



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

**Media Release – For Immediate Distribution**

## **InSphero Inc. Streamlines 3D Microtissue Business Operations in US**

***US subsidiary of Swiss-based InSphero AG initiates direct product shipments, progresses towards full-scale US bioproduction capabilities.***

**Schlieren, Switzerland – February 1, 2016** – [InSphero AG](http://www.insphero.com), the leading supplier of easy-to-use solutions for production, culture, and assessment of organotypic 3D cell culture models, has begun accepting orders and processing all financial transactions with customers in North America through its US subsidiary InSphero Inc. The February 1 transition immediately streamlines access to InSphero products and services for US-based customers, and positions the company to initiate operations in its recently completed US bioproduction laboratory in Brunswick, Maine.

Stewart Hunt, General Manager of InSphero Inc., states the transition, “Will dramatically simplify order fulfillment for customers in our largest market by allowing them to order directly from a US source. This eliminates customs processing, additional handling fees, as well as potential delays encountered by those currently using shipment brokers to manage direct international purchases.”

The transition accompanies another milestone for InSphero Inc., which has assembled an expert team of 11 US-based employees over the past year. Says Hunt, “We’ve moved quickly, adding key scientific staff to lead the laboratory build-out, technology transfer, and initiate product manufacturing in our state-of-the-art Brunswick facility. Concurrently we have established an experienced team of sales managers, field application scientists, marketing, and product management professionals to provide responsive education and support to our North American customers.”

For more information about InSphero, visit [www.insphero.com](http://www.insphero.com).

###

### **InSphero contacts**

Dr. Randy Strube  
Director of Global Marketing  
Phone +1 800-779-7558 ext. 102  
[randy.strube@insphero.com](mailto:randy.strube@insphero.com)

Stewart Hunt  
General Manager, InSphero Inc.  
Phone +1 800-779-7558 ext. 101  
[stewart.hunt@insphero.com](mailto:stewart.hunt@insphero.com)



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

## About InSphero

InSphero provides superior biological relevance to *in vitro* testing with its easy-to-use solutions for production, culture and assessment of more organotypic 3D cell culture models. The company's patented 3D cell culture platforms and methods enable large-scale, reproducible production of a broad range of assay-ready 3D InSight™ Microtissues derived from liver, pancreas, tumor, heart, brain and skin. These models and contract research services utilizing them help to identify promising drugs and toxic liabilities with greater predictivity at early development stages, enabling better pre-clinical decision making, saving development cost, and shortening time to market. InSphero technologies drive significant findings in [peer-reviewed journals](#), through collaborative projects such as [EU Body on a Chip](#) and [HeCaToS](#), and have gained validation in the world's largest government institutions and pharmaceutical, chemical and cosmetics companies. This 3D know-how is also being applied in the diagnostics field to aid development of personalized chemotherapeutic strategies for the treatment of cancer.

Founded in 2009, the privately held company is headquartered in Schlieren, Switzerland with subsidiaries in the United States (Brunswick, ME) and Waldshut, Germany. It has been recognized for its scientific and commercial achievements with a number of national and international [awards](#).

Follow us on [Twitter](#) and [www.insphero.com](http://www.insphero.com).