




# **iCell<sup>®</sup> Mesenchymal Stem Cells Prototype User's Guide**

DOCUMENT #DX21838



## Trademarks

iCell is a registered trademark, 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.

THE iCELL MESENCHYMAL STEM CELLS ARE PROVIDED “AS IS” TO YOU. TO THE FULLEST EXTENT PERMITTED BY APPLICABLE LAW, CDI MAKES NO, AND DISCLAIMS ALL, REPRESENTATIONS, WARRANTIES, CONDITIONS OR COVENANTS, EXPRESS OR IMPLIED, WITH RESPECT TO ANY PRODUCT REFERENCED HEREIN OR PERFORMANCE OF ANY SERVICES REFERENCED HEREIN, INCLUDING BUT NOT LIMITED TO, 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 Mesenchymal Stem Cells as described herein. You assume all risk in connection with your use of iCell Mesenchymal Stem Cells.

## Conditions of Use

iCell Mesenchymal Stem Cells 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 Mesenchymal Stem Cells.

## Origin

iCell Mesenchymal Stem Cells are manufactured in the United States of America.

## Copyright Notice

© 2016 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.

## Revision History

Document #DX21838

Version 0.1: November 2016

# 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
Required Equipment and Consumables.....	2
Technical Support and Training.....	3
Workflow Diagram .....	4
<b>Chapter 2. Handling and Storage</b> .....	<b>5</b>
<b>Chapter 3. Preparing Media</b> .....	<b>6</b>
<b>Chapter 4. Thawing iCell Mesenchymal Stem Cells</b> .....	<b>7</b>
<b>Chapter 5. Plating iCell Mesenchymal Stem Cells</b> .....	<b>9</b>
<b>Chapter 6. Maintaining iCell Mesenchymal Stem Cells</b> .....	<b>10</b>
<b>Appendices</b> .....	<b>12</b>
Appendix A. Intellectual Property Rights, Use Restrictions, and Limited License...12	
Appendix B. Product Provided “AS IS” .....	12
Appendix C. Limited Liability.....	13

## Before You Begin

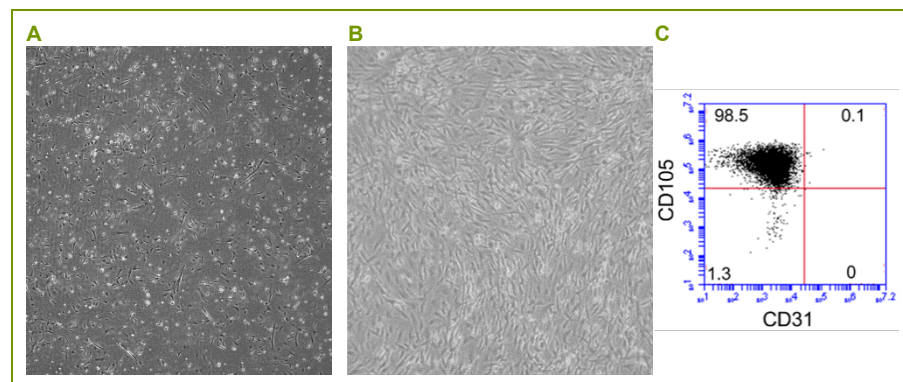
Notes

- Immediately transfer the frozen vials to liquid nitrogen storage.
- Read this entire iCell® Mesenchymal Stem Cells Prototype User's Guide before handling or using iCell Mesenchymal Stem Cells.
- iCell Mesenchymal Stem Cells 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 Mesenchymal Stem Cells 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 Mesenchymal Stem Cells.

## Chapter 1. Introduction

Cellular Dynamics International's (CDI) iCell Mesenchymal Stem Cells are a highly pure population of human mesenchymal stem cells derived from induced pluripotent stem (iPS) cells using CDI's proprietary differentiation protocols. iCell Mesenchymal Stem Cells exhibit characteristic gene and protein expression and multilineage differentiation potential (i.e. adipocyte, chondrocyte, and osteocyte). These cells provide a reliable source of human mesenchymal stem cells suitable for use in targeted drug discovery, toxicity testing, and other life science research.

When thawed and plated according to this User's Guide, iCell Mesenchymal Stem Cells can be maintained in culture for up to 5 passages. These cells are suitable for an array of assays including but not limited to flow cytometry, qPCR, and immunomodulatory assays.



