

Culture of Human Liver Organoids

Using Cultrex™ UltiMatrix Basement Membrane Extract

This protocol provides a procedure for subculturing normal human liver organoids. This protocol was modified from the submerged method described in Huch, M. *et al.* (2015) *Cell*. **160**: 299. The protocol provided below is intended to culture organoids from normal human liver tissues using Cultrex™ UltiMatrix Reduced Growth Factor (RGF) Basement Membrane Extract as a scaffold. The majority of reagents used in this protocol were sourced from the Bio-Techne brands of R&D Systems™ and Tocris Bioscience™.

TABLE // 01

Materials Needed for Human Liver Organoid Culture

Product Name	Supplier	Catalog #
Cultrex Organoid Harvesting Solution	R&D Systems	3700-100-01
Cultrex UltiMatrix Reduced Growth Factor Basement Membrane Extract	R&D Systems	BME001-05
Advanced DMEM/F-12 Cell Culture Medium	Thermo Fisher	12634010
GlutaMAX™	Thermo Fisher	35050061
HEPES	Tocris Bioscience	3173
Penicillin/Streptomycin	Various	Various
N21-MAX Supplement	R&D Systems	AR008
N-2 MAX Supplement	R&D Systems	AR009
N-Acetylcysteine	Tocris Bioscience	7874
Gastrin I (Human)	Tocris Bioscience	3006
Nicotinamide	Tocris Bioscience	4106
Y-27632 dihydrochloride (Rho Kinase inhibitor)	Tocris Bioscience	1254
A 83-01 (ALK5 inhibitor)	Tocris Bioscience	2939

Product Name	Supplier	Catalog #
Forskolin	Tocris Bioscience	1099
DAPT	Tocris Bioscience	2634
Dexamethasone	Tocris Bioscience	1126
Recombinant Human EGF	R&D Systems	236-EG
Recombinant Human R-Spondin 1	R&D Systems	4645-RS
Recombinant Human Noggin	R&D Systems	6057-NG
Recombinant Human FGF-10	R&D Systems	345-FG
Recombinant Human FGF-19	R&D Systems	969-FG
Recombinant Human BMP-7	R&D Systems	354-BP
Recombinant Human HGF	R&D Systems	294-HG
Recombinant Human Wnt-3a	R&D Systems	5036-WN

Other Required Reagents

- Distilled (DW) or deionized water (DI)
- Phosphate buffered saline (PBS)

Reagent Preparation

Use aseptic technique at all times during this protocol. This protocol is optimized for human liver organoids. Organoids from other tissues may have different culture requirements.

1. Thaw Cultrex UltiMatrix RGF Basement Membrane Extract (BME) on ice for four hours or overnight at 2 - 8 °C (on ice in the refrigerator).
2. Prepare Liver Organoid Initiation Medium (Table 2), Liver Organoid Expansion Medium (Table 3), and Liver Organoid Differentiation Medium (Table 4).
3. Sterile filter the media.

TABLE // 02

Human Liver Organoid Initiation Medium

Reagent Name	[FINAL]
Advanced DMEM/F-12 Cell Culture Medium	NA
N21-MAX Supplement	1X
GlutaMAX™	2 mM
HEPES	10 mM
Penicillin/Streptomycin	1X
N-2 MAX Supplement	1X
Nicotinamide	10 mM
A 83-01	5 µM
N-Acetylcysteine	1.25 mM
Recombinant Human FGF-10	100 ng/mL
Recombinant Human R-Spondin 1	0.5 µg/mL
Gastrin I Human	10 nM
Recombinant Human BMP-7	25 ng/mL
Recombinant Human EGF	50 ng/mL
Recombinant Human HGF	25 ng/mL
Recombinant Human Noggin	25 ng/mL
Recombinant Human Wnt-3a	100 ng/mL
Forskolin	10 µM
Y-27632 dihydrochloride	10 µM

TABLE // 03

Human Liver Organoid Expansion Medium

Reagent Name	[FINAL]
Advanced DMEM/F-12 Cell Culture Medium	NA
N21-MAX Supplement	1X
GlutaMAX™	2 mM
HEPES	10 mM
Penicillin/Streptomycin	1X
N-2 MAX Supplement	1X
Nicotinamide	10 mM
A 83-01	5 µM
N-Acetylcysteine	1.25 mM
Recombinant Human FGF-10	100 ng/mL
Recombinant Human R-Spondin 1	0.5 µg/mL
Gastrin I Human	10 nM
Recombinant Human BMP-7	25 ng/mL
Recombinant Human EGF	50 ng/mL
Recombinant Human HGF	25 ng/mL
Forskolin	10 µM

TABLE // 04

Human Liver Organoid Differentiation Medium

Reagent Name	[FINAL]
Advanced DMEM/F-12 Cell Culture Medium	NA
N21-MAX Supplement	1X
GlutaMAX™	2 mM
HEPES	10 mM
Penicillin/Streptomycin	1X
N-2 MAX Supplement	1X
A 83-01	0.5 µM
Recombinant Human FGF-19	100 ng/mL
Gastrin I Human	10 nM
Recombinant Human BMP-7	25 ng/mL
Recombinant Human EGF	50 ng/mL
Recombinant Human HGF	25 ng/mL
DAPT	10 µM
Dexamethasone	30 µM

Methods for Expanding and Differentiating Human Liver Organoids

Expansion

1. Prepare a suspension of isolated and dissociated human liver tissues as detailed in Huch, M. *et al.* (2015) *Cell* **160**: 299.
2. Resuspend the cell pellet in Cultrex UltiMatrix RGF Basement Membrane Extract and aliquot into wells by dispensing 50 μL of the Cultrex UltiMatrix RGF Basement Membrane Extract liver cell mixture in the center of each well of a 24-well plate to create domes.

