



DISTRIBUTED BY



# **iCell® Hepatocytes 2.0**

## **User's Guide**

Document ID: X1007



## Trademarks

iCell and MyCell are registered trademarks, and Cellular Dynamics and the  logo are trademarks of Cellular Dynamics International, Inc. and may not be used without the express written permission of Cellular Dynamics International, Inc. (CDI).

All other brands, product names, company names, trademarks, and service marks are the properties of their respective owners.

## Restrictions and Liabilities

This document is provided “as is.” CDI assumes no responsibility for any typographical, technical, or other inaccuracies in this document. CDI reserves the right to periodically change information that is contained in this document; however, CDI makes no commitment to provide any such changes, updates, enhancements, or other additions to this document to you in a timely manner or at all.

OTHER THAN THE EXPRESS LIMITED WARRANTY SET FORTH IN THIS USER’S GUIDE, CDI MAKES NO REPRESENTATIONS, WARRANTIES, CONDITIONS OR COVENANTS, EITHER EXPRESS OR IMPLIED (INCLUDING WITHOUT LIMITATION, ANY EXPRESS OR IMPLIED WARRANTIES OR CONDITIONS OF FITNESS FOR A PARTICULAR PURPOSE, NON-INFRINGEMENT, MERCHANTABILITY, DURABILITY, TITLE, OR RELATED TO THE PERFORMANCE OR NON-PERFORMANCE OF ANY PRODUCT REFERENCED HEREIN OR PERFORMANCE OF ANY SERVICES REFERENCED HEREIN).

This document might contain references to third-party sources of information, hardware or software, products or services and/or third-party websites (collectively the “Third-Party Information”). CDI does not control, and is not responsible for, any Third-Party Information, including, without limitation the content, accuracy, copyright compliance, compatibility, performance, trustworthiness, legality, decency, links, or any other aspect of Third-Party Information. The inclusion of Third-Party Information in this document does not imply endorsement by CDI of the Third-Party Information or the third party in any way.

CDI does not in any way guarantee or represent that you will obtain satisfactory results from using iCell Hepatocytes 2.0 as described herein. The only warranties provided to you are included in the Limited Warranty enclosed with this guide. You assume all risk in connection with your use of iCell Hepatocytes 2.0.

## Conditions of Use

iCell Hepatocytes 2.0 are for life science research use only and subject to the use restrictions contained in Appendix A. You are responsible for understanding and performing the protocols described within this guide. CDI does not guarantee any results you may achieve. These protocols are provided as CDI’s recommendations based on its use and experience with iCell Hepatocytes 2.0.

## Origin

iCell Hepatocytes 2.0 are manufactured in the United States of America.

## Copyright Notice

© 2018 Cellular Dynamics International, Inc. All rights reserved. This document may not be reproduced, distributed, modified or publicly displayed without the express written permission of Cellular Dynamics International, Inc.

Copyright © 2018 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, and Scientists Helping Scientists are trademarks of STEMCELL Technologies Canada Inc. All other trademarks are the property of their respective holders. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

## Revision History

Document ID: X1007  
Version 1.0: July 2018

# Table of Contents

<b>Before You Begin</b> .....	<b>ii</b>
<b>Chapter 1. Introduction</b> .....	<b>1</b>
Components Supplied by Cellular Dynamics.....	2
Technical Support, Knowledge Base, and Training.....	3
Workflow Diagram .....	4
<b>Chapter 2. Handling and Storage</b> .....	<b>5</b>
Handling iCell Hepatocytes 2.0.....	5
Handling iCell Hepatocytes 2.0 Medium Supplement.....	5
<b>Chapter 3. Preparing Cell Culture Surfaces</b> .....	<b>6</b>
Recommended Vessel: Using the Pre-coated Cell Culture Vessel .....	6
Alternative Vessel: Preparing the Collagen I Cell Culture Vessel.....	6
<b>Chapter 4. Preparing Media</b> .....	<b>7</b>
<b>Chapter 5. Thawing iCell Hepatocytes 2.0</b> .....	<b>8</b>
<b>Chapter 6. Plating iCell Hepatocytes 2.0</b> .....	<b>9</b>
<b>Chapter 7. Maintaining iCell Hepatocytes 2.0</b> .....	<b>10</b>
<b>Appendices</b> .....	<b>11</b>
Appendix A. Intellectual Property Rights, Use Restrictions, and Limited License .....	11
Appendix B. Limited Warranty .....	11
Appendix C. Limited Liability.....	13

