



Our mission is to create a healthier society by enabling life science companies to invent novel compounds in a safer, more efficient, faster, and animal-free way. As the pioneer in engineering and industry-scale production of assay-ready 3D microtissues, InSphero provides solutions that help pharmaceutical, cosmetics, and chemical companies accelerate safety and efficacy testing, and bring better drugs to market.

InSphero engineered the Akura™ Flow microphysiological discovery platform to be compatible with our complete portfolio of 3D InSight™ microtissues including disease models for NAFLD/NASH and diabetes, **3D tumor models engineered for Immuno-Oncology applications** and liver models for predictive liver toxicology. It opens the door to a near unlimited variety of pre-clinical efficacy and toxicity testing applications in a multi-tissue, microfluidic assay format.

InSphero is headquartered in Schlieren, Switzerland, a biotech hub of Zurich. We are looking for a highly motivated individual to join our team as

#### **Postdoctoral Researcher in Immuno-oncology Organ-on-chip Solutions (2-year contract)**

In the frame of a **partnering project between F. Hoffmann-La Roche** in Basel and InSphero, you will contribute to advancing immune-competent organ-on-chip solutions to best support immuno-oncology drug development. Based on our scalable and flexible Akura™ technology and 3D InSight™ microtissue platforms, you will be developing industry-relevant applications for personalized drug development strategies in the oncology field. You will be working in a highly interdisciplinary group at the interface of bioengineering, 3D cell culture and advanced analytical methods and gain unique insights in industry-relevant assay development.

Starting date: as soon as possible. Place of work: Schlieren, reporting to the Head of Technology and Platforms Group.

#### **Your Profile**

- PhD in Cell biology, Immunology or Oncology with deep knowledge in cell/tissue-based assays (especially in immuno-oncology) and basic know-how in advanced in-vitro assays (incl. microfluidics)
- Minimum 2 years post-PhD experience
- Solid scientific track record (incl. publications)
- Laboratory experience in primary cell culture techniques, ideally 3D tissue culture, immune cell assays
- Microscopy skills (especially confocal microscopy) and image processing
- Basic knowledge in scripting (data analysis) and statistics
- Strong communication and presentation skills, networker, team player
- Professional English is a must, additional languages are a plus
- Applicants must be either Swiss citizens, EU-nationals or have the legal right to live and work in Switzerland

We are looking forward receiving your application (resume, cover letter, diplomas, job references) by e-mail at [hr@insphero.com](mailto:hr@insphero.com).



**InSphero AG**  
Wagistrasse 27a  
CH-8952 Schlieren  
Switzerland  
[www.insphero.com](http://www.insphero.com)