

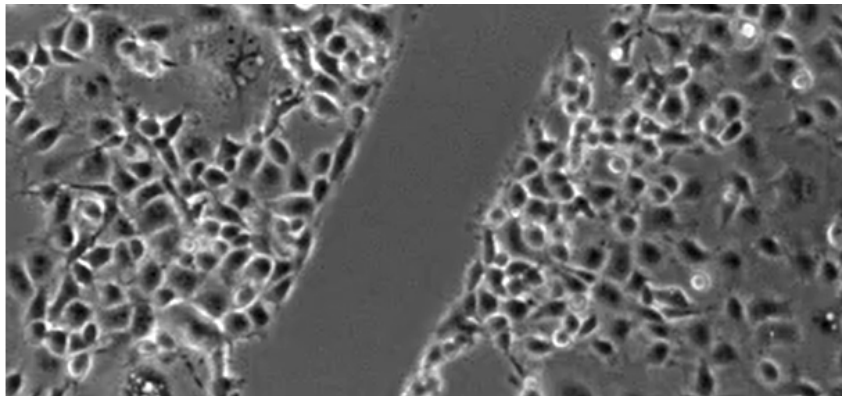


Stencell

Q&A

Q/ What is the smallest design you can craft out of the foil?

The smallest feature that has been tested was 250 μm .



COS-7 cells colonizing a gap induced by the removal of a PDMS stencil made of 2 rectangular chambers of 10 mm * 3 mm separated by 250 μm .

Image credit: © Pierre-Olivier Strale

Q/ Are Stencells sterile?

No, however Stencells can be treated with UV under the hood and/or washed with ethanol.

Q/ How long can I store the Stencells?

Stencell can be stored in the protective foil for up to 2 years. Do not remove the protective foil. If you wish to store the assembled Stencells, avoid dust deposition.

Q/ Can I design my own Stencell?

Yes. We have put up a process to be able to answer any design requirement very fast. [Let us](#)

[know about your project](#), we'll be pleased to have this discussion with you!

Q/ Can I stack two (or more) Stencells on top of each other?

Sure, it has been performed. See [Pasturel et al. Adv Biosystems 2020](#)

Q/ What supports are compatible with Stencells compatible with plastic Petri dishes as well?

Stencells have been successfully used with:

- Microscope slide
- Plastic Petri dishes
- Transwell
- Polyacrylamide gel

Q/ Are Stencells reusable?

No, Stencells are single-use consumables.

February 24, 2022