website creation."



Singapore-headquartered conversation and social collaboration platform [Unified Inbox](#) announced it has acquired Concord, Mass.-based [SocialGrow](#). As part of the acquisition, SocialGrow's Co-founder and Chief Marketing Officer **Ken Herron** will join Unified Inbox as the company's head of marketing. (More on that [here](#).)

Mark Roskey joined Lexington, Mass.-based high definition diagnostics company [Quanterix](#) as its vice president and general manager of applications and reagents. Roskey previously worked in a similar VP/GM role at [PerkinElmer](#), and boasts more than 25 years of experience in the life science industry.

[RAMP](#) added two new executives to its team. The end-to-end, cloud-based video content platform named **Bob Orlando** and **Erik Herz** as its chief financial officer and vice president of business development, respectively. Shared RAMP CEO Tom Wilde in a statement:

We are thrilled to welcome Bob and Erik to the RAMP team. Bob brings extensive experience building high tech companies and will help us grow our market presence as we add to our roster of marquee customers for our media and enterprise video solutions. Erik has been a highly-respected innovator in video technology solutions for years and will enable us to add to our growing list of enterprise video solutions and engage with businesses who want to use video to communicate with customers, prospects, and employees.

Biopharmaceutical company [Akebia Therapeutics](#) recently [announced](#) the expansion of its management team, including three key additions: **Brad Maroni**, senior vice president and chief medical officer; **Mark De Rosch**, vice president of regulatory affairs; and **Tamara Dillon**, vice president of human resources. The Cambridge company is still looking to hire more senior talent, however, to bolster its growth. Interested? Go [here](#).

Streetwise Media announced a hire of its own this week. Kyle Alspach will be leaving *The Boston Globe* to join Streetwise as our new national tech editor. (More on that [here](#).)

Bonus opportunity: [The New England Venture Capital Association](#) is looking to bring on a full-time staff member to manage its new student outreach program. The program is meant to ensure the state's top talent stays in Boston post-graduation. Click [here](#) to learn more.