



Organoid media formulation #2

Components required

Item	Manufacturer	Catalog #	Storage
Organoid Growth Kit 1C	ATCC	ACS-7102	-20°C or below
L-Glutamine	ATCC	30-2214	-20°C or below
DMSO	ATCC	4-X	2-8°C
Advanced DMEM:F12	Thermo Fisher Scientific	12634028	2-8°C
HEPES	Thermo Fisher Scientific	15630080	2-8°C
B-27 Supplement	Thermo Fisher Scientific	17504-044	-20°C or below
HA-R-Spondin1-Fc 293T (RSPO1) Conditioned Media	For each 250 mL of complete media, 25 mL of RSPO1 conditioned media is required. Refer to vendors instructions to prepare conditioned medium from Trevigen Cultrex® HA-R-Spondin1-Fc 293T Cells (Trevigen Cat # 3710-001-01). The protocol for cell culture and conditioned medium generation is available at: https://trevigen.com/docs/protocol/protocol_3710-001-01.pdf		
Refer to manufacturer documentation for expiration dates and safe handling information.			

Complete 1X growth medium preparation procedure (makes ~250 mL)

1. Thaw B-27 and L-Glutamine on ice or in a refrigerator at 2-8°C. Aliquot stock bottles into working volumes and store at -20°C or below. Avoid multiple freeze/thaw cycles. Thaw DMSO at ambient temperature. Place Organoid Growth Kit at ambient temperature.
2. Prepare supplemented basal medium. Aseptically combine the following components in a sterile 250 mL bottle.

Item	Volume
Advanced DMEM:F12	215.0 mL
HEPES	2.5 mL
L-Glutamine	2.5 mL
B-27	5.0 mL
Total volume	225.0 mL

3. Briefly centrifuge the vials in the Organoid Growth Kit to ensure the material is at the bottom of the vial.
4. Aseptically reconstitute the individual kit components in the indicated buffer. After adding buffer to each vial, incubate for 15 minutes at room temperature. Mix by repeated pipetting. If the N-Acetyl Cysteine is difficult to dissolve, periodic vortexing and incubation in a 37°C water bath for 10-20 minutes can help the material enter solution.



Item	Catalog #	Buffer	Volume of buffer
Noggin	ACS-7201	Supplemented basal medium	1.0 mL
Nicotinamide	ACS-7214	Supplemented basal medium	2.5 mL
N-Acetyl-Cysteine	ACS-7215	Supplemented basal medium	1.0 mL
FGF-10	ACS-7205	Supplemented basal medium	1.0 mL
FGF-7	ACS-7207	Supplemented basal medium	1.0 mL
A 83-01	ACS-7210	DMSO	0.1 mL
SB 202190	ACS-7212	DMSO	0.1 mL

Note: Once reconstituted components should be used immediately. Do not store reconstituted components.

5. Aseptically combine the reconstituted kit components, conditioned media and supplemented basal media.

Item	Volume
RSPO1 Conditioned Media	25.0 mL
Supplemented basal media	218.5 mL
Noggin	1.0 mL
Nicotinamide	2.5 mL
N-Acetyl-Cysteine	1.0 mL
FGF-10	1.0 mL
FGF-7	1.0 mL
A 83-01	0.1 mL
SB 202190	0.1 mL
Total volume	~250.0 mL

6. Aseptically filter the complete growth medium through an 0.22 μ M PES bottle-top filter unit.
7. (Optional) Place the supplied sticker on the final collection bottle to indicate media preparation is complete. Label with an expiration date 4 weeks from date of preparation.

Notes

- Once prepared, store complete medium at 2-8°C in the dark.
- Complete medium expires after 4 weeks or at the expiration date of any of the components, whichever comes first.
- Do not freeze complete medium and avoid extended light exposure.

Warranty and Disclaimer

ATCC® products are warranted for 30 days from the date of shipment, and is valid only if the product is stored and used according to the information included on this product information sheet. This product is intended for laboratory research purposes only. It is not intended for use in humans. While ATCC uses reasonable efforts to include accurate and up-to-date information on this product sheet, ATCC makes no warranties or representations as to its accuracy.



ATCC does not warrant that such information has been confirmed to be accurate. ATCC is not liable for any damages or injuries arising from receipt and/or use of this product.