

**Media Release – For Immediate Distribution****InSphero and Hamilton Announce Collaboration to Improve the Production and Handling of Spheroids Using Advanced Liquid Handling Technology**

*Both partners have a proven track record for being a successful entrepreneur and a technical disruptor at the intersection of 3D in vitro technologies and automation*

**Schlieren, Switzerland – May 19, 2021** InSphero AG, the pioneer of 3D cell-based assay technology and [Hamilton](#), an innovation-driven leading supplier of automated liquid handling equipment, today announced the companies have entered into a collaboration to improve the production and handling of 3D microtissues (spheroids).

“InSphero develops microtissues using the most advanced 3D cell culture technologies to produce highly predictive, organotypic *in vitro* models that are ready to use in critical applications including drug discovery, efficacy and safety,” said Wolfgang Moritz, PhD, InSphero co-founder. “Through this collaboration, we will leverage Hamilton’s breakthrough MagPip™ liquid and cell handling technology to optimize the processes used to produce and manipulate these microtissues.”

With the ability to rapidly and accurately transfer liquid volumes in a sub-microliter range, the MagPip™ liquid handling system allows for precise and gentle handling of even the most delicate microtissues without harming their integrity.

“The combination of InSphero’s advanced 3D in vitro models and our MagPip™ technology on the Microlab VANTAGE™ will allow us to explore new ways to not only produce microtissues, but potentially expand how they are used in a wide range of applications,” said Patrick Pfenninger, Senior Product Manager, Hamilton. “We expect MagPip enable new opportunities in microtissue sorting and storage, as well as in the areas of downstream processing for drug efficacy and safety testing.”

View this [video](#) to see how the MagPip™ system safely and gently handles the InSphero microtissues.

For more information about InSphero 3D microtissues, visit [www.insphero.com/products](http://www.insphero.com/products).

###

**About InSphero**

InSphero is the pioneer of industrial-grade, 3D-cell-based assay solutions and scaffold-free 3D organ-on-a-chip technology. Through partnerships, InSphero supports pharmaceutical and biotechnology researchers in successful decision-making by accurately rebuilding the human physiology *in vitro*. Its robust and precisely engineered suite of 3D InSight™ human tissue platforms are used by major pharmaceutical companies worldwide to increase efficiency in drug discovery and safety testing. The company specializes in liver toxicology, metabolic diseases (e.g., T1 & T2 diabetes and NAFLD & NASH liver disease), and oncology (with a focus on immuno-oncology and PDX models). The scalable Akura™ technology underlying the company’s 3D InSight™ Discovery and Safety Platforms includes 96 and 384-well plate formats and the Akura™ Flow organ-on-a-chip system to drive efficient innovation throughout all phases of drug development.

Learn more at [www.insphero.com](http://www.insphero.com) and follow us on [Twitter](#) and [LinkedIn](#).

### **InSphero Contact**

Dr. Frank Junker

Chief Business Officer

Phone +41 44 5150490

[frank.junker@insphero.com](mailto:frank.junker@insphero.com)

### **Images**



### **About Hamilton**

Hamilton Bonaduz AG is a leading global manufacturer that provides automated workstations for liquid handling and laboratory automation technology to the scientific community. With a focus on innovative design, Hamilton products integrate patented liquid handling technologies into a portfolio that includes liquid handling platforms, standard application-based solutions, small appliances, consumables, and OEM liquid handling solutions. Hamilton products are known for their advances in life sciences, clinical diagnostics, forensics and biotechnology, offering reliability, performance and flexibility. Hamilton ensures an ongoing commitment to quality and uses state-of-the-art production facilities in Reno, Nevada and Bonaduz, Switzerland, and has achieved a worldwide ISO 9001 certification. Hamilton is privately owned and headquartered in Reno, Nevada; Franklin, Massachusetts; and Bonaduz, Switzerland, along with subsidiaries around the world.  
[www.hamiltoncompany.com/robotics](http://www.hamiltoncompany.com/robotics)

### **Hamilton Contact**

Dr. Nikolai Kraut

Product Manager Cellomics

Phone: +41 58 610 13 90

[nkraut@hamilton.ch](mailto:nkraut@hamilton.ch)

### **Images**

