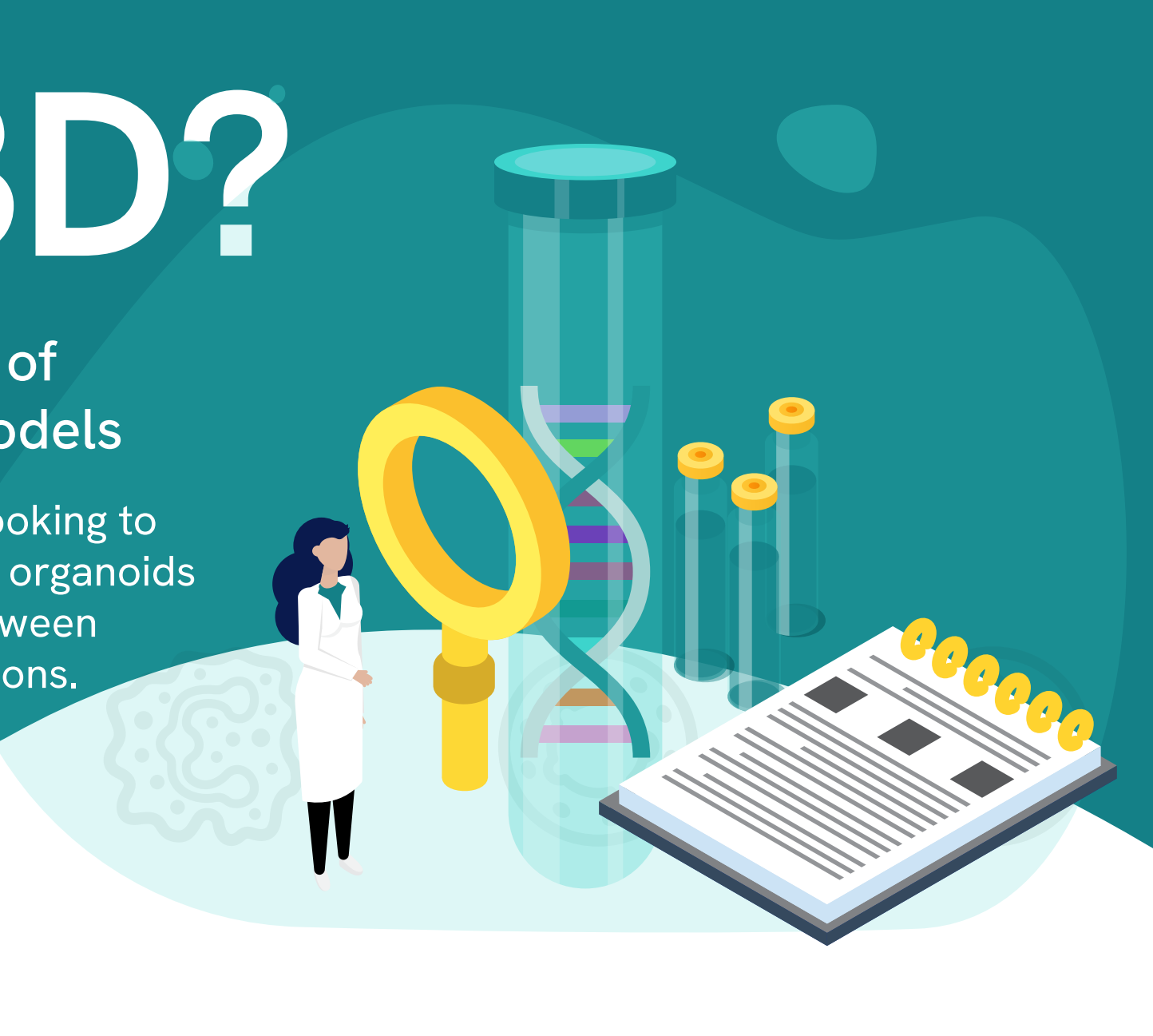



# Why 3D?

## Stripping Complexity Out of Multi-Dimensional Cell Models


More than ever, researchers are looking to 3D cell cultures, microtissues, and organoids to bridge the translational gap between 2D cell cultures and *in vivo* conditions.





