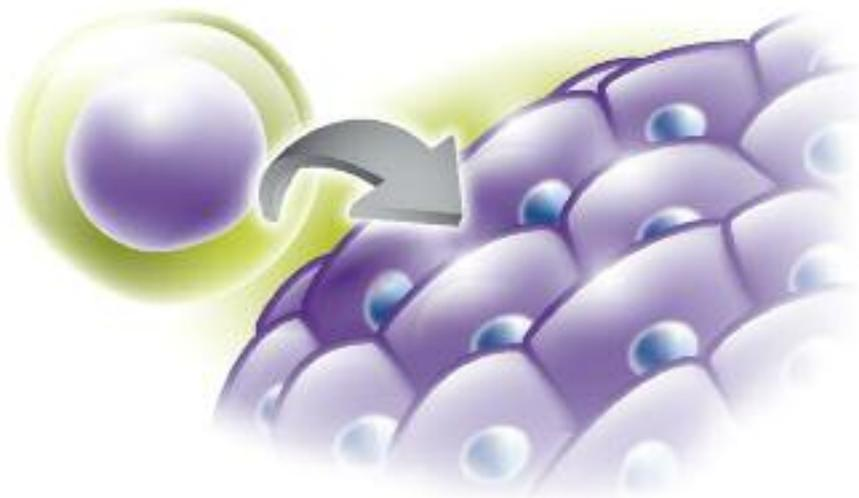


Stem cell differentiation protocol

abcam[®]
discover more

Human embryonic stem cells to hepatic endoderm



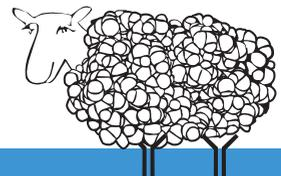
The ability to produce a homogenous population of functional cells in large quantities is one of the major challenges for stem cell biology. This is a robust 17 day differentiation protocol to direct human embryonic stem cells (hESCs) towards hepatocyte-like cells.

Edited protocol reproduced with kind permission of: David Hay and the Journal of Visualized Experiments

Taken from:

Robust Generation of Hepatocyte-like Cells from Human Embryonic Stem Cell Populations

Claire N. Medine, Baltasar Lucendo-Villarin, Wenli Zhou, Christopher C. West, David C. Hay. *J. Vis. Exp.* 2011 (56):e2969



Discover more at abcam.com/protocols

Section 1

Initial preparation of all chemical stocks and coating of culture plasticware

All steps to be carried out in a tissue culture hood under aseptic conditions.

1. Preparation of human basic Fibroblast Growth Factor (hbFGF)

- 1.1. Prepare 10% BSA solution in PBS and filter through a 0.22 μm filter.
- 1.2. From the 10% BSA solution prepare a 0.2% BSA solution.
- 1.3. Add 10 ml 0.2% BSA solution/100 μg hbFGF.
- 1.4. Pre-wet a 0.22 μm filter by filtering 5 ml 10% BSA solution through the filter. Discard the 10 ml BSA wash.
- 1.5. Filter the hbFGF through the pre-washed filter.
- 1.6. Aliquot the hbFGF in sterile Eppendorf tubes and store at -20°C .

2. Preparation of Human Activin A Stock Solution

- 2.1. Add 1 ml of 0.2% BSA into a syringe and pre-wet the filter.
- 2.2. Dilute the Activin A in 0.2% BSA to a stock concentration of 100 $\mu\text{g}/\text{ml}$.
- 2.3. Filter the Activin A solution and aliquot in sterile Eppendorf tubes, and store at -20°C .

3. Preparation of Mouse Wnt3a Stock Solution

- 3.1. Add 200 μl of PBS to a 2 μg vial of Wnt3a to a stock concentration of 10 $\mu\text{g}/\text{ml}$.
- 3.2. Aliquot in sterile Eppendorf tubes and store at -20°C .

4. Preparation of human Hepatocyte Growth Factor (HGF) Stock Solution (1000X)

- 4.1. Dilute the HGF in PBS to a stock concentration of 10 $\mu\text{g}/\text{ml}$.
- 4.2. Filter the HGF solution and aliquot in sterile Eppendorf tubes, and store at -20°C .

5. Preparation of Oncostatin M (OSM) Stock Solution (1000X)

- 5.1. Dilute the OSM in PBS to a stock concentration of 20 $\mu\text{g}/\text{ml}$.
- 5.2. Filter the OSM solution and aliquot in sterile Eppendorf tubes, and store at -20°C .

6. Coating of culture plasticware with Matrigel

- 6.1. Thaw the 10 ml stock bottle of Matrigel overnight at 4°C on ice and then add 10 ml of KO-DMEM (Knock Out-DMEM). Mix well using chilled pipettes and store 1 ml aliquots at -20°C .
- 6.2. Thaw an aliquot of Matrigel at 4°C for at least 2 hours or overnight to avoid the formation of a gel.
- 6.3. Add 5 ml of cold KO-DMEM to the Matrigel, mix well with a pipette.
- 6.4. Make up to 15 ml with cold KO-DMEM and mix using a pipette.
- 6.5. Add Matrigel to the plate or flask to be coated (see table below)

Recommended volumes of Matrigel for coating typical plasticware for hESC culture:

| Plate / Flask | Volume / Well or Flask |
|-------------------------|------------------------|
| 12 well plate | 0.5 ml per well |
| 6 well plate | 1 ml per well |
| 25cm ² plate | 2 ml per flask |

- 6.6. Incubate the coated plate or flask overnight at 4°C or room temperature for 1 hour before use.
- 6.7. Plates or flasks which have been coated with Matrigel can be stored at 4°C for up to 1 week. They should be clearly labelled with the date they were coated. Discard any plates or flasks not used within one week.
 - (a). Before use allow the coated culture container to come up to room temperature inside a tissue culture hood.
 - (b). Immediately prior to use aspirate the Matrigel and add the cell suspension to the well or flask.

