

iCell® Retinal Pigment Epithelial Cells

Handling and Storage

Upon receipt, immediately transfer the cryovial to liquid nitrogen storage.

Preparing Cell Culture Surfaces

For best results, use vitronectin-coated vessels. Alternatively, tissue culture-treated vessels can be directly plated with cells in serum-containing medium.

1. Prepare 2.5 µg/ml vitronectin. For example to prepare 1 ml, combine:
 - 990 µl CellAdhere Dilution Buffer (STEMCELL Technologies, # 07183)
 - 10 µl of 250 µg/ml Vitronectin XF (STEMCELL Technologies, # 07180)
2. Coat vessel by adding the recommended coating volume per well (**Table 1**).
3. Incubate for ≥1 hour at room temperature.
4. Remove coating solution immediately before plating the cells (no rinse needed).

Preparing the Medium

Cells can be plated and cultured in either serum-free or serum-containing medium.

1. Prepare medium (**Table 2**); sterile filter using a 0.2 µm PES filter unit.
2. Store medium at 4°C for up to 2 weeks.

Thawing the Cells

1. Warm 25 ml of medium to room temperature.
2. Dispense 8 ml of medium into sterile 15 ml centrifuge tube.
3. Thaw the cryovial in a 37°C water bath for 3 minutes; clean with 70% ethanol.
4. Transfer the cells to the centrifuge tube containing the 8 ml of medium.
5. Rinse the cryovial with 1 ml of medium and transfer to centrifuge tube.
6. Centrifuge the cells at 300 x g (~1,000 rpm) for 5 minutes; discard the supernatant.

Plating the Cells

1. Check the Certificate of Analysis to obtain the number of expected cells.
2. Resuspend the cells at ~0.5 x 10⁶ cells/ml.
3. Add the cells to vessel using the recommended culture volume per well (**Table 1**).
4. Incubate the cells at 37°C, 5% CO₂.

Replacing the Medium

Feed the cells every 2 days by replacing the culture volume (**Table 1**) with an aliquot warmed to room temperature.

For most applications, the cells should be cultured for a minimum of 21 - 28 days (**Figure 2**).

Note: The cells are for LIFE SCIENCE RESEARCH USE ONLY. See www.cellulardynamics.com/product-warranty/ for USE RESTRICTIONS applicable to the cells and other terms and conditions related to the cells and their use.

Contacting Technical Support

Email: support@cellulardynamics.com

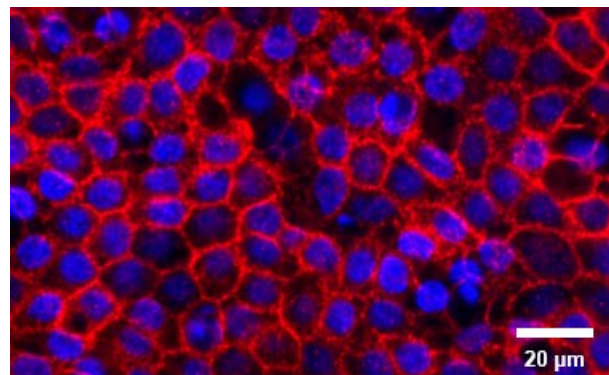


Figure 1: iCell Retinal Pigment Epithelial Cells were cultured for 31 days in serum-free medium: mature RPE marker BEST1 (red) and nuclei (blue).

Culture Vessel	Surface Area (cm ²)	Coating Volume (ml)	Culture Volume (ml)	Cell Number (cells)
6-well	9.5	2	3	1.5 x 10 ⁶
12-well	3.8	0.8	1.2	0.6 x 10 ⁶
12-well Transwell	1.12	0.24	0.35	0.175 x 10 ⁶
24-well	1.9	0.4	0.6	0.30 x 10 ⁶
48-well	0.95	0.2	0.3	0.15 x 10 ⁶
96-well	0.32	0.07	0.1	0.05 x 10 ⁶

Table 1: Cell Culture Volumes and Measures (per well)

Component	Volume (ml)	Final Concentration
MEM alpha ThermoFisher, # 12571-063	93.3	91.3%
KnockOut SR* ThermoFisher, # 10828-028	5	5%
N-2 Supplement ThermoFisher, # 17502-048	1	1%
Hydrocortisone, 50 µM Sigma, # H6909	0.11	55 nM
Taurine Sigma, # T0625 → prep 50 mg/ml**	0.5	250 µg/ml
Triiodo-L-thyronine (T ₃) Sigma, # T5516 → prep 20 µg/ml** Dilute 1:1,000 immediately before use	0.07	14 pg/ml
Gentamicin, 50 mg/ml ThermoFisher, # 15750-060 → optional	0.05	25 µg/ml

Table 2: Medium Preparation (adapted from JoVE 45, e2032)

* Alternatively, 5% fetal bovine serum can be used.

