



Gri3D[®] Hydrogel Microcavity Plate - Plastic **Technical Specifications**



SPEC015b, October 2025

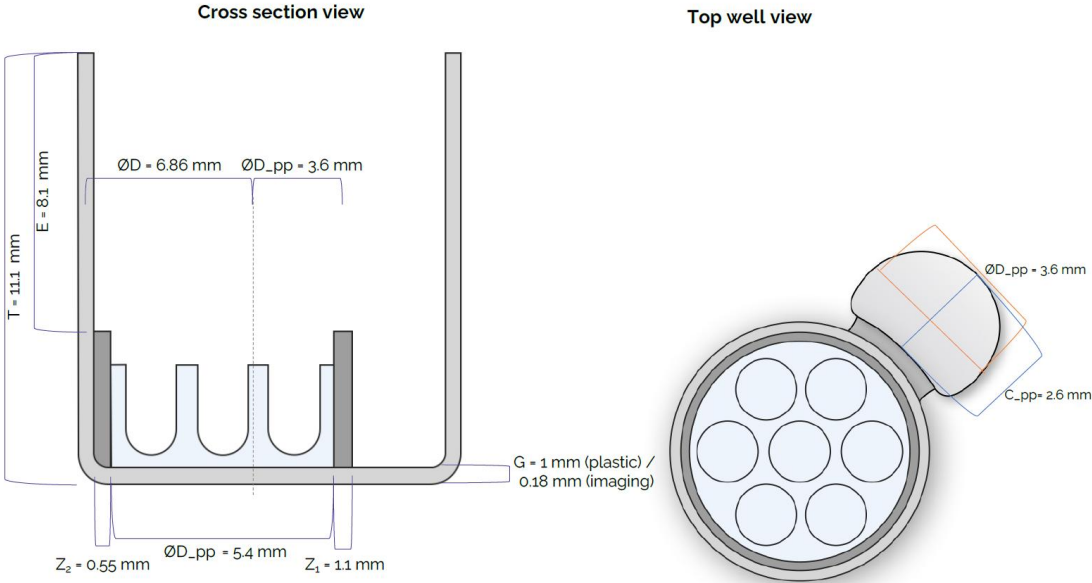
Gri3D® Hydrogel Microplate Specifications

Article number:	GRI3D-96P-S-96-400/500/600/800
Product name:	Gri3D® Hydrogel Microcavity Plate - Plastic
Pack size:	1x Gri3D® Microplate with lid, sterile, 1 mm plastic bottom
Plate & lid material:	Polystyrene, transparent
Hydrogel microcavities:	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	1 mm
∅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	2.35 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



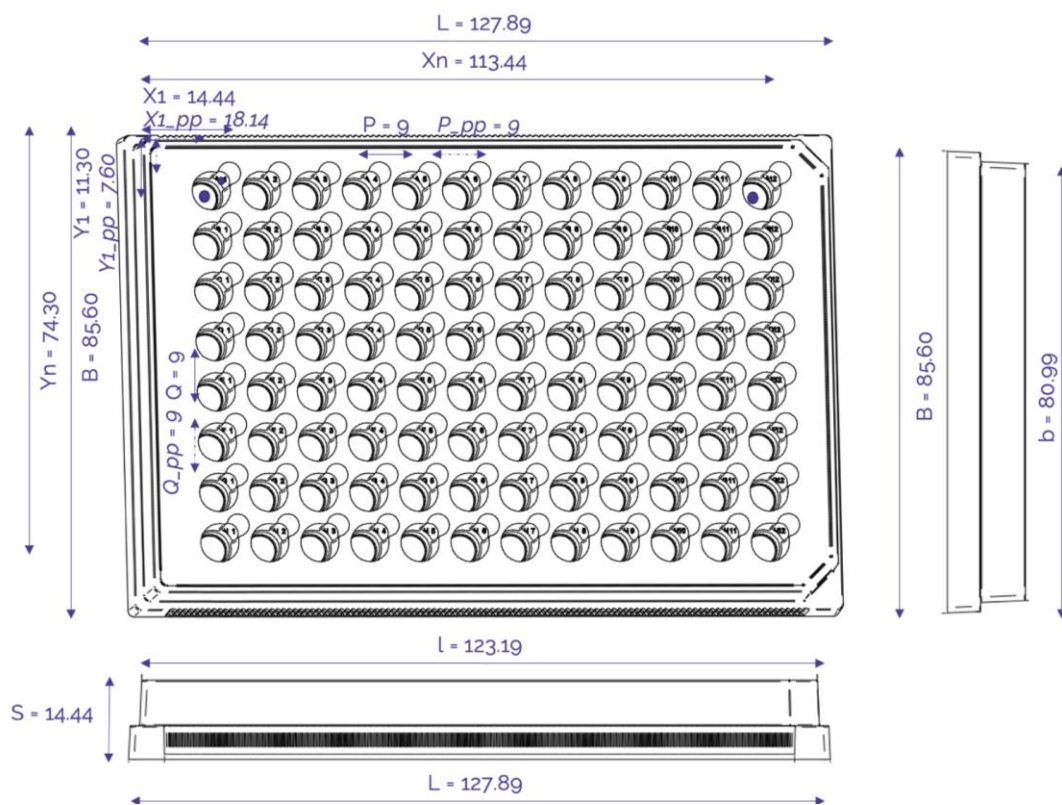
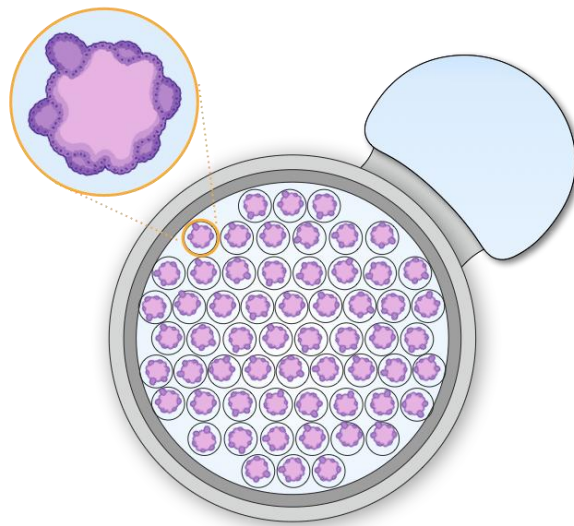


Figure 1: Technical specifications of the Gri3D® Hydrogel Microplate in mm.

A



B



*Figure 2: **A.** Angled view of Gri3D® Hydrogel Microplate. **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

Sign up for latest news and updates at
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.