



Welcome to the October edition of our newsletter!

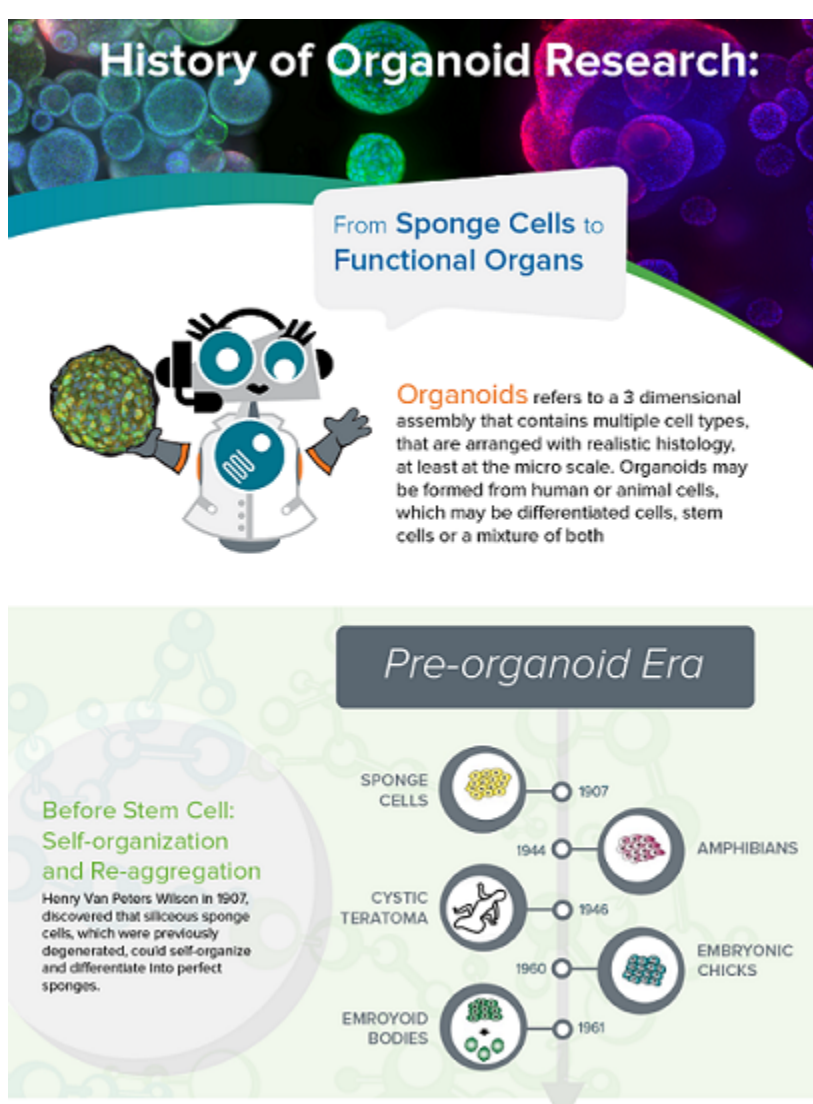
Bringing you our latest discoveries in high-throughput screening, genomic and cellular analysis, colony selection, and microplate detection. Visit us at: moleculardevices.com

NEW
History of Organoid Research

From Sponge Cells to Functional Organs

Interestingly, the idea of organoids is not new. Today's organoid technology is the product of decades of research. In fact, the foundations of the concept go back to the beginning of the 20th century.

Here is a brief rundown of the history of organoids and how Molecular Devices found itself a spot on the timeline with its extensive organoid research solutions.



[Learn More](#)

Organoids for disease modeling and *in vitro* drug screening

Leverage automation and advanced high-content imaging to increase throughput and data from your organoid assays

Organoids are an essential tool for disease modeling and assessment of compound effects. Automated imaging and analysis of organoids are equally as important for quantitative assessment of organoid phenotypic changes and to increase experiment throughput. Read our new application note to learn about an integrated system that allows automated monitoring, maintenance, and characterization of the growth and differentiation of organoids and stem cells, and a workflow to test the effects of various compounds.



[Download the Research](#)



Phenotypic screening with iPSC-derived cardiomyocytes and neurons

Identify therapeutic effects and adverse responses to compounds earlier in the drug discovery process

In this eBook, we use both imaging and calcium oscillation analysis to develop profiles of compounds in iPSC-derived cardiomyocytes such as HERG blockers, β -adrenergic agonists, and environmental toxins. iPSC-derived neuronal cultures were evaluated with neuromodulators as well as environmental toxins.

[Download eBook](#)

Assess virus neutralization with a rapid, HTS-friendly assay

Discover how you can identify SARS-CoV-2 neutralizing antibodies in one hour!

The worldwide COVID-19 pandemic caused by the SARS-CoV-2 virus has necessitated the fast-tracked development of many research tools for understanding this virus's pathogenesis, as well as vaccine discovery and development.

As neutralizing antibodies are key biomarkers of immune response and vaccine efficacy, levels of neutralizing antibodies in patient serum samples are an important parameter to be able to monitor efficiently.



[Download Application Note](#)



Let's connect, live and in person!

We are looking forward to connecting at SynBioBeta, now January 25-27th 2022.

Molecular Devices is a proud sponsor of [SynBioBeta](#). After this landmark year for synthetic biology applications, we cannot wait to see what our community has been creating.

Can't wait for 2022?
Check out our virtual session from SynBioBeta 2020.

Join our team and highlighted QPix® Microbial Colony Picker users, as they explore diverse applications across gene editing, strain engineering, and protein engineering.

[Watch the Session](#)

EVENTS

SBI[®] (Society of Biomolecular Imaging and Informatics) 2021
October 5-6, 2021 | Virtual

TEDD Annual Meeting 2021
October 14, 2021 | Wädenswil, Switzerland

ILMAC
October 19-21, 2021 | Basel, Switzerland

SLAS 2021 Building Biology in 3D Symposium - Hybrid
October 26-27, 2021 | Jupiter, FL

Discovery UK
October 26-27, 2021 | London, UK

Society for Neuroscience (SfN) 50th Annual Meeting 2021
November 8-11, 2021 | Virtual
November 13-16, 2021 | Chicago, IL

Cell Bio Virtual 2021
December 1-10, 2021 | Virtual

Antibody Engineering & Therapeutics 2021
December 12-16, 2021 | San Diego, CA

JOIN OUR TEAM

Ready to reach your full potential? Click [here](#) to see what life is like at Molecular Devices.