**Figure 1: iCell Mesenchymal Stem Cells Represent a Highly Pure Population of Human Mesenchymal Stem Cells**

*These images show iCell Mesenchymal Stem Cells at (A) day 1 and (B) day 5 post-plating. iCell Mesenchymal Stem Cells form a confluent monolayer in culture in approximately 5 days post-plating and (C) show the appropriate human mesenchymal stem cell markers as demonstrated by flow cytometry.*

## Components Supplied by Cellular Dynamics

Notes

Item	Catalog Number
iCell Mesenchymal Stem Cells Prototype <sup>1</sup>	MSC-301-010-001-PT
iCell Mesenchymal Stem Cells Prototype User's Guide <sup>1</sup>	
Certificate of Testing <sup>2</sup>	
Certificate of Origin If required for shipping purposes	
<sup>1</sup> Safety Data Sheets and User's Guide available online at <a href="http://www.cellulardynamics.com/lit/">www.cellulardynamics.com/lit/</a> <sup>2</sup> Also available online at <a href="http://www.cellulardynamics.com/cot/">www.cellulardynamics.com/cot/</a>	

## Required Equipment and Consumables

**Note:** As required for the intended use, see the following iCell Mesenchymal Stem Cells Prototype Application Protocols for assay-specific equipment and consumables before thawing cells:

- Modeling Adipocyte Differentiation
- Modeling Chondrocyte Differentiation
- Modeling Osteocyte Differentiation

These Application Protocols are available online at [www.cellulardynamics.com/lit/](http://www.cellulardynamics.com/lit/).

Item	Vendor	Catalog Number
<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*	Multiple Vendors	
Liquid Nitrogen Storage Unit	Multiple Vendors	
Pipettors	Multiple Vendors	
Tabletop Centrifuge	Multiple Vendors	
<b>Consumables</b>		
1-Thioglycerol (MTG)	Sigma	M6145
15 ml Centrifuge Tubes	Multiple Vendors	
6-well Cell Culture Plates	Multiple Vendors	
10 cm Cell Culture Dish	Multiple Vendors	
Ascorbic Acid	Multiple Vendors	
B-27 Supplement, Minus Vitamin A	Life Technologies	12587
Bovine Serum Albumin (BSA)	Sigma	A1470
bFGF from Zebrafish	Multiple Vendors	
Dulbecco's Phosphate Buffered Saline without Ca <sup>2+</sup> and Mg <sup>2+</sup> (D-PBS)	Life Technologies	14190, 14040
Fetal Bovine Serum (FBS)	Multiple Vendors	
GlutaMAX Supplement	Life Technologies	35050
Ham's F12 Medium	Corning	10-080-CV
Iscove's Modified Dulbecco's Medium (IMDM)	Life Technologies	12440
N-2 Supplement	Life Technologies	17505
PDGF-BB	Multiple Vendors	
Penicillin/Streptomycin	Life Technologies	15140

## Notes

Item	Vendor	Catalog Number
PES Filter Unit, 0.2 µm, 500 ml	Multiple Vendors	
Trypan Blue	Life Technologies	15250
TrypLE Select Enzyme, No Phenol Red (TrypLE)	Life Technologies	12563

\* Ensure the automated cell counter is appropriately calibrated before use.

---

## Technical Support and Training

CDI's Technical Support Scientists have the necessary laboratory and analytical experience to respond to your inquiries. In addition, in-lab training may be available upon request.

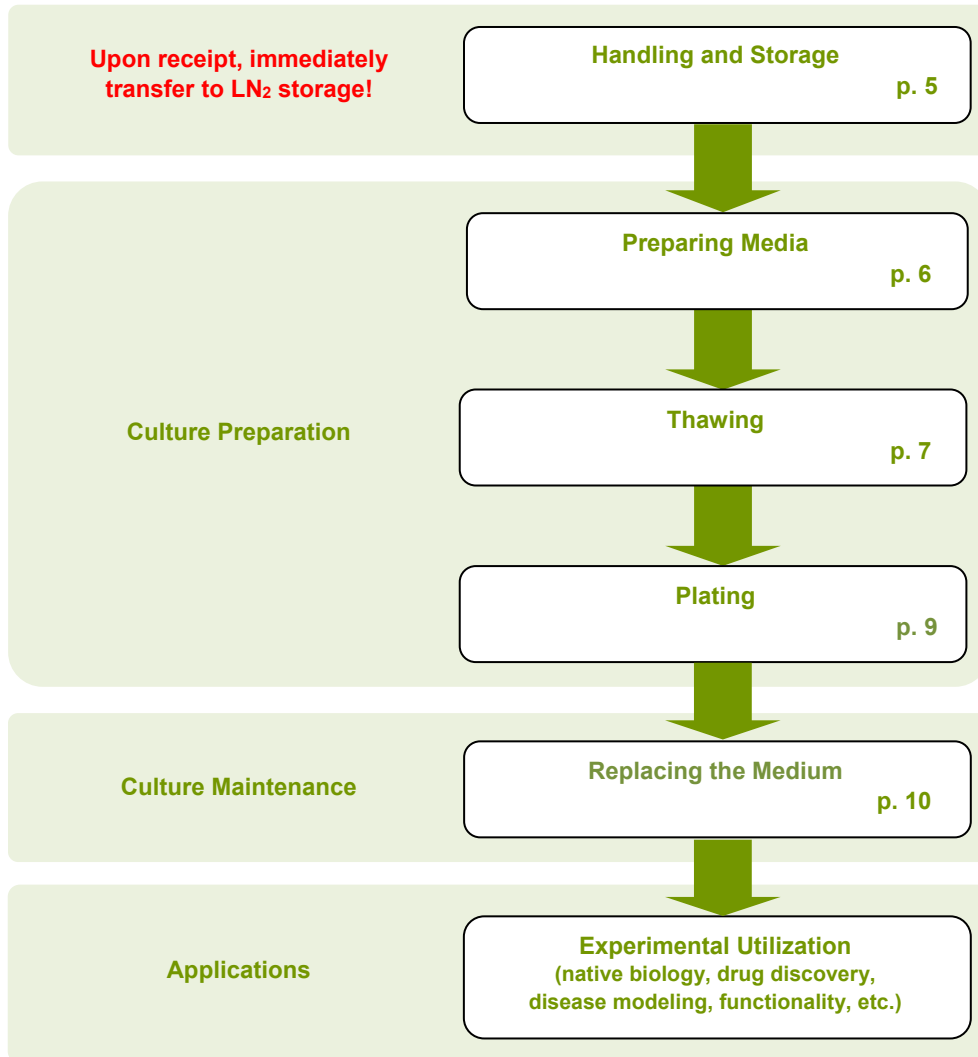
**Telephone** (877) 320-6688 (US toll-free) / (608) 310-5100 x5  
Monday - Friday, 8:30 am - 5:00 pm US Central Time