## Before You Begin

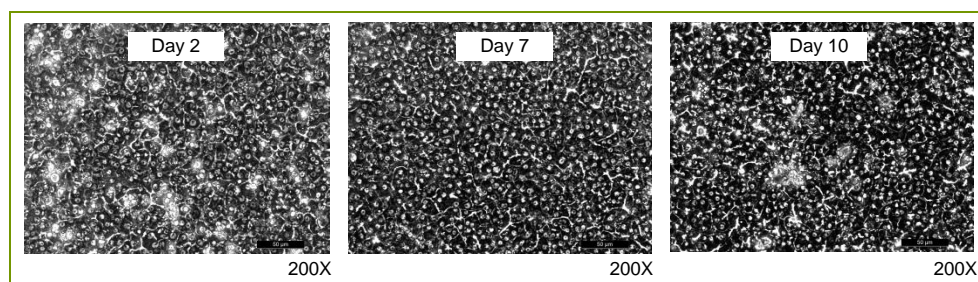
- Immediately transfer the frozen vial(s) to liquid nitrogen storage.
- Read this entire iCell® Hepatocytes 2.0 User's Guide before handling or using iCell Hepatocytes 2.0.
- iCell Hepatocytes 2.0 are for life science research use only. See Appendix A for more information and other restrictions.
- A Safety Data Sheet (SDS) for dimethyl sulfoxide (DMSO), in which iCell Hepatocytes 2.0 are frozen, is available online at [www.cellulardynamics.com/lit/](http://www.cellulardynamics.com/lit/) or on request from Cellular Dynamics International. Only technically qualified individuals experienced in handling DMSO and human biological materials should access, use, or handle iCell Hepatocytes 2.0.

Notes

## Chapter 1. Introduction

Cellular Dynamics International's (CDI) iCell Hepatocytes 2.0 are a highly pure population of human hepatocytes derived from induced pluripotent stem (iPS) cells using CDI's proprietary differentiation and purification protocols. iCell Hepatocytes 2.0 provide, for the first time, access to commercial quantities of high quality, high purity human liver cells for preclinical drug discovery, hepatotoxicity, and disease research.

When handled and maintained as recommended in this User's Guide, iCell Hepatocytes 2.0 exhibit a mature hepatocyte morphology, characteristic gene and protein expression, and active metabolism. iCell Hepatocytes 2.0 maintain their morphology and function for at least 14 days in culture making them amenable to a variety of assay endpoints.



**Figure 1: iCell Hepatocytes 2.0 Exhibit Mature Hepatocyte Morphology**

*These images show iCell Hepatocytes 2.0, 01434 at days 2, 7, and 10 post-plating. The cells display an adherent monolayer and cobblestone morphology within 48 hours and remain viable and adherent for an extended period in culture ( $\geq 14$  days). Some debris in the culture is expected but will decrease over time with medium replacements.*

## Components Supplied by Cellular Dynamics

Notes

Item	Catalog Number	
iCell Hepatocytes 2.0 Kit, 01434	R1027	
•iCell Hepatocytes 2.0, 01434 <sup>1, 2, 3</sup>	•C1026 ( $\geq 10.0 \times 10^6$ viable cells)	
•iCell Hepatocytes 2.0 Medium Supplement <sup>1, 3</sup>	•2 x M1024 (3 ml)	
•iCell Hepatocytes 2.0 User's Guide <sup>3</sup>	•X1007	
iCell Hepatocytes 2.0 Kit, 01279	R1025	R1104
•iCell Hepatocytes 2.0 Kit, 01279 <sup>2, 3</sup>	•C1023 ( $\geq 10.0 \times 10^6$ viable cells)	•C1090 ( $\geq 3.0 \times 10^6$ viable cells)
•iCell Hepatocytes 2.0 Medium Supplement <sup>1, 3</sup>	•2 x M1024 (3 ml)	•2 x M1024 (3 ml)
•iCell Hepatocytes 2.0 User's Guide <sup>3</sup>	•X1007	•X1007
Certificate of Analysis <sup>4</sup>		
Certificate of Origin		
If required for shipping purposes		

1 These products were formerly known by these names and/or catalog numbers:

iCell Hepatocytes 2.0, 01434 = iCell Hepatocytes 2.0 (Cat. No. PHC-100-020-001)

iCell Hepatocytes 2.0 Medium Supplement = iCell Hepatocytes 2.0 Medium Supplement (Cat. No. HCS-100-021-001)

**Note:** You may receive product labeled with its former name and/or catalog number until current stock is depleted. There were no changes in the manufacture of the cells or supplement.