**Advantages**


- More physiologically relevant
- Closely mimic cell-cell and cell-extracellular matrix interactions
- Help for a better understanding of biology and thus to identify new targets
- Improved lead validation and characterization for effective and safe drugs




**Landscape**

The need for more biologically relevant disease models has spurred a steady **8% annual growth in research** on 3D cell models since 2013\*.

\* Source: Google Scholar



Publications (Total)



**Applications**

3D cell models can improve and accelerate different steps along the drug discovery workflow.

Model disease *in vitro* to better understand the underlying biology

Identify new targets using more physiologically relevant models

Screen drug candidates in high-throughput format




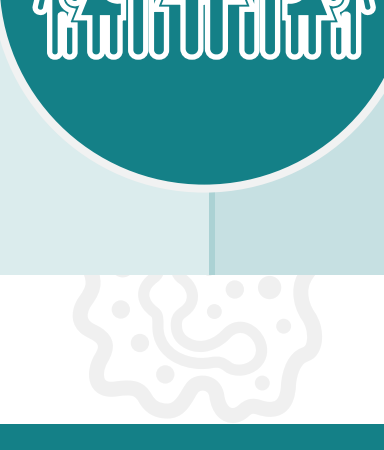
Perform efficacy and toxicity profiling *in vitro* for lead optimization & selection

Preclinical use of 3D models for

- Regenerative Medicine
- Transplantation
- Precision medicine

## A precise distinction

The difference between 2D and 3D

2D Cell Model		3D Cell Model
<p><b>Lacks <i>in vivo</i> imitation</b></p> <p>Doesn't mimic <i>in vivo</i> - like structures or molecular mechanisms</p>		<p><b><i>In vivo</i> imitation</b></p> <p><i>In vivo</i> - like architecture and stable expression of relevant genes and proteins</p>
<p><b>Limited cell interactions</b></p> <p>Limited cell-cell interactions or cell-extracellular environment interactions</p>		<p><b>Cell interactions</b></p> <p>Interactions of cell-cell and cell-extracellular microenvironment</p>
<p><b>Simple culture</b></p> <p>High performance and reproducibility, short culturing times</p>		<p><b>Complex culture</b></p> <p>Challenging to define proper seeding methods and conditions, lower reproducibility, long culture formation times</p>
<p><b>Changed &amp; uniform characteristics</b></p> <p>Changed morphology, loss of diverse phenotype, and polarity</p>		<p><b>Preserved &amp; diverse characteristics</b></p> <p>Preserved morphology, diverse phenotypes, and polarity</p>