**Fax** (608) 310-5101

**Email** [support@cellulardynamics.com](mailto:support@cellulardynamics.com)

## Workflow Diagram

Notes





## Chapter 2. Handling and Storage

iCell Mesenchymal Stem Cells are provided as cryopreserved single-cell suspensions in 1.5 ml cryovials. Upon receipt, directly transfer the cryobox containing iCell Mesenchymal Stem Cells to the vapor phase of liquid nitrogen storage dewar. CDI strongly recommends transferring the entire cryobox into the storage rack to avoid transferring individual vials.



*It is critical to maintain cryopreserved iCell Mesenchymal Stem Cells at a stable temperature. Minimize exposure of cryopreserved iCell Mesenchymal Stem Cells to ambient temperature when transferring vials to liquid nitrogen storage.*

## Chapter 3. Preparing Media

Notes

**Note:** *iCell Mesenchymal Stem Cells will plate and function optimally on tissue culture treated vessels.*

Prepare and store Plating Medium and Maintenance Medium for iCell Mesenchymal Stem Cells as follows:

Plating Medium Component	Amount (ml)	Final Concentration
FBS	5	10%
IMDM	45	90%

- 1 Filter Plating Medium using a 0.2  $\mu$ m PES filter unit.
- 2 Store Plating Medium at 4°C for up to 2 weeks.

Maintenance Medium Component	Amount (ml)	Final Concentration
Ascorbic Acid, 20 mg/ml	2.5	50 $\mu$ g/ml
B-27 Supplement, Minus Vitamin A, 50X	10	1%
bFGF, 1 mg/ml	0.05	50 ng/ml
BSA, 10% in D-PBS	5	0.05%
GlutaMAX Supplement	10	1%
Ham's F-12 Medium	250	25%
IMDM	750	75%
MTG, 11.5 M	0.04	450 $\mu$ M
N-2 Supplement, 100X	5	0.5%
PDGF-BB, 100 $\mu$ g/ml	0.5	50 ng/ml
Penicillin/Streptomycin	10	1%

- 1 Prepare 10% BSA in D-PBS. Filter 10% BSA using a 0.2  $\mu$ m PES filter unit. Store remaining 10% BSA at 4°C for up to 1 month.
- 2 Filter Maintenance Medium using a 0.2  $\mu$ m PES filter unit.
- 3 Store Maintenance Medium at 4°C for up to 2 weeks.

## Chapter 4. Thawing iCell Mesenchymal Stem Cells

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

1. Equilibrate Plating Medium and Maintenance Medium at room temperature before thawing iCell Mesenchymal Stem Cells.
2. Remove the iCell Mesenchymal Stem Cells 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 and gently swirl (avoid submerging the cap) until the cell suspension is fully thawed. Use of a floating microcentrifuge tube rack is recommended.
4. When the cell suspension is fully thawed, immediately remove the cryovial from the water bath, spray with 70% ethanol, and place in a biological safety cabinet.
5. Gently transfer the iCell Mesenchymal Stem Cells cryovial contents to a sterile 15 ml centrifuge tube using a 1 ml pipettor.



