

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com info@eaglebio.com



EAGLE
BIOSCIENCES



Stampwell Q&A

Q/ What should I use the rectangular stamp for?

This shape is perfect for the imaging of your zebrafish or medaka embryos.

Q/ What should I use the V-shape 300μ stamp for?

This shape is perfect for the imaging of your medium-sized rounded 3D objects like medium-sized spheroids and organoids. If you need to work on bigger objects, you might use the V-shape 500μ. If you need to work on smaller objects, [contact us](#) and we will study an optimized option for you.

Q/ What should I use the V-shape 500μ stamp for?

This shape is perfect for the imaging of your large size rounded 3D objects like big spheroids and organoids. If you need to work on smaller objects, you might use the V-shape 300μ. If you want to work on bigger objects, [just tell us](#) and we will study an optimized option for you.

Q/ What should I use the U-shape stamp for?

This shape is perfect to easily grow spheroids or organoids in a reproducible, fast, and cheap way. If you need to work on more objects, or if you want that your stamp is compatible with other culture supports, [just tell us](#) and we will study an optimized option for you.

Q/ Which microscopes are compatible with Stampwell?

You can use any kind of microscope (inverted, upright, with/without deeping lens, confocal, spinning disk, two-photon excited, stereomicroscope, macroconfocal, videomicroscope, ...). The only parameters that matter are the objective field of view and its numerical aperture / working distance.

Q/ How to make sure that my samples fit the wells?

Please contact us, we will be happy to discuss your specific needs and provide you advice on the best shapes.

Q/ What hydrogels are compatible with Stampwell ?

- 1-2% agarose
- Collagen-1
- Gelrite
- PDMS

Q/ My cells can't grow without Matrigel. How to do so ?

We recommend to prepare, on ice with coolen reagents and consumables, 1 mL of a mix containing:

- 5 - 20% (v/v) of Matrigel
- Culture medium
- Cells

Q/ What kind of tools do I need to use Stampwell?

All you need is:

- a Stampwell
- hydrogel (agarose, alginate, phytagel, methylcellulose, PDMS ...),
- 35 mm petri dish, with or without glass bottom,
- a pipetman
- a cutter blade
- a non-wounding spatula
- and a microscope :-)

May 24, 2022

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com info@eaglebio.com



EAGLE
BIOSCIENCES