# Human HepatiCult™ Protocol Walkthrough

### Lecture 3

### Presenter

Kiran Bhullar Scientist, Scientific Support, Epithelial Cell Biology

**Note:** The images in these slides depict <u>representative</u> morphologies. Users may see some variation in their own cultures compared to these images. Deviation from the morphology shown here does not necessarily indicate that an experiment will fail.

Magnifications (2X, 4X, 10X, 40X) refer to the objective magnification used.





# **Learning Objectives**

After this session, you should be able to:



- Apply confidently the key techniques and methodological considerations for hepatic organoid cultures
- Describe morphological features of cultures through different stages of the workflow

Property of StEM



### **Outline**

- Human HepatiCult<sup>™</sup> Expansion Workflow 1.
- Property of STERMCELL Technologies Human HepatiCult<sup>™</sup> Differentiation Workflow 2.



### Section 1 | Human HepatiCult™ Expansion Workflow



# HepatiCult<sup>™</sup> Workflow: Initiation from Tissue-Derived Hepatic Ducts



# **Media Preparation**



### Complete HepatiCult<sup>™</sup> Organoid Initiation Medium (OIM)

- HepatiCult<sup>™</sup> Organoid Basal Medium + Organoid Supplement + HepatiCult<sup>™</sup> Organoid Growth Supplement + Y-27632 (final concentration 10 µM) + antibiotics (e.g. final concentration 50 µg/mL gentamicin)
- Thaw the supplements overnight (O/N) at 2 8°C. If not used immediately, aliquot and store at -20°C. After thawing, use immediately. Don't re-freeze
- Complete HepatiCult<sup>™</sup> OIM can be stored at 2 8°C for up to 2 weeks
- Warm the complete media to room temperature (RT) before use

### Complete HepatiCult<sup>™</sup> Organoid Growth Medium (OGM)

- HepatiCult<sup>™</sup> Organoid Basal Medium + HepatiCult<sup>™</sup> Organoid Growth Supplement + antibiotics (e.g. final concentration 50 µg/mL gentamicin)
- Thaw the supplement O/N at 2 8°C. If not used immediately, aliquot and store at -20°C. After thawing, use immediately. Don't re-freeze
- Complete HepatiCult<sup>™</sup> OGM can be stored at 2 8°C for up to 2 weeks
- Warm the complete media to room temperature (RT) before use



# **Tissue Sourcing for Liver**

### Considerations



- Tissue Type Normal vs Diseased
  - Fresh, unprocessed, normal (ideally non-fibrotic/cirrhotic) liver tissue is preferred
  - Cadaveric tissue and biopsy samples/resections can be used for organoid establishment
  - If working with diseased samples, digestion protocol will likely be different depending on the biomechanical properties and cellular composition of the tissue
  - Some diseased samples might be easier to grow than others

### • Sources

- Researcher/institutional collaborations and affiliations
- Biobanks or commercial sources (e.g. BiolVT)

### • Tissue Processing

- Tissue should be processed fresh, whenever possible
- Tissue can be cryopreserved for a later use with CryoStor® CS10
- Liver samples/biopsies are usually large, it may not be feasible to freeze down the entire section of tissue
  - Split up the tissue into small enough pieces for fitting into a cryovial for cryopreservation
- Organoid yield is not correlated to tissue sample size we have used tissue weighing <0.5g</li>
- Flash frozen tissue has not been tested



# **Hepatic Tissue Sources**

Hepatic Ti	ssue S	ources			dies	
	BiolVT	Discovery Life Sciences	US Biolabs Custom Collections	IIAM (Custom Request)	National Disease Research Interchange	ATCC
Normal Tissue	Yes	Yes	Yes	Yes	Yes	No
Cancer Tissue	Yes	Yes	Yes	Maybe	Yes	Yes
Other Diseases	No	No	<ul> <li>Alcoholic Steatohepatitis (ASH)</li> <li>Liver Fibrosis</li> <li>Nonalcoholic Steatohepatitis (NASH)</li> <li>Liver Cirrhosis</li> </ul>	<ul> <li>Nonalcoholic Steatohepatitis (NASH)</li> <li>Hepatitis</li> </ul>	Yes	No
	05	08				MCELL



## **Tissue Digestion**

### Before starting the digestion, prepare the following solutions:

**a. Wash Solution** = Fetal Bovine Serum (FBS) + DMEM/F-12 with 15 mM HEPES (*Store on ice*)

**b. Tissue Dissociation Cocktail** = Collagenase Type IV + DNase I Solution (1 mg/mL) + DMEM/F-12 with 15 mM HEPES (*Warm to 37*°C *before use*)

c. TrypLE<sup>™</sup> + BSA in water (Store on ice)

**d. AdvDMEM + DNase I** = DMEM/F-12 + BSA + DNase I Solution (1 mg/mL) + Y-27632 (*Store on ice*) Washing and mincing tissue

Digest the tissue into ductal material DNase I = Break up the DNA from dead cells, makes samples less 'sticky'

Hepatic ducts can be quite large in size, TrypLE<sup>™</sup> breaks them into more manageable sizes (not necessarily single cells), BSA reduces stickiness