*Avoid repeated pipetting of the thawed iCell Mesenchymal Stem Cells suspension.*

6. Rinse the empty iCell Mesenchymal Stem Cells cryovial with 1 ml of room temperature Plating Medium to recover residual cells from the cryovial. Transfer the 1 ml of Plating Medium rinse drop-wise to the 15 ml centrifuge tube containing the iCell Mesenchymal Stem Cells suspension. Gently swirl the tube while adding the medium to mix the solution completely and minimize the osmotic shock on the thawed cells.



*Drop-wise addition of Plating Medium to the cell suspension is critical to minimize osmotic shock and ensure maximum viability and subsequent attachment of the cells.*

7. Slowly add 8 ml of Plating Medium to the 15 ml centrifuge tube. Gently swirl the tube while adding the Plating Medium.



*It is critical to add the 8 ml of Plating Medium slowly to ensure maximum viability and attachment of the cells once plated.*

8. Gently mix the contents of the 15 ml centrifuge tube by pipetting the cell suspension once. Avoid repeated pipetting, vigorous shaking, or vortexing of the cell suspension.

9. Centrifuge the cell suspension at 400 x g for 5 minutes at room temperature.
10. Carefully aspirate the supernatant, taking care not to disturb the cell pellet.
11. Slowly add 5 ml of Maintenance Medium to resuspend the cell pellet.

**Note:** Thaw no more than 3 vials of iCell Mesenchymal Stem Cells at one time. Once thawed, combine the contents of the cryovials after adding the rinse and final volume of Plating Medium. For example, if pooling 3 cryovials, add each 1 ml of rinse and 8 ml of additional Plating Medium. Once thawed and diluted to the desired density, cells can be pooled together for plating.

Notes

## Chapter 5. Plating iCell Mesenchymal Stem Cells

The following procedure describes how to plate iCell Mesenchymal Stem Cells at 35,000 viable cells/cm<sup>2</sup> into a 6-well cell culture plate. Scale volumes appropriately for other cell culture vessel formats.

1. Invert the thawed iCell Mesenchymal Stem Cells suspension 2 - 3 times to ensure an even cell distribution before performing the cell count.
2. Remove a sample of cells to confirm the viability of the cells using a hemocytometer (using trypan blue exclusion to identify viable cells) or an automated cell counter.
3. Dilute the cell suspension in room temperature Maintenance Medium to 1.68 x 10<sup>5</sup> cells/ml).

Culture Vessel	Surface Area (cm <sup>2</sup> )	Maintenance Medium (ml)	Cell Number (3.5 x 10 <sup>4</sup> cells/cm <sup>2</sup> )
6-well Cell Culture Plate	9.6	2	3.36 x 10 <sup>5</sup>
10 cm Cell Culture Dish	57	10	1.995 x 10 <sup>6</sup>

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

*All volumes and measures are per well, if applicable.*

4. Invert the cell suspension 2 - 3 times. Immediately dispense 2 ml/well of cell suspension into the cell culture vessel (~336,000 viable cells/well).
5. Culture iCell Mesenchymal Stem Cells in a cell culture incubator at 37°C, 5% CO<sub>2</sub>.

## Chapter 6. Maintaining iCell Mesenchymal Stem Cells

iCell Mesenchymal Stem Cells are cryopreserved at high purity. The cells retain their high purity when maintained and cultured as recommended.

1. Equilibrate Plating Medium and Maintenance Medium at room temperature before use.
2. 48 hours post-plating iCell Mesenchymal Stem Cells, aspirate the spent medium and replace with the appropriate volume of fresh Maintenance Medium.
3. Replace the Maintenance Medium every other day.
4. Culture iCell Mesenchymal Stem Cells in a cell culture incubator at 37°C, 5% CO<sub>2</sub>.
5. Passage iCell Mesenchymal Stem Cells at 80% - 90% confluency. The following instructions apply to one well of a 6-well cell culture plate. Scale volumes appropriately for other vessel formats.

**Note:** *iCell Mesenchymal Stem Cells will plate and function optimally on tissue culture treated vessels.*

- a. Aspirate the spent medium from the 80% - 90% confluent cell culture.
- b. Add 1 ml/well of TrypLE.
- c. Incubate at 37°C until cells detach from the plate (approximately 5 - 7 minutes).
- d. Tap the vessel gently to dislodge cells from the plate. Gently pipette up and down once to dislodge cells. Transfer the cell suspension to a 15 ml centrifuge tube. Microscopically check the plate to confirm that cells have been removed.



*Do not remove cells from the plate by scraping. If cells are not harvested according to these instructions, there may be a loss of viability, poor recovery, or loss of cell surface markers.*

- e. Rinse each well with 2 ml of Plating Medium. Transfer the 2 