Note: The Cultrex UltiMatrix RGF BME domes should not touch the sides of the well.
3. Incubate the plate in the cell culture incubator for 15 minutes to polymerize the Cultrex UltiMatrix RGF Basement Membrane Extract.
4. Add the appropriate volume ($\sim 500 \mu\text{L}/\text{well}$) of Liver Organoid Initiation Medium (Table 2).
5. Return the plate containing the organoid cultures to the cell culture incubator to promote growth.
6. After 3 days, aspirate the culture medium from each well and add fresh Liver Organoid Expansion Medium (Table 3) every other day ($\sim 500 \mu\text{L}/\text{well}$).

Note: Medium should be gently aspirated from and pipetted into the corner of the well away from the Cultrex UltiMatrix RGF BME/ organoids to prevent their disruption.
7. Passage organoids (See Passage Liver Organoids section) or start differentiation after 7-10 days.

Differentiation

1. Following 7-10 days of expansion, aspirate the Liver Organoid Expansion Medium, wash once with PBS, then add an equal volume ($\sim 500 \mu\text{L}/\text{well}$) of Liver Organoid Differentiation Medium (Table 4). Incubate for an additional 11-13 days.
2. The culture medium should be aspirated from each well and replaced with fresh Liver Organoid Differentiation Medium every other day.

Note: Medium should be gently aspirated from and pipetted into the corner of the well away from the Cultrex UltiMatrix RGF BME/ organoids to prevent their disruption.

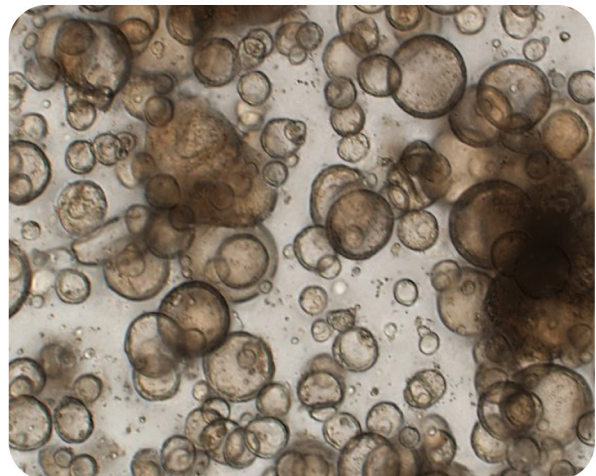
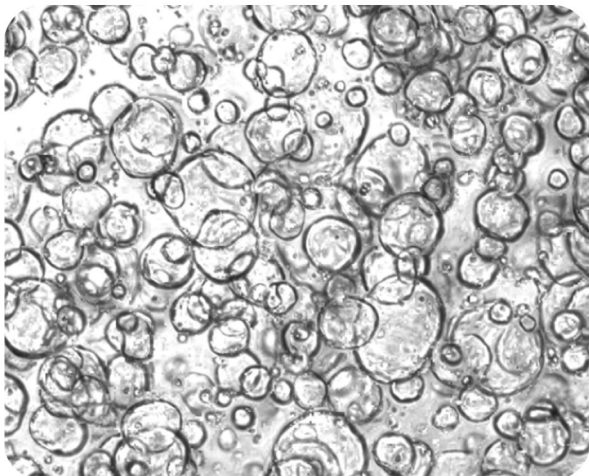


Figure 1. Undifferentiated and Differentiated Human Liver Organoids. Representative brightfield images of undifferentiated (left) and differentiated (right) human liver organoids that were cultured using [Cultrex UltiMatrix RGF Basement Membrane Extract](#) (R&D Systems, Catalog # BME001-05) and the reagents listed in this protocol.