Dilutes the TrypLE<sup>™</sup> enzymatic reaction, washing/resuspending pellet at the end of digestion



# Critical Steps: Tissue Digestion

### Key Points for Tissue Dissociation:

- Prior to enzymatic digestion  $\rightarrow$  keep everything cold (on ice)
- Digestion  $\rightarrow$  water bath (37°C)
  - Repeat the digestion steps (at least 4 times)
- echnologies Assess the size of the pellet to decide if additional digestion might be required
  - If digestion is complete, there should be no tissue pieces, only white floating hepatic ducts (looks like fibres)
  - Repeat additional digestion cycle if tissue pieces still remain. Keep repeating the digestion until tissue pieces have been completely dissociated into hepatic ducts, and there are no visible tissue pieces
- Optional filtration step  $\rightarrow$  using a 100 µm reversible cell strainer to separate the large ductal material from the smaller ductal material
- Optional ACK lysis step  $\rightarrow$  if there is RBC contamination, a thin layer of RBCs is visible
  - If the pellet looks red, perform lysis. If the pellet looks clear, ACK lysis can be skipped 0
- After digestion is complete  $\rightarrow$  keep ductal material cold (on ice)

ropert





#### **Important Points:**

- Perform a viable cell count from the liver digestion
- Use 15 mL falcon tubes to get a more "compact" pellet
- Leave 5 10 µL of supernatant behind to avoid aspirating off the pellet
- Important to keep the pelleted cells/tube "cold"
- Be careful to **not introduce bubbles** when resuspending pellet in Matrigel®

- Use pre-cooled pipette tips to minimize premature solidification of Matrigel®
- Only dispense to the first stop to **avoid bubbles** on the top of the dome.
- Don't pipette directly onto the domes when adding media
- Passage cultures before the lumen turns dark and organoids collapse (up to 2 weeks)



# **Before Passaging**

- Place a 24-well tissue culture-treated plate in a 37°C incubator for at least 1 hour. Plates can be left overnight in a 37°C incubator
- Thaw ~40 µL of Corning® Matrigel® on ice for each dome to be seeded. Keep Matrigel® on ice when handling to prevent it from solidifying
- Prepare the complete OGM media as per the Technical Manual. If already prepared, leave an aliquot at RT
- Assess the morphology of hepatic organoids



# Hepatic Organoid Initiation: Morphology









# **Optimal Hepatic Organoid Fragment Sizes for Passaging**

Recommended fragment sizes for passaging

# (10 µl droplet)

Additional trituration recommended









30 μm, triturate further 30 μm, usually single cells 30-100 μm, desired size

### Fragment generation and counting

- Generate uniformly sized fragments (recommended = 30 - 100 μm)
- Use cell strainers to size-filter fragments
- Properly mix fragment suspensions before sampling - via stirring or gentle vortexing
- Dilute fragment suspensions for counts if needed - important for accurate counts
- Fragment size can influence organoid size and culture uniformity
  - Larger fragments will yield organoids earlier than smaller fragments/single cells



# Representative Organoid Morphology: Post-thaw Time Course



00 um





Cultures passaged again on Day 6



### **Important Points:**

- Bubbles may form when seeding domes (red arrow), but should not impact organoid growth
- Average organoid size is influenced by the organoid density in the dome
- Later in the passage, organoids may collapse (blue box) and fuse (green arrow)
- Culture is ready to be passaged on Day 5 6 (the recommended density is 1000 fragments/well)



All images taken at 2X magnification

# Representative Organoid Morphology: Day 8 in HepatiCult™ OGM



Single cells

"Organoids" outside of the dome "Large" organoids visible (overgrown)

### **Important Points:**

- Optimal passaging day can vary for donors and cultures this culture should have been passaged 1-2 days earlier
- Average organoid size is influenced by the organoid density in the dome - larger organoid size likely related to a less densely seeded dome
- Organoids can 'push out' of the Matrigel® dome edges (red arrows)
- Collapsed organoids can 'shed' or extrude cells in addition to deflating (blue arrows)



# **Identifying When to Passage Hepatic Organoids**

D7/8+ D1 - 4 D4 - 8 Too early to passage Ready to passage Should have been passaged 1-2 days earlier

#### When to passage

Monitor hepatic progenitor organoids daily; they should be passaged before the lumen turns dark and organoids collapse (generally 7 - 10 days)



Ċ.

18

# **Summary**

- Hepatic progenitor organoids can be established from fresh or frozen liver tissues
- Hepatic organoids can be monitored daily; they should be passaged before the lumen turns dark and organoids collapse
- Cultures require passaging every 7 10 days
- Routine culture maintenance is recommended using fragment generation and counting (recommended = 30 - 100 μm)



### Section 2 | Human HepatiCult<sup>™</sup> Differentiation Workflow





#### 

# **Media Preparation**

### Complete HepatiCult<sup>™</sup> Organoid Differentiation Medium (ODM)

- HepatiCult<sup>™</sup> Organoid Basal Medium + HepatiCult<sup>™</sup> Organoid Differentiation Supplement + dexamethasone (final concentration 3 µM) + antibiotics (e.g. final concentration 50 µg/mL gentamicin)
- Thaw the supplement O/N at 2 8°C. If not used immediately, aliquot and store at -20°C. After thawing, use immediately. Don't re-freeze.