Section 2

Preparation of media for differentiation of hESCs to hepatic endoderm

All media preparation should be carried out in a tissue culture hood under aseptic conditions.

1. Preparation of RPMI:B27 priming medium for endoderm differentiation

- 1.1. For RPMI-B27 medium, mix RPMI 1640 (500 ml) and B27 (50x, 10 ml).
- 1.2. Swirl to mix components.
- 1.3. Add all components to a filter unit and filter under vacuum, store at 4°C.

2. Preparation of SR-DMSO medium for hepatocyte differentiation

- 2.1. For SR-DMSO medium, mix 80% KO-DMEM, 20% Knock Out Serum Replacement medium (KO-SR), 0.5% L-glutamine, 1% non-essential amino acids, 0.1 mM β -Mercaptoethanol and 1% DMSO.
- 2.2. Filter the solution under vacuum, store at 4°C and aliquot and store at -20°C if required.
- 2.3. Use 4 ml per well of a 6-well plate, and 6 ml per T25 flask.

3. Preparation of L15 maturation medium for hepatocyte maturation

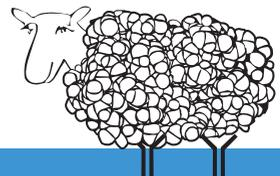
- 3.1. For L-15 medium, mix 500 ml Leibovitz L-15 medium, tryptose phosphate broth (final concentration 8.3%), heat inactivated foetal bovine serum (final concentration 8.3%), 10 μ M hydrocortisone 21-hemisuccinate, 1 μ M Insulin (bovine pancreas), 1% L-Glutamine, 0.2% ascorbic acid.
- 3.2. Filter the solution under vacuum, store at 4°C and aliquot and store at -20°C if required.

4. Preparation of final RPMI:B27 priming medium

- 4.1. Dispense the required volume of priming medium for the experiment (1 ml per well of a 6-well plate, and 2 ml per T25 flask).
- 4.2. Add Activin A to a final concentration of 100 ng/ml.
- 4.3. Add recombinant Wnt3a to a final concentration of 50 ng/ml.
- 4.4. Mix well and the media is now ready for use.
- 4.5. This final media should be made up fresh each day.

5. Preparation of final L-15 maturation medium

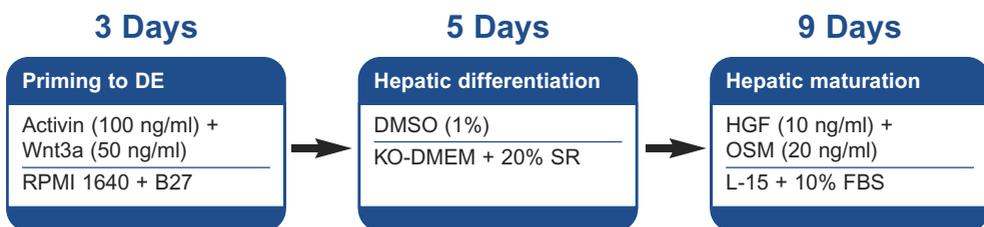
- 5.1. Dispense the required volume of L-15 medium for the experiment (4 ml per well of a 6-well plate, and 6 ml per T25 flask).
- 5.2. Add HGF to a final concentration of 10 ng/ml.
- 5.3. Add OSM to a final concentration of 20 ng/ml.
- 5.4. Mix well and the media is now ready for use.
- 5.5. This final media should be made up fresh each day.



Section 3

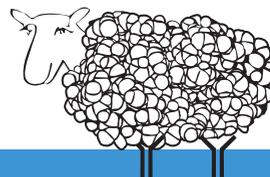
Priming hESCs to definitive endoderm

1. Culture hESCs (H1, H9 and RCM-1) and propagate on Matrigel coated plates with mouse embryonic fibroblast MEF-CM (mouse embryonic fibroblast culture media) supplemented with bFGF.
2. Initiate hepatic differentiation when hESCs reach a confluency level of approximately 30%-60% (depending on the hESC line) by replacing the MEF-CM with priming medium (RPMI 1640-B27 supplemented with 100 ng/ml Activin A and 50 ng/ml Wnt3a).
3. The cells are cultured in priming medium for 3 days (changing the medium every 24 hours), and final priming medium with Activin A and Wnt3a is made up fresh each day.
4. After 72 hours in priming medium, change the medium to the second differentiation medium (SR-DMSO) for 5 days (changing the medium every 48 hours).
5. At day 8, culture the cells in maturation and maintenance medium (L-15) supplemented with 10 ng/ml HGF and 20 ng/ml OSM for 9 days (changing medium every 48 hours). Maturation and maintenance medium with HGF and OSM is made up fresh each day.
6. The cells gradually exhibit morphological changes from a spiky/triangular shape to a characteristic liver morphology displaying a polygonal appearance.
7. Schematic for hepato-cellular differentiation:



Notes:

1. All priming, differentiation and maturation media are filtered under vacuum before use.
2. Priming, differentiation and maturation media are stored at 4°C for no longer than 2 weeks. Assess how much medium is required for the experiment and aliquot the remaining media and store at -20°C for future use.
3. Matrigel is made up as per the manufacturer's instructions; 1 ml aliquots can be stored at -20°C until use.
4. Growth factors once made up and aliquoted can be stored at -20°C and when thawed can be stored at 4°C for no longer than 2 weeks.



Discover more at abcam.com/protocols