2 iCell Hepatocytes 2.0, 01434 and iCell Hepatocytes 2.0, 01279 were derived from apparently healthy, normal donors. iCell Hepatocytes 2.0, 01434 are exclusive to CDI.

3 Safety Data Sheets and User's Guide available online: [www.fujifilmcdi.com/lit/](http://www.fujifilmcdi.com/lit/)

4 Available online: [www.fujifilmcdi.com/coa/](http://www.fujifilmcdi.com/coa/)

## Required Equipment and Consumables

Item	Vendor	Catalog Number(s)
<b>Equipment</b>		
37°C Water Bath	Multiple Vendors	
Biological Safety Cabinet with UV Lamp	Multiple Vendors	
Cell Culture Incubator	Multiple Vendors	
Hemocytometer or Automated Cell Counter <sup>1</sup>	Multiple Vendors	
Liquid Nitrogen Storage Unit	Multiple Vendors	
Pipettors	Multiple Vendors	
Tabletop Centrifuge	Multiple Vendors	
<b>Consumables</b>		
B27 Supplement, 50X	Thermo Fisher Scientific	17504
Collagen I Multiwell Plates, Biocoat <sup>2</sup>	Becton Dickinson	354408 (24-well) 354407 (96-well)
Conical Tubes, 15 ml, Falcon (Centrifuge Tubes) <sup>3</sup>	STEMCELL Technologies	38009
Dexamethasone	Thermo Fisher Scientific	ICN19456125
Gentamicin	Thermo Fisher Scientific	15750
Oncostatin M	R&D Systems	295-OM

## Notes

Item	Vendor	Catalog Number(s)
RPMI	Thermo Fisher Scientific	11875
Serological Pipettes, 5, 10, 25 ml <sup>3</sup>	STEMCELL Technologies	38003, 38004, 38005
Trypan Blue <sup>3</sup>	STEMCELL Technologies	07050

1 Ensure the automated cell counter is appropriately calibrated before use.

2 CDI recommends using pre-coated cell culture vessels for culturing iCell Hepatocytes 2.0. However, instructions for manually coating cell culture vessels are provided in Chapter 3, Preparing Cell Culture Surfaces.

3 Similar products are available from multiple vendors.

---

## Product and Scientific Support

Please contact STEMCELL's Product and Scientific Support for any technical inquiries or questions related to plating, media, cell culture and general assay methods.