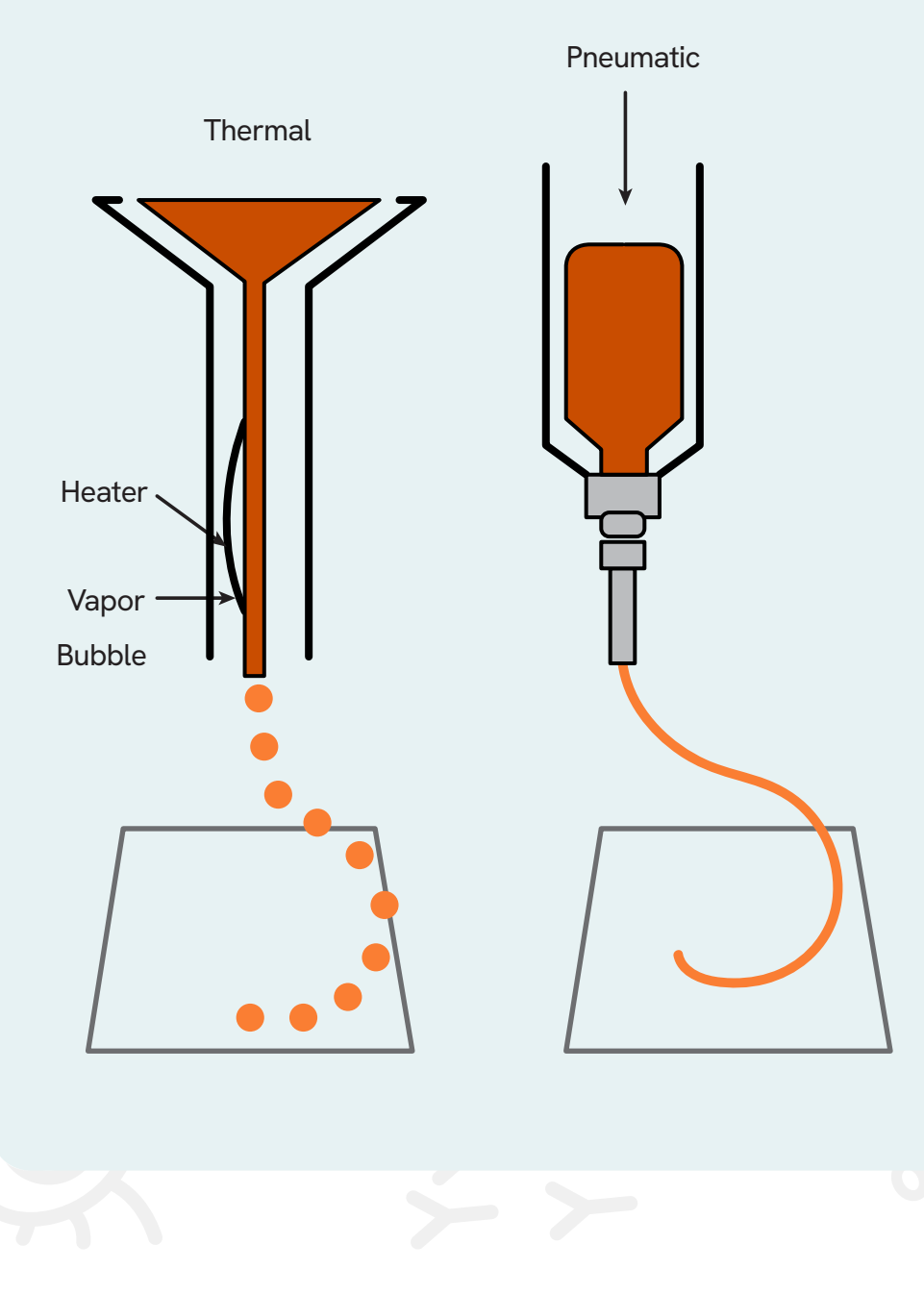
## Types of cell models

Structure	Spheroid	Organoids	Organ-on-a-chip
Cell Type	<ul style="list-style-type: none"> <li>• Cell line monoculture or co-culture</li> </ul>	<ul style="list-style-type: none"> <li>• Multiple organ-specific cell types (ESC, iPSC or patient derived)</li> </ul>	<ul style="list-style-type: none"> <li>• Multiple organ-specific cell types (primary cells, iPSCs)</li> </ul>
Assembly method	<ul style="list-style-type: none"> <li>• Cell aggregation</li> </ul>	<ul style="list-style-type: none"> <li>• Self-assembly</li> <li>• Differentiation</li> </ul>	<ul style="list-style-type: none"> <li>• Self-assembly</li> <li>• Differentiation</li> </ul>
Advantages	<ul style="list-style-type: none"> <li>• Easy-to-use protocol</li> <li>• High reproducibility</li> <li>• Co-culture ability</li> </ul>	<ul style="list-style-type: none"> <li>• Multiple cell lineages</li> <li>• <i>In vivo</i> - like complexity and architecture</li> </ul>	<ul style="list-style-type: none"> <li>• <i>In vivo</i> - like architecture, microenvironment, and flow</li> <li>• Transport of soluble factors, nutrients, and oxygen</li> <li>• Good imaging compatibility</li> </ul>
Disadvantages	<ul style="list-style-type: none"> <li>• Simplified architecture</li> <li>• Lack vasculature</li> <li>• Only partial tissue components</li> <li>• Limited imaging depth</li> </ul>	<ul style="list-style-type: none"> <li>• High variability</li> <li>• Potential lack of vasculature</li> <li>• Limited imaging depth</li> </ul>	<ul style="list-style-type: none"> <li>• High variability</li> <li>• Lower throughput</li> <li>• Careful handling is required, e.g. air bubbles</li> </ul>