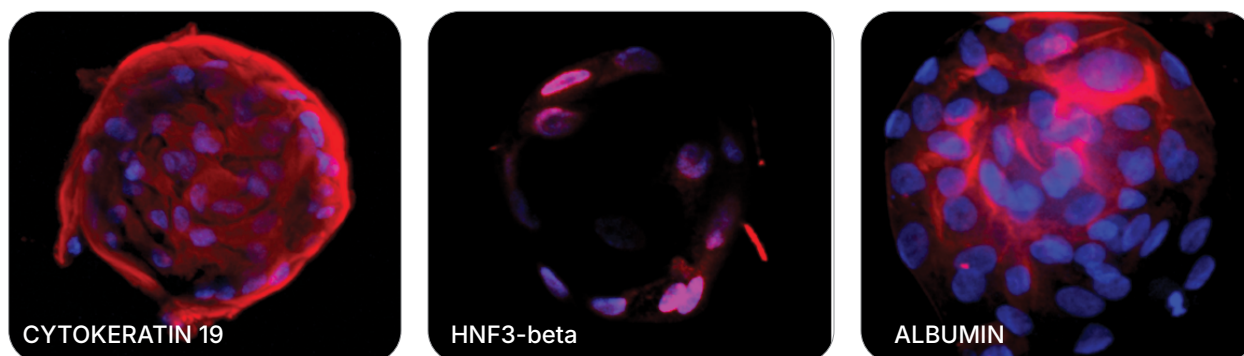


Figure 2. Characterization of Differentiated Human Liver Organoids. Adult stem cell-derived liver organoids were cultured using [Cultrex UltiMatrix RGF Basement Membrane Extract](#) (R&D Systems, Catalog # BME001-05) and the other reagents listed in this protocol. Differentiated human liver organoids were stained for proteins characteristic of differentiated hepatocytes using a [Sheep Anti-Human Cytokeratin 19 Polyclonal Antibody](#) (R&D Systems, Catalog # AF3506), a [Goat Anti-Human HNF3-beta Polyclonal Antibody](#) (R&D Systems, Catalog # AF2400), and a [Mouse Anti-Human Serum Albumin Monoclonal Antibody](#) (R&D Systems, Catalog # MAB1455). Organoids were counterstained with [DAPI](#) (Tocris, Catalog # 5748).

Passaging Liver Organoids

1. View liver organoids under the microscope. Each well should contain approximately 500 organoids for optimal growth. Organoid cultures exhibiting rapid growth may be split 1:4 during passaging, while slow growing cultures may benefit from a 1:1 split. Make this determination prior to harvesting to estimate reagent needs prior to starting.

Note: Organoid density is important for optimal growth; too many organoids will strain culture resources, while too few organoids lack paracrine signaling necessary to sustain growth.
2. Transfer the 24-well plate containing liver organoids from the cell culture incubator to the cell culture hood.
3. Aspirate the medium without disturbing the Cultrex UltiMatrix RGF BME-contained organoids at the bottom of the well.
4. Gently wash each well with 10 volumes of cold (2-8 °C) PBS (Table 5). Be careful not to disrupt the basement membrane matrix-containing organoids.
5. Aspirate the PBS, and add 10 volumes of cold (2-8 °C) Cultrex Organoid Harvesting Solution to each well (Table 5).
6. Incubate the plate at 2-8 °C or on ice for 30–90 minutes with moderate shaking. This incubation

is complete when the basement membrane matrix dome is no longer visible at the bottom of the well and the organoids are seen floating at the bottom of the well.

Note: Dislodging the dome with a cell scraper or pipet may accelerate this process.

7. Once the matrix depolymerizes, transfer the contents of the well into a tube on ice. Single wells may be transferred to a microtube while multiple domes may necessitate a 15 mL or 50 mL conical tube.
8. Centrifuge the tube at 500 x g for 5 minutes at 2-8 °C in a swinging bucket rotor to pellet the organoids. Aspirate the supernatant.
9. Wash the organoids with 10 volumes of cold (2-8 °C) PBS, and repeat centrifugation at 500 x g for 5 minutes at 2- 8 °C in a swinging bucket rotor to pellet the organoids. Aspirate the PBS. Add fresh ice-cold Liver Organoid Expansion Medium.
10. Pipet up and down three times with a serological pipette to mix the organoids.

TABLE // 05

Suggested Working Volumes of PBS and Cultrex Organoid Harvesting Solution

Plate Type	Volume of Basement Membrane Matrix	Volume of PBS and Organoid Harvesting Solution
96-well plate	5 µL	50 µL
48-well plate	25 µL	250 µL
24-well plate	50 µL	500 µL

- Centrifuge the tube at 500 × g at room temperature for 3 minutes.
- Aspirate the medium, but be careful not to disturb the organoid pellet.
- Resuspend the organoids in Cultrex UltiMatrix RGF Basement Membrane Extract, and dispense 50 µL of the mixture in the center of each well of a 24-well plate to form domes. Follow the density/splitting ratios recommended in Passaging Protocol Step 1.

Note: The Cultrex UltiMatrix RGF BME-contained organoids should not touch the sides of the well.

- Incubate the plate in the cell culture incubator for 15 minutes to polymerize the Cultrex UltiMatrix RGF Basement Membrane Extract.
- Add 500 µL of Liver Organoid Initiation Medium per well.

Note: Medium should be gently pipetted into the corner of the well away from the Cultrex UltiMatrix RGF BME/organoids to prevent their disruption.

- Return the plate containing organoid cultures to the cell culture incubator to promote organoid growth. Follow Liver Organoid Expansion Protocol.



Learn More

Scan the QR Code or Visit:

rndsystems.com/products/organoid-and-3d-culture-reagents



View Our Full Cultrex UltiMatrix Protocol Video

Scan the QR Code or Visit:

rndsystems.com/UltiMatrix

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners. 8553015712_0225