



# Gri3D® Hydrogel Microcavity Plate

## - Imaging

## Technical Specifications



SPEC015a, October 2025

## Gri3D® Hydrogel Microplate Specifications

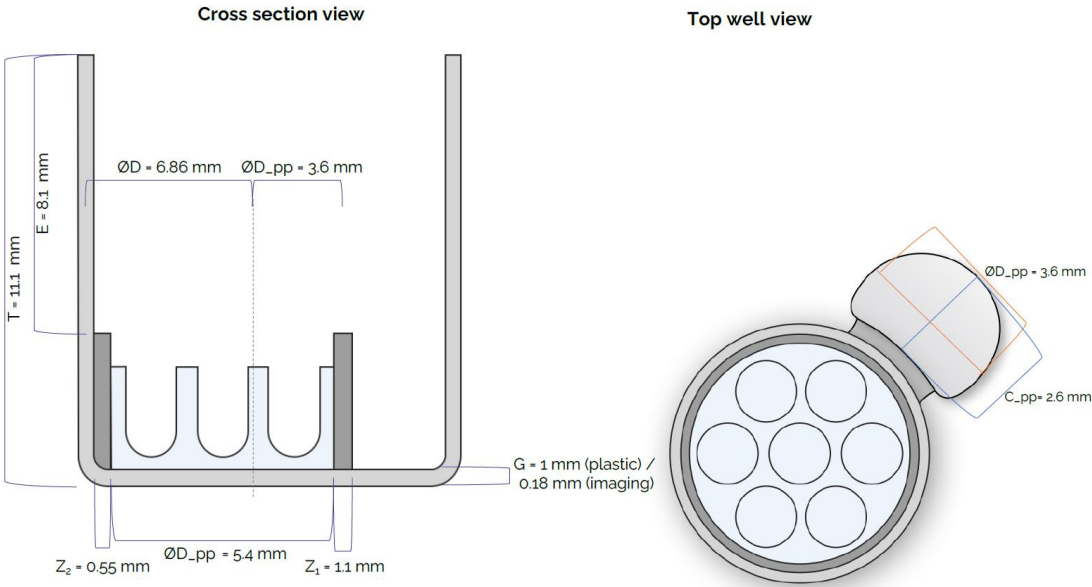
---

<b>Article number:</b>	GRI3D-96IBI-S-96-400/500/600/800
<b>Product name:</b>	Gri3D® Hydrogel Microcavity Plate - Imaging
<b>Pack size:</b>	1x Gri3D® Microplate with lid, sterile, 180 µm imaging bottom
<b>Plate &amp; lid material:</b>	COP (Cyclo-olefin polymer), transparent
<b>Hydrogel microcavities:</b>	Cell-repellent PEG-based hydrogels promote uniform organoid aggregation and culture within precisely engineered high-aspect ratio microcavities for several weeks up to several months.

### Plate dimensions:

Parameter:	Description:	Measurements:
L	Length (with skirt)	127.89 mm
l	Length (without skirt)	123.19 mm
1	Left well position A1	14.44 mm
X1_pp	Left pipetting port position A1	18.14 mm
Y1	Upper well position A1	11.30 mm
Y1_pp	Upper pipetting port position A1	7.60 mm
P	Horizontal well to well distance	9.00 mm
Q	Vertical well to well distance	9.00 mm
Q_pp	Vertical pipetting port to pipetting port distance	9.00 mm
B	Width (with skirt)	85.60 mm
b	Width (without skirt)	80.99 mm
Xn	Right well position	113.44 mm
Yn	Lower well position	74.30 mm
S	Plate without lid	14.45 mm
R	Skirt	3.35 mm
G	Thickness bottom	180 µm
∅D	Diameter top	6.86 mm
∅d	Diameter ground	5.40 mm
∅D_pp	Diameter pipetting port	3.6 mm
C_pp	Cross section pipetting port	2.6 mm
T	Depth	11.10 mm
B	Bottom from plate ground	3.35 mm

C	Distance bottom side well to plate support	3.17 mm
E	Depth to seeding ring	8.10 mm
Z1	Width wall seeding ring to pipetting port	1.10 mm
Z2	Width wall seeding ring	0.55 mm