## Common seeding methods

Different seeding methods, each with unique characteristics, must be selected based on the assay type and research questions at hand.

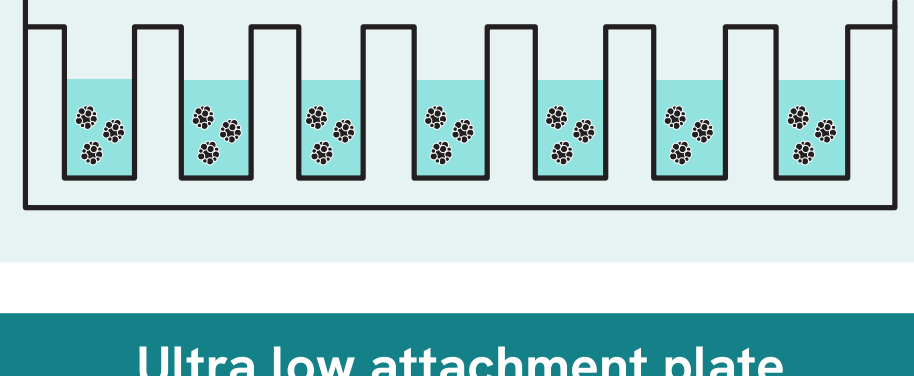
**Bioprinter**



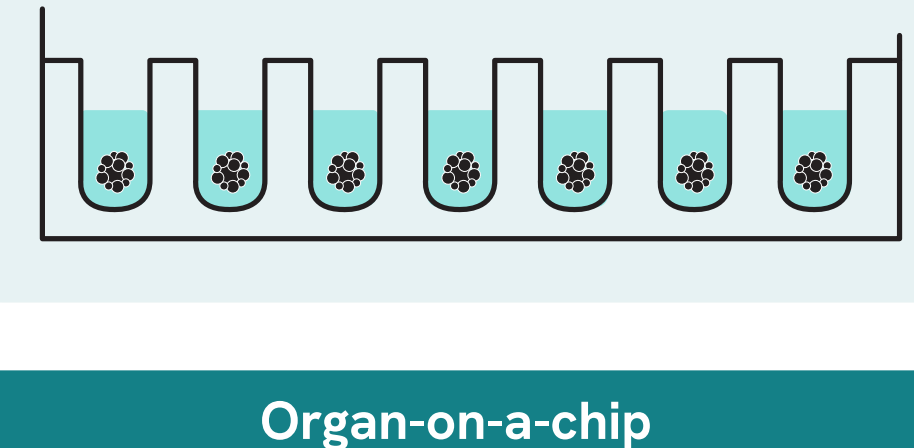
Thermal: Heater, Vapor Bubble

Pneumatic

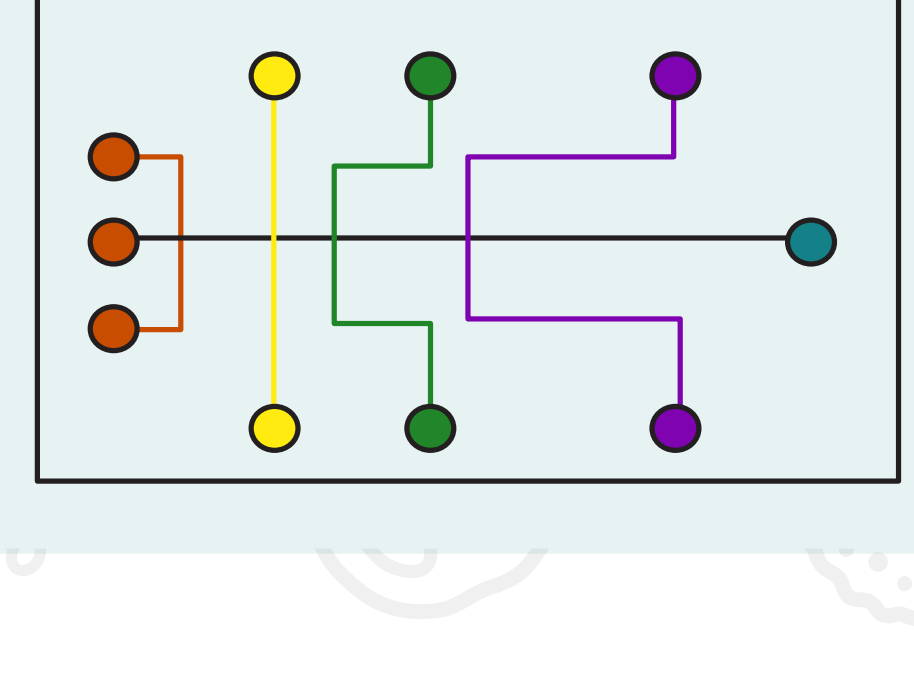
**Biological hydrogels**



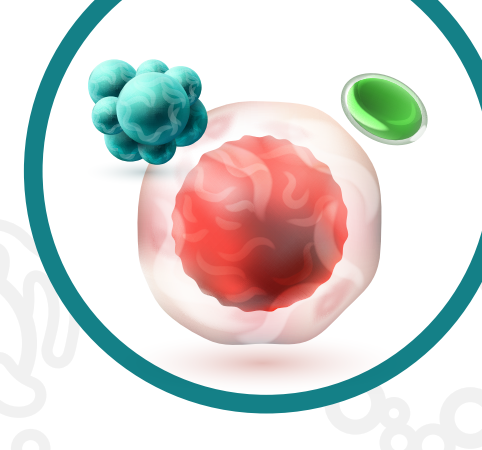
**Ultra low attachment plate**




**Organ-on-a-chip**



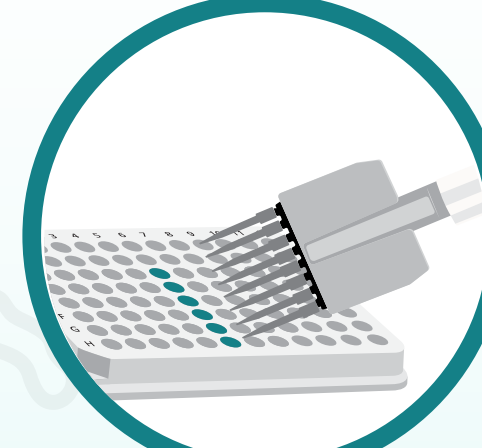
## The 3D cell culture workflow




**Selection of cell model and growth method**



**Grow 3D cell model**




**Treatment**




**Sample preparation**

- Staining and Clearing
- Preparing Detection Assay
- Nucleic Acid Isolation & NGS Library Prep



**Read out**

- High Content Imaging
- Multimode Plate Readers
- NGS



**Analysis**

## Providing confidence and ease in 3D cell models

What to look for in technologies to help make it a seamless workflow.



Reliability



Reproducibility



Speed



Biologically relevant results



Analyze data in context

For more information, visit us at [www.revivity.com](http://www.revivity.com)

