

Media Release

InSphero Named #1 Swiss Startup for 2014

Schlieren, Switzerland, September 25, 2014 – Leading supplier of 3D microtissues for in vitro toxicity and efficacy testing earns top honors in annual Swiss startup competition.

InSphero AG, the leading provider of assay-ready 3D microtissues and 3D-focused contract research services, was recognized as the #1 company in the [Top 100 Swiss Startups Rankings](#) for 2014, per results published in Switzerland on September 18th. The honor marks the fourth consecutive year that InSphero has made the Top 10 in the rankings, but is the first time the company has been awarded the #1 overall startup designation across all industries. Founded in 2009 in Zurich, Switzerland, InSphero now has commercial offices in Germany and the United States, and counts the world's top 15 pharmaceutical and cosmetics companies amongst its customers. Already in 2014, the company has added novel [3D pancreatic microislets](#) to its 3D InSight™ microtissue line, along with a [spheroid imager](#), and 3D-optimized [cytotoxicity assay kit](#) incorporating technology from Promega.

Dr. Jan Lichtenberg, Chief Executive Officer and co-founder of InSphero says the award is affirmation that the company's effort to expand the reach of their customer base along with the breadth of their product portfolio are paying off. "We started with the mission to make 3D cell culture a robust and easy to use tool for more predictive drug testing. Today, all major pharma companies have InSphero products in-house. We are proud that the panel has recognized this success."

InSphero, a spin-off of the ETH Zurich and the University of Zurich, placed fifth overall in the competition in 2013. The Swiss Startups Rankings are determined by a panel of 100 Swiss startup experts, each of whom individually names and ranks ten promising startup companies. Ten points are awarded to the top nominee on each list, down to one point for the tenth, and the startups with the highest combined points enter the Top 100 rankings.

For more information about InSphero, visit www.insphero.com.

InSphero contact

Dr. Randy Strube, Director of Global Marketing, phone +1 800-779-7558 ext. 102,
randy.strube@insphero.com, www.insphero.com

Dr. Jan Lichtenberg, CEO and Co-Founder, phone +41 44 5150490,
jan.lichtenberg@insphero.com, www.insphero.com

About InSphero

InSphero is a leading supplier of organotypic, biological in vitro 3D microtissues for highly predictive drug testing. The company, headquartered in Zurich, Switzerland, with subsidiaries in the USA and in Germany, currently counts all of the top ten global pharmaceutical and cosmetics companies as customers. InSphero 3D InSight™ Microtissues enable more biologically relevant in vitro applications in efficacy and toxicology. The spin-off company of the Swiss Federal Institute



InSphero AG
Wagistrasse 27
CH-8952 Schlieren
Switzerland
Tel: +41 (44) 515049-0
Fax: +41 (44) 515049-1
www.insphero.com

of Technology (ETH) Zurich and the University Zurich has been recognized for its scientific and commercial achievements with a number of national and international awards, including being named the #1 Swiss Startup for 2014. InSphero is certified to the ISO 9001:2008 standard for its Quality Management System.

Follow us on   and www.insphero.com