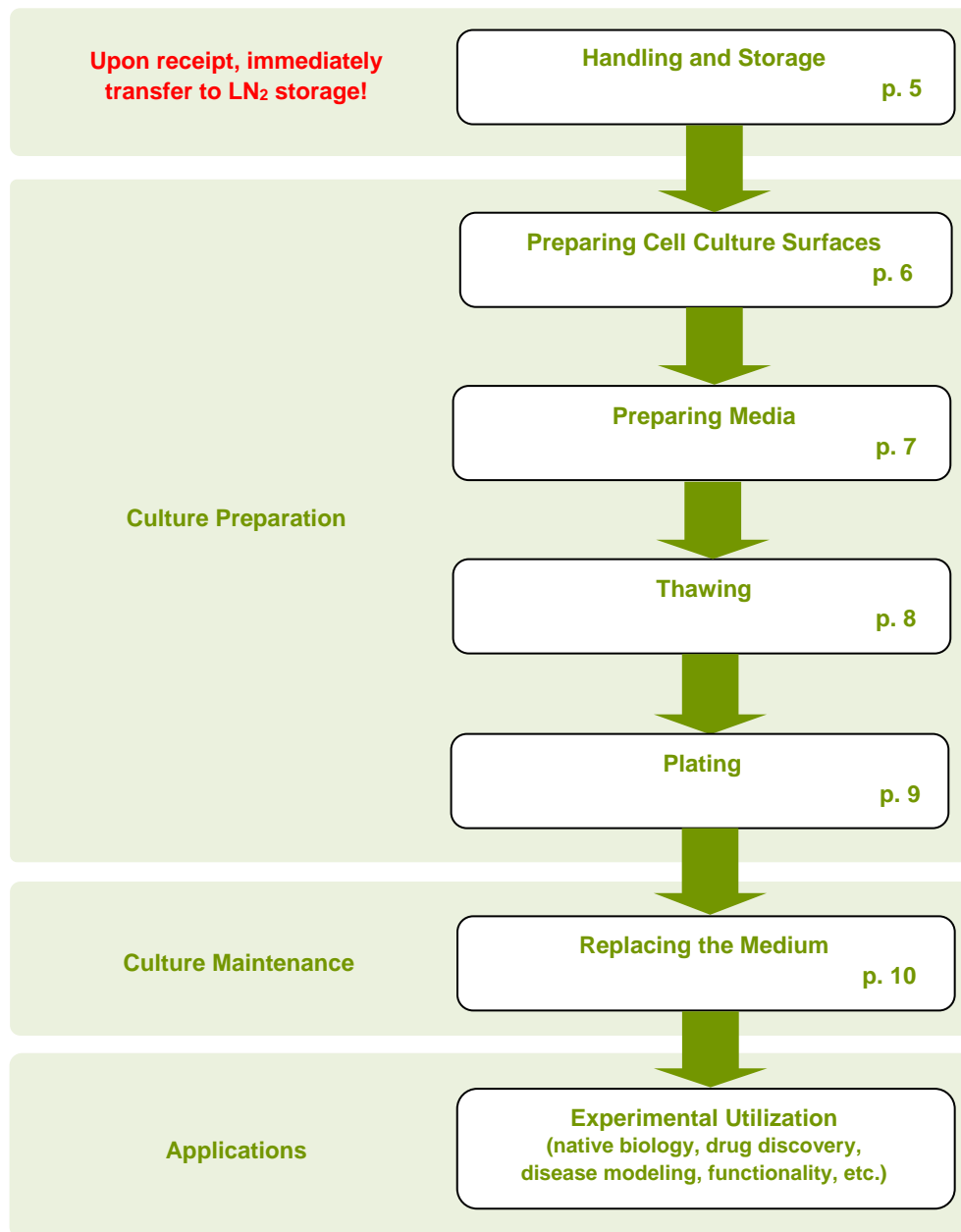
**Telephone** 1 800 667 0322

**Email** [techsupport@stemcell.com](mailto:techsupport@stemcell.com)

**Website** [www.stemcell.com](http://www.stemcell.com)

## Workflow Diagram

Notes





## Chapter 2. Handling and Storage

---

### Handling iCell Hepatocytes 2.0

iCell Hepatocytes 2.0 are provided as cryopreserved single-cell suspensions in 1.5 ml cryovials. Upon receipt, directly transfer the individual vials containing iCell Hepatocytes 2.0 to the vapor phase of a liquid nitrogen storage dewar.

**Note:** Do not store iCell Hepatocytes 2.0 on dry ice or at  $-80^{\circ}\text{C}$  as this may impact their performance post-thawing.



It is critical to maintain cryopreserved iCell Hepatocytes 2.0 at a stable temperature. Minimize exposure of cryopreserved iCell Hepatocytes 2.0 to ambient temperature when transferring vials to liquid nitrogen storage.

---

### Handling iCell Hepatocytes 2.0 Medium Supplement

iCell Hepatocytes 2.0 Medium Supplement is provided in 3 ml tubes and shipped frozen on dry ice. Upon receipt, store iCell Hepatocytes 2.0 Medium Supplement at  $-80^{\circ}\text{C}$  until ready for use.

## Chapter 3. Preparing Cell Culture Surfaces

iCell Hepatocytes 2.0 will plate and function optimally on collagen I-coated cell culture vessels. You can purchase pre-coated vessels (recommended) or coat them manually. Regardless of the coating methodology, have the plating surfaces ready before thawing iCell Hepatocytes 2.0.

### Recommended Vessel: Using the Pre-coated Cell Culture Vessel

See Required Equipment and Consumables section in Chapter 1, Introduction, for ordering information for recommended pre-coated cell culture vessels.

