



Dr. Nick Seay
COO, Cellular Dynamics International (USA)

Biography

Nick Seay formerly practiced law in the area of biotechnology, intellectual property and related business transactions. He currently holds the title of Chief Technology Officer, managing the research activities, at Cellular Dynamics International Inc. (CDI), a privately held stem cell company founded in 2005 by the scientific team of Dr. James Thomson, Dr. Craig January and Dr. Tim Kamp in partnership with Tactics II Ventures, LLP. Thomson, January, and Kamp are internationally renowned for their research accomplishments related to stem cells and to human heart function. Cellular Dynamics is making a suite of cellular products from pluripotent cells for drug testing, beginning with tests for cardiac risk and toxicity. CDI also intends to be a leader in newly developing technology of induced pluripotent stem cells, which are stem cells made from samples of a living person's skin.