100165

- Complete HepatiCult<sup>™</sup> ODM can be stored at 2 8°C for up to 2 weeks
- Complete HepatiCult™ ODM is light-sensitive; minimize exposure to light
  - To minimize exposure to the light, on days when the medium is required, aliquot the required volume into a tube, and warm to RT before use
  - Minimize repeated exposure of the stock bottle to the light
  - The medium should ideally be warmed and used in a BSC without the lights on, and if being stored in a fridge that is opened frequently, wrapping in foil is a good suggestion





#### **Important Points:**

- Perform at least one passage after initiating organoids
- For the fragment count, calculate the volume required to transfer 2000 fragments/well
- Seed additional domes in complete HepatiCult™ OGM for use as progenitor organoid controls/references in downstream analyses
- Perform a full media change on Day 3 with the complete HepatiCult™ OGM
- On Day 5 post-seeding, proceed with the differentiation step with HepatiCult™ ODM





### **Important Points:**

- Aliquot the required volume of ODM and warm to RT
- Perform complete media change on Day 5 post-seeding
  - Aspirate off OGM from well. Wash domes by adding 750 µl of RT DMEM/F-12 + 15 mM HEPES
- Perform full media change every 3 days using complete RT ODM
- Full media change is performed on Days 8, 11, and 14
  - For assays investigating secretory functions using the spent media, perform a final media change on Day 13 (i.e. at least 48 hours before final harvest on Day 15)
  - ο If domes are loose, perform a partial media change (e.g 500 μl instead of 750 μl)
- On Day 15, harvest/process domes for downstream characterization/functional assays
- Harvest not recommended later than Day 15



# **Representative Morphology: Differentiation Time Course**

### HepatiCult™ OGM

Day 2 (OGM)



HepatiCult<sup>™</sup> Differentiation Medium (ODM)

Day 5 (ODM Day 0)



Day 10 (ODM Day 5)

Day 7 (ODM Day 2)



Day 15 (ODM Day 10)





# Representative Morphology: Differentiation Day 1-30



### Important Points:

00 um

- Organoids begin to undergo morphology changes rapidly when differentiated
- Thickened epithelia (white arrows), compacted organoid size, and shift from cystic to dense morphologies are evident as early as differentiation day 1
- Onset and extent of morphological changes can be donor dependent



# Representative Morphology: Differentiation Day 5









### **Important Points:**

- Organoid morphologies continue to change over the course of the differentiation
- Thickened epithelia, compacted organoid size (blue arrows), and shift from cystic to dense morphologies are more evident
- Cell debris (red arrow) and organoids that appear collapsed are often observed and do not negatively affect differentiation outcome
- Differentiating organoid morphologies are heterogeneous and may vary by donor
- A further 3-5 days of differentiation is recommended to generate organoids with mature hepatic functionality



# **Differentiation Optimization**

### Seeding density:

- Recommended seeding densities to test: 1000 2500 fragments per dome
- Optimize organoid differentiation protocol parameters for every donor and assay of interest

### Number of domes for differentiation assay:

- Recommended testing ranges: 3 6 wells per donor per differentiation
- Intracellular protein- and nucleic acid-based analyses require more differentiated organoid material
- Assays investigating enzymatic activity or secretory functions (spent media) can be performed on fewer wells
- Application dependent: Important to seed at appropriate density for desired downstream application

### Duration of culture:

- Recommended testing range: Culture in complete HepatiCult™ OGM for 3 5 days
- Organoids should be differentiated once they are expanding, but before they start collapsing or deflating
- 14 15 days is the upper limit for differentiation duration



20100182

# Differentiation Optimization - Dome Disintegration

Technical Matrigel® dome handling and media change tips:

- Ensure protein content of Matrigel® lots being used is ≥8 mg/mL
- Ensure plates are being sufficiently pre-warmed at 37°C
- Remove as much of the supernatant from pellet fragments as possible when seeding domes (i.e keep Matrigel® content as close to 100% as possible for dome stability)
- Ensure medium is warmed before performing media changes on the domes
- Ensure dome is not disturbed during media changes.
  - Tilt the plate during media changes to visualize where the dome is
  - Perform slow media changes, removing and adding media to the wall of the well
  - For domes that look a little more loose, perform partial media changes, i.e. replace ~75% of the media in the well.
     Loose domes can still be used for analyses if media changes are performed carefully



# **Summary**

- Protocol optimization recommended for every donor to determine:
  - Seeding densities for passaging
  - Passaging frequency before organoids collapse

roperty

- Differentiation set-up
- A total of 1 2 passages is recommended prior to starting differentiation, functional applications and biobanking

echnologies

- Donor-to-donor variability observed in functional assays may be reflective of patient heterogeneity
  - Important to optimize experimental parameters for each donor and readout of interest