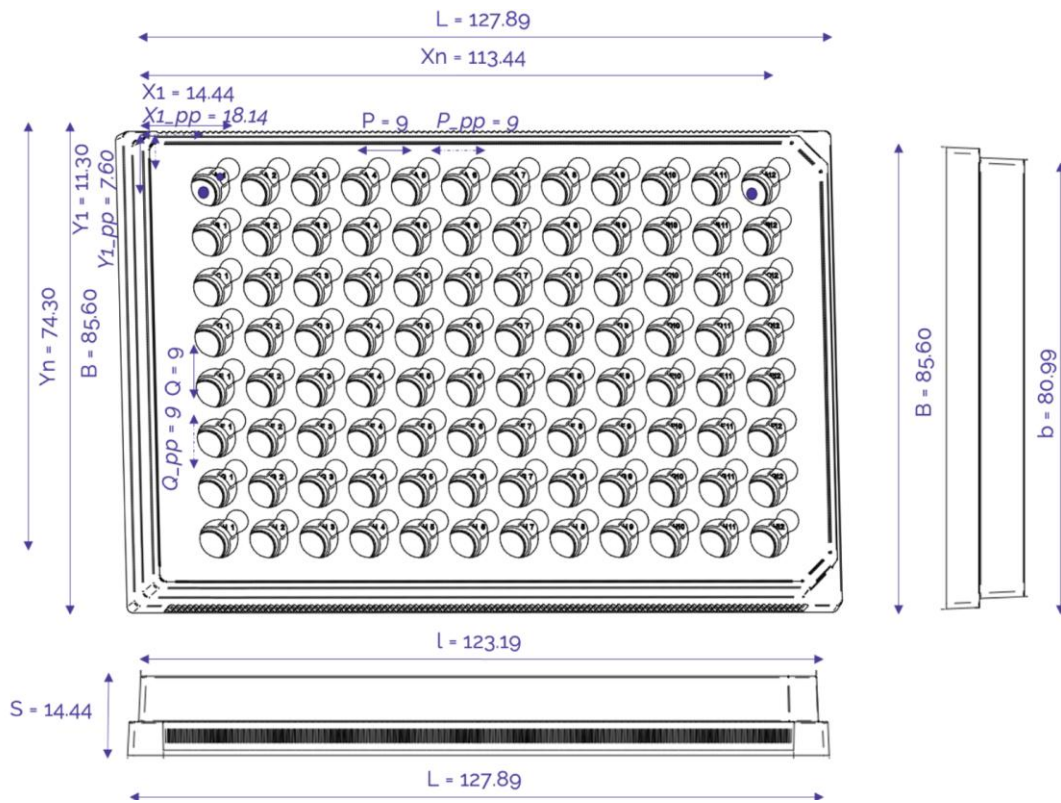
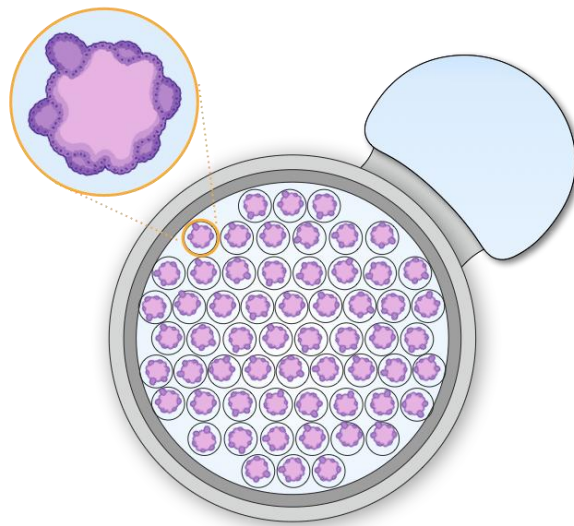


Figure 1: Technical specifications of the Gri3D® Hydrogel Microplate in mm.

**A**



**B**



*Figure 2: **A.** Angled view of Gri3D® Hydrogel Microcavity Plate. **B.** Image of one well containing microcavities, and pipetting port.*



InSphero AG  
Schlieren, Switzerland

InSphero Inc.  
Brunswick, ME, USA

If you have more questions, please refer to  
the FAQs section on [shop.insphero.com](https://shop.insphero.com)

Sign up for latest news and updates at  
[insphero.com/newsletter](https://insphero.com/newsletter)

Follow us on LinkedIn and X

InSphero is ISO 9001:2015 certified

All rights reserved, © 2025 InSphero AG.  
3D InSight, Akura and InFloat are  
trademarks of InSphero AG. For life  
science research only. Not for use in  
diagnostic procedures.