### Alternative Vessel: Preparing the Collagen I Cell Culture Vessel

1. Prepare 0.02 M acetic acid (Sigma, Cat. No. 320099) in sterile Milli-Q (or equivalent) water.



Observe safety precautions for handling concentrated acids. Wear eye protection and impervious gloves. Use caution and always add acid slowly to water, never the reverse.

2. Dilute 3 mg/ml collagen I solution (Gibco, Cat. No. A1048301) in 0.02 M acetic acid to a final concentration of 100 µg/ml.
3. Select the cell culture vessel appropriate for your experimental use. Use the volumes specified in the table below in the following coating procedure. Scale volumes appropriately for other vessel formats.

Culture Vessel	Volume of 100 µg/ml Collagen I Solution (ml)	Volume of Water Rinse (ml)
6-well Cell Culture Plate	1.0	2.0
12-well Cell Culture Plate	0.4	0.8
24-well Cell Culture Plate	0.2	0.4
48-well Cell Culture Plate	0.1	0.2
96-well Cell Culture Plate	0.04	0.08

**Table 1: Summary of Useful Volumes**

All volumes are **per well**.

4. Add the collagen I solution to each well of the vessel(s).
5. Tap the vessel(s) to distribute liquid evenly.
6. Incubate the vessel(s) at room temperature for at least 1 hour.
7. After incubation, completely aspirate the collagen I solution from each well. Rinse each well 3 times with sterile Milli-Q water and aspirate completely.

**Note:** If necessary, air-dry the vessels coated with collagen I and store at 4°C for up to 7 days before use. Equilibrate the vessel(s) at room temperature before plating iCell Hepatocytes 2.0.

## Chapter 4. Preparing Media

Prepare and store the Plating Medium and Maintenance Medium for iCell Hepatocytes 2.0 as follows:

Plating Medium <sup>1, 2, 3, 4</sup>		
Component	Amount (ml)	Final Concentration
RPMI	72	96%
B27 Supplement, 50X	1.5	1X
Oncostatin M, 10 µg/ml	0.15	20 ng/ml
Dexamethasone, 5 mM	0.0015	0.1 µM
Gentamicin	0.0375	25 µg/ml
iCell Hepatocytes 2.0 Medium Supplement	1.5	1X

- 1 Prepare the Plating Medium on the day of thaw to ensure optimal viability of cells upon plating.
- 2 Prepare stock solutions of 10 µg/ml oncostatin M and 5 mM dexamethasone. Follow the manufacturer's instructions for storage and reconstitution.
- 3 Prepare 10 ml aliquots of Plating Medium in 15 ml centrifuge tubes. Up to 2 vials of iCell Hepatocytes 2.0 can be thawed into 1 tube of 10 ml Plating Medium.
- 4 Store the Plating Medium at 4°C for up to 1 week. Do not store at -20°C.

Maintenance Medium*		
Component	Amount (ml)	Final Concentration
RPMI	72	96%
B27 Supplement, 50X	1.5	1X
Dexamethasone, 5 mM	0.0015	0.1 µM
Gentamicin	0.0375	25 µg/ml
iCell Hepatocytes 2.0 Medium Supplement	1.5	1X

- \* Store the Maintenance Medium at 4°C for up to 1 week. Do not store at -20°C.

## Chapter 5. Thawing iCell Hepatocytes 2.0

Maintain iCell Hepatocytes 2.0 in liquid nitrogen until immediately before thawing to ensure maximal performance of the cells. Complete the following steps of the thawing procedure in a time-efficient manner to facilitate optimal iCell Hepatocytes 2.0 viability and performance.

**Note:** Thaw no more than 2 vials of iCell Hepatocytes 2.0 at one time.

1. Equilibrate a 10 ml aliquot of Plating Medium in a 37°C water bath before thawing iCell Hepatocytes 2.0.
2. Remove the iCell Hepatocytes 2.0 cryovial from the liquid nitrogen storage tank.

**Note:** If necessary, place cryovials on dry ice for up to 10 minutes before thawing.

3. Immerse the cryovial in a 37°C water bath for **exactly 3 minutes** (avoid submerging the cap) holding the tube stationary (no swirling). Use of a floating microcentrifuge tube rack is recommended.



Precise timing is critical to maximizing viable cell recovery and attachment.

