

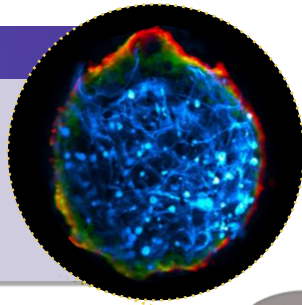
BLOOD-BRAIN BARRIER

Cells: human primary astrocytes, pericytes, and endothelial cells - 1:1:1 ratio

3000 cells per microcavity



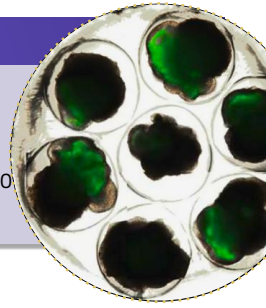
Time of culture: 3-5 days
Gri3D format: Gri3D® 96WP 600 µm
[Simonneau et al., 2021](#)



RETINA

Cells: mouse embryonic stem cells,
300 cells per microcavity

Time of culture: 26 days
Gri3D format: Gri3D® 96WP 1600
[Decembrini et al., 2020](#)



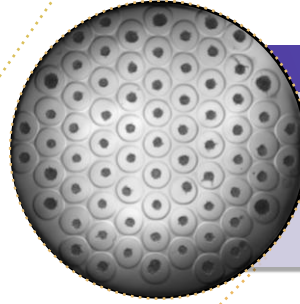
MELANOMA

Cells: mouse primary melanocytes (seeded fresh from biopsy)

500 cells per microcavity

Time of culture: 1-2 weeks

Gri3D format: Gri3D® 96WP 500 µm



LIVER

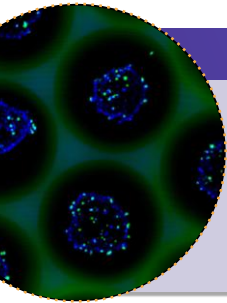
Cells: iPSC-derived hepatocytes, endothelial cells, and stellate cells - 10:3:1 ratio

4200 cells per microcavity

Time of culture: up to 1 month

Gri3D format: Gri3D® 96WP 500 µm

[Meseguer-Ripolles et al., 2021](#)



PANCREAS

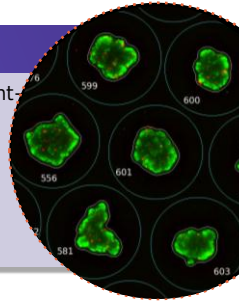
Cells: human pancreatic adenocarcinoma patient-derived organoids

200 cells per microcavity

Time of culture: 1 week

Gri3D format: Gri3D® 96WP 500 µm

[Roch et al., ISSCR 2022 - Poster](#)



INTESTINE

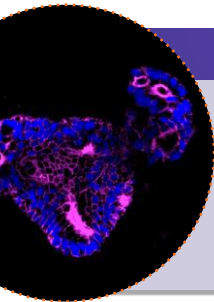
Cells: human adult stem cell-derived intestinal organoids

200 cells per microcavity

Time of culture: 7 days

Gri3D format: Gri3D® 96WP 500 µm

[Brandenberg et al., 2020](#)



(IMMUNO)-ONCOLOGY

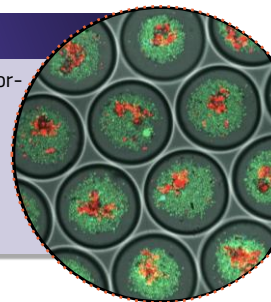
Cells: human colorectal cancer derived, and tumor-infiltrating lymphocytes at different ratios

200 cells per microcavity

Time of culture: 1 week

Gri3D format: Gri3D® 96WP 500 µm

[Dutta et al., 2021](#)



GASTRULOID

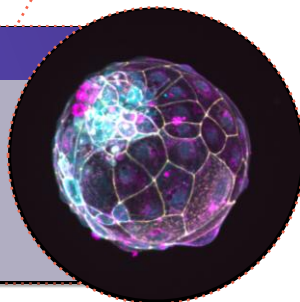
Cells: human pluripotent stem cells

45 cells per microcavity

Time of culture: 3-4 days

Gri3D format: Gri3D® 96WP 200 µm

[Khoei et al., 2023](#)



BONE MARROW

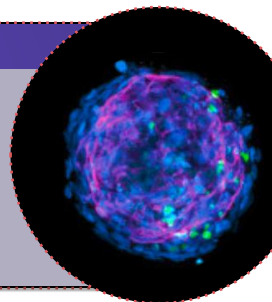
Cells: HUVECs* + mesenchymal stem cells - 75:25 ratio, addition of HSPC** at different ratios

500 cells per microcavity

Time of culture: 1 week

Gri3D format: Gri3D® 96WP 400 µm

[Giger et al., 2022](#)



*Human umbilical vein endothelial cells / ** Hematopoietic stem and progenitor cells