** Follow manufacturer's guidelines; stock solution may be aliquoted and frozen.

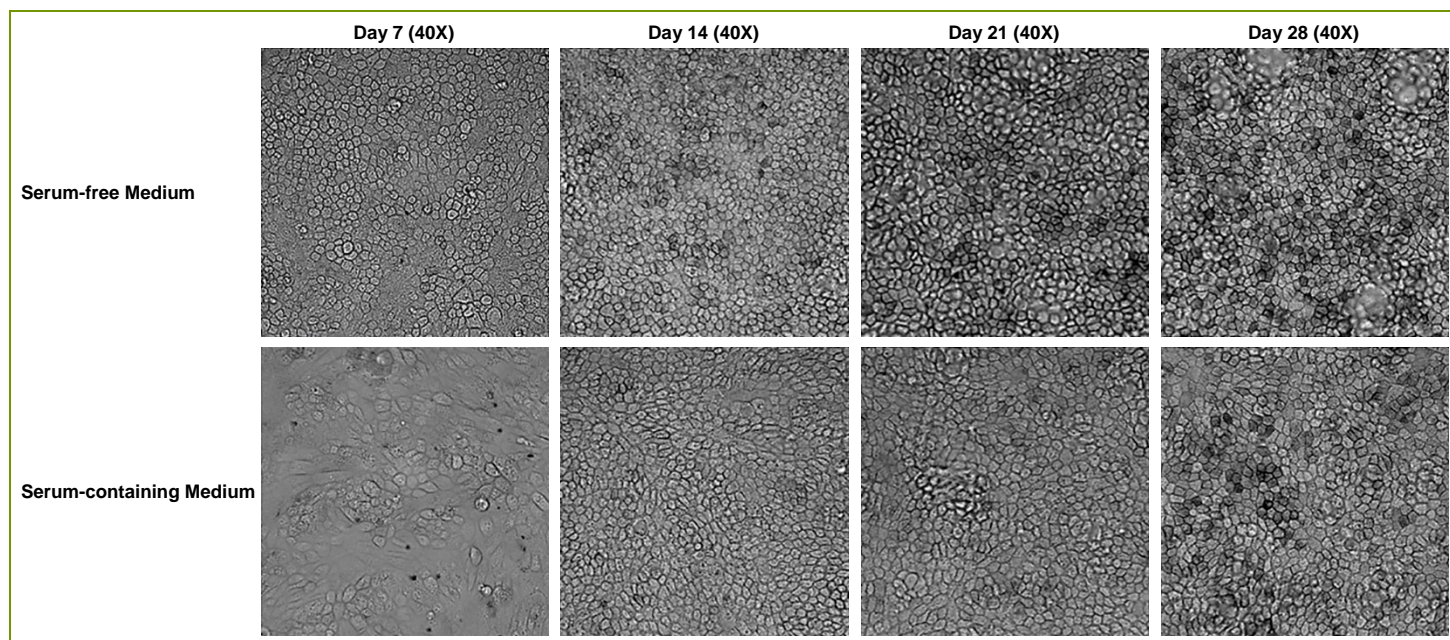



Figure 2: iCell Retinal Pigment Epithelial Cells can be cultured under serum-free or serum-containing conditions to form a tight monolayer with polygonal cell morphology that becomes increasingly pigmented with time in culture. Images were taken with a 40X objective.

Conditions of Use

The cells are for LIFE SCIENCE RESEARCH USE ONLY. See www.cellulardynamics.com/product-warranty/ for USE RESTRICTIONS applicable to the cells and other terms and conditions related to the cells and their use.

Trademarks

iCell and MyCell are registered trademarks, and Cellular Dynamics and the  logo are trademarks of Cellular Dynamics International, Inc. All other brands, product names, company names, trademarks, and service marks are the properties of their respective owners.

Copyright Notice

© 2017 Cellular Dynamics International, Inc. All rights reserved. This document may not be reproduced, distributed, modified or publicly displayed without the prior express written permission of Cellular Dynamics International, Inc.

Revision History

Document ID: X1015

Version 1.2: November 2017