



InSphero is an innovative biotech company in a rapidly growing sector helping its customers in pharmaceutical and chemical industry to make their compound discovery and development safer and more efficient. InSphero has partnered with leading pharmaceutical and biotech companies to develop and produce highly functional, organotypic, three-dimensional cell-based microtissues in the areas of toxicology, liver diseases, diabetes and oncology. With manufacturing sites in both Switzerland and Brunswick, ME, InSphero is well positioned to be a global solution provider in drug discovery and development. We are looking for a highly motivated individual to join our team in the Brunswick Bio-production unit. Starting date: as soon as possible.

### **3D Cell Culture Lab Technician (100%)**

#### **Main responsibilities**

- Routine production of 3D Insight™ microtissues and cell culture media for commercial sales and internal use following established SOPs
- Thorough documentation of production processes in manufacturing checklists
- Isolation of primary cells
- Support of product development and maintenance (validation of new cell lots, testing of media batches, evaluation of new cell sources)
- General laboratory maintenance (lab cleaning, equipment monitoring, preparation of reagent aliquots, lab inventory)

#### **Profile**

- BSc/MSc in Biology or other Life Sciences with experience in cell culture technologies and animal handling/simple surgery
- Ability to work under aseptic conditions
- Good hand and eye coordination to perform biochemical assays and use technical equipment with accuracy
- Experience in working according to standard operating procedures (ISO 9001 certification)
- Good skills with Microsoft Office Applications
- Flexible with regards to working hours
- Resilient and reliable team player but also able to work independently with high attention to detail
- Authorization to work in the U.S. for any employer (candidates must be U.S. citizens or Permanent Resident Aliens)

We are looking for a highly reliable personality who assumes responsibility. We are offering an international and dynamic work environment, which will allow you to participate in shaping InSphero's future. We are looking forward to receiving your application by e-mail at [hr@insphero.com](mailto:hr@insphero.com).



**InSphero Inc.**  
TechPlace  
74 Orion Street  
Brunswick, ME 04011  
USA  
[www.insphero.com](http://www.insphero.com)