4. Immediately remove the cryovial from the water bath, spray with 70% ethanol, and place in a biological safety cabinet.
5. Gently transfer the iCell Hepatocytes 2.0 cryovial contents into the 15 ml centrifuge tube containing 10 ml of 37°C Plating Medium.



Avoid repeated pipetting of the thawed iCell Hepatocytes 2.0 cell suspension.

6. Rinse the empty iCell Hepatocytes 2.0 cryovial with 1 ml of 37°C Plating Medium to recover any residual cells from the vial. Transfer the 1 ml of Plating Medium rinse from the cryovial to the 15 ml centrifuge tube containing the iCell Hepatocytes 2.0 cell suspension. Invert the tube slowly to mix the cell suspension.
7. Centrifuge the cell suspension at 200 x g for 3 minutes at room temperature.
8. Carefully aspirate the supernatant, taking care not to disturb the cell pellet.
9. Slowly add 2 ml of 37°C Plating Medium with a wide-bore pipette to resuspend the cell pellet. Avoid vigorous shaking or vortexing of the cell suspension.

**Note:** Avoid repeated pipetting of the thawed iCell Hepatocytes 2.0 to ensure maximum viability and cell attachment.

**Note:** Thaw no more than 2 vials of iCell Hepatocytes 2.0 at one time. Once thawed, pour the contents of the 2 vials into the same 15 ml centrifuge tube containing 10 ml of 37°C Plating Medium and rinse each vial before inverting the tube.

## Chapter 6. Plating iCell Hepatocytes 2.0

The recommended plating density of iCell Hepatocytes 2.0 is  $3 \times 10^5$  viable cells/cm<sup>2</sup>. Adjust the cell concentration in Plating Medium as needed.

1. Obtain the number of viable cells/vial and viability from the Certificate of Analysis.
2. Remove a sample of cells to confirm viability using a hemocytometer (using trypan blue exclusion to identify viable cells) or an automated cell counter.

**Note:** Use a hemocytometer until validation of the automated cell counter can be completed as cell debris may cause altered viability values as calculated on automated counters.

3. Calculate the final volume of Plating Medium needed to obtain a desired cell plating density (i.e.  $1 \times 10^6$  cells/ml) using the number of viable cells/vial from the Certificate of Analysis.
4. Dispense the cell suspension into the appropriate cell culture vessel(s). Use the volumes specified in the table below for the selected vessel format. For 24-well cell culture plates or larger, gently shake the plates to distribute the cells evenly. Do not shake 96-well cell culture plates.

Culture Vessel	Surface Area (cm <sup>2</sup> )	Plating Volume (ml)	Cell Number ( $3 \times 10^5$ cells/cm <sup>2</sup> )
6-well Cell Culture Plate	9.5	2.9	$2.9 \times 10^6$
12-well Cell Culture Plate	3.8	1.1	$1.1 \times 10^6$
24-well Cell Culture Plate	1.9	0.57	$5.7 \times 10^5$
48-well Cell Culture Plate	0.95	0.29	$2.9 \times 10^5$
96-well Cell Culture Plate	0.32	0.096	$0.96 \times 10^5$

**Table 1: Summary of Recommended Volumes and Measures**

All volumes and measures are **per well**.

5. Culture iCell Hepatocytes 2.0 in a cell culture incubator at 37°C, 5% CO<sub>2</sub> for 3 - 4 hours.
6. During incubation, equilibrate an aliquot Plating Medium at room temperature.
7. Following incubation, shake 24-well cell culture plates in the diagonal planes 4 times to remove dead cells and debris. Do not shake 96-well cell culture plates.
8. Aspirate the Plating Medium from the plate using a pipettor and replace with room temperature Plating Medium.
 

**Note:** Do not aspirate the medium from more than 12 wells of any vessel format at one time to prevent wells from drying.
9. Culture iCell Hepatocytes 2.0 in a cell culture incubator at 37°C, 5% CO<sub>2</sub> overnight.

## Chapter 7. Maintaining iCell Hepatocytes 2.0

Notes

iCell Hepatocytes 2.0 are shipped cryopreserved at high purity. The hepatocytes preserve a high purity for at least 14 days after thawing if plated and maintained as recommended.



*Plating Medium and Maintenance Medium are stable for 1 week when stored at 4°C.*

1. Equilibrate the Plating Medium at room temperature before use.
2. Approximately 24 hours post-plating iCell Hepatocytes 2.0, aspirate the spent medium and replace (100% exchange) with the appropriate volume of Plating Medium. Recommended volumes are as follows:
  - **6-well cell culture plate:** 3.0 ml/well
  - **12-well cell culture plate:** 1.0 ml/well
  - **24-well cell culture plate:** 0.6 ml/well
  - **48-well cell culture plate:** 300  $\mu$ l/well
  - **96-well cell culture plate:** 100  $\mu$ l/well
3. Replace 100% of the spent Plating Medium daily until day 5 post-plating.
4. At day 5 post-plating, equilibrate the Maintenance Medium at room temperature before use.
5. Aspirate the spent Plating Medium and replace (100% exchange) with the appropriate volume of Maintenance Medium. Recommended volumes are the same as in step 2.
6. Replace 100% of Maintenance Medium every 2 days thereafter.
7. Culture iCell Hepatocytes 2.0 in a cell culture incubator at 37°C, 5% CO<sub>2</sub>.

## Appendices

---

### Appendix A. Intellectual Property Rights, Use Restrictions, and Limited License

A. **OWNERSHIP.** The Products are covered by pending patents and patents: [cellulardynamics.com/about-us/patents/](http://cellulardynamics.com/about-us/patents/). Customer has a limited license to use the Products for internal research purposes for the sole benefit of the Customer, subject to the use restrictions included in subsection B of this Appendix A. Customer acknowledges and agrees that the receipt or purchase of the Products by Customer shall not be construed as a transfer of any title or the grant of any rights in or to the intellectual property embodied in the Products owned or licensed by Cellular Dynamics. In particular, no right or license to make, have made, offer to sell, or sell the Products, to modify or reproduce the Product or any part thereof, or to use the Products in combination with any other product(s), except product(s) provided or expressly licensed to Customer by Cellular Dynamics for such use, is implied or conveyed by the sale or transfer of Products to Customer.

B. **USE RESTRICTIONS.** The Products are licensed for internal research purposes only, and may not be used for any other purpose. The Products must be used in accordance with this User's Guide to which Customer, by ordering and accepting the Products, agrees. Customer shall not make, have made, offer to sell, or sell the Products. Unless expressly permitted in a contract executed between Customer and Cellular Dynamics, Customer shall not use the Products (or any modifications Customer makes to the Products or any cells derived, developed or expanded from the Products) in (i) the manufacture of any products, or (ii) any services for a third party. Customer may not transfer the Products (or any modifications Customer makes to the Products or any cells derived, developed or expanded from the Products) to any third party without Cellular Dynamics' prior written consent. Customer shall not reverse engineer the Products. Customer shall not use the Products, components or modifications thereof, or any cells derived, developed or expanded therefrom, in humans, in clinical trials, for diagnostic purposes involving human subjects, or for any investigational or other therapeutic use. Customer shall not use the Products directly or indirectly to derive or make any human gamete or gamete precursor cell. Customer shall use the Products in accordance with all applicable laws and regulations and any applicable institutional review board approved protocol and/or privacy office approval. Customer is not entitled to receive any data or information from Cellular Dynamics that directly identifies the donor of the biological materials from which the Products indirectly are derived or were made. Customer shall not attempt in any way to determine the identity of the donor of the biological materials from which the Products indirectly are derived or was made.

---

### Appendix B. Limited Warranty

A. During the Warranty Period (as defined below) and subject to subsection F of this Appendix B, Cellular Dynamics warrants that its Products conform to the specifications contained in the Certificate of Analysis for the Product shipped to Customer. Customer's sole and exclusive remedy (and Cellular Dynamics' sole and exclusive liability) with respect to any defective Products shall be replacement of the defective Products by Cellular Dynamics pursuant to this Appendix B.

B. Under no circumstances shall Cellular Dynamics' liability to Customer exceed the amount paid by Customer for the Products to Cellular Dynamics. Cellular Dynamics will bear all reasonable shipping costs if the Products are replaced pursuant to this warranty. For clarity, this warranty automatically shall be void, and any claims under it invalid, (i) if Customer's use of the Products is other than solely in accordance with this User's Guide and Cellular Dynamics' Terms and Conditions (or such other written agreement between Cellular Dynamics and Customer under which the Products are sold or transferred to Customer) or for a purpose or in a manner other than that for which the Products were designed; or (ii) if Customer fails to follow this User's Guide for the use, storage, and handling of the Products

however such failure is caused; or (iii) if Customer fails to comply with any of the provisions of Appendix A in this User's Guide; or (iv) if there is any abuse, other misuse or neglect of the Products by Customer or to the extent of any damage or loss of the Products by events or occurrences beyond a person's (e.g., Cellular Dynamics') control including without limitation, accident, fire, vandalism and natural disasters (acts of God). This warranty applies only to Customer and not to third parties. This warranty is not assignable.

C. TO THE FULLEST EXTENT PERMITTED BY APPLICABLE LAW, CELLULAR DYNAMICS DISCLAIMS ALL REPRESENTATIONS, AND WARRANTIES, EXPRESS OR IMPLIED, WITH RESPECT TO THE PRODUCTS, INCLUDING BUT NOT LIMITED TO, ANY IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE OR NON-INFRINGEMENT, AND CUSTOMER WAIVES ALL RIGHTS AND REMEDIES, WITH RESPECT TO ANY DEFECTIVE PRODUCTS OTHER THAN THE EXPRESS WARRANTY AND REMEDY THEREFOR STATED ABOVE IN THIS APPENDIX B.

D. Within five (5) business days of thawing the Product but prior to the expiration date of the Product as listed on the Certificate of Analysis and/or Product's label (the "Warranty Period"), Customer must notify Cellular Dynamics in writing of any nonconformity of the Products, describing the nonconformity in detail. Customer's failure to properly notify Cellular Dynamics in the Warranty Period voids the limited warranty set forth above in this Appendix B.

E. Customers who believe they have a warranty claim should call Cellular Dynamics' Technical Support line at (608) 310-5100 ext. 3 or email at support@cellulardynamics.com to request a replacement Product based on a breach of the limited warranty set forth above in this Appendix B. Any action by Customer for Cellular Dynamics' breach of this limited warranty, for which Customer has given timely and proper notice of such breach during the Warranty Period and otherwise in accordance with this Appendix B, must be commenced by Customer within 18 months following the date of such breach.

F. Cellular Dynamics makes no warranty of any kind or nature, neither express nor implied, for any product sold together with, or as a part of, the Products (e.g., an accessory accompanying a Product or a discrete component part of a Product that is a kit) that is not manufactured by Cellular Dynamics. Any such accessory to or part of the Products shall have the warranty, if any, that is offered and granted (and, for clarity, extended by its terms to Customer) by the manufacturer of such other accessory or component product accessories.

G. Customer acknowledges and agrees that Cellular Dynamics may fill Customer's order with any number of units of Products. Such units may be more units than Customer ordered. Customer will not be charged extra for any adjustments made by Cellular Dynamics. The number of cells in a unit is determined by the Product's Certificate of Analysis. The number of cells that are contained in a unit accounts for both viability and plating efficiency percentages. Because this may vary from lot to lot, Cellular Dynamics reserves the right to fill the order with that number of units which is sufficient to fill Customer's order and such adjustments shall not constitute a breach of the limited warranty set forth herein.



## Appendix C. Limited Liability

TO THE FULLEST EXTENT PERMITTED UNDER APPLICABLE LAW, CELLULAR DYNAMICS SHALL NOT HAVE ANY LIABILITY FOR INCIDENTAL, COMPENSATORY, PUNITIVE, CONSEQUENTIAL, INDIRECT, SPECIAL OR OTHER SIMILAR DAMAGES, HOWEVER CAUSED AND REGARDLESS OF FORM OF ACTION WHETHER IN CONTRACT, TORT (INCLUDING NEGLIGENCE), STRICT PRODUCT LIABILITY OR OTHERWISE, EVEN IF CELLULAR DYNAMICS HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES. NOTWITHSTANDING ANY OTHER TERM OR IMPLICATION TO THE CONTRARY, UNDER NO CIRCUMSTANCES SHALL CELLULAR DYNAMICS' LIABILITY TO CUSTOMER EXCEED THE AMOUNT PAID BY CUSTOMER FOR THE PRODUCTS TO CELLULAR DYNAMICS.

© 2018 Cellular Dynamics International, Inc. All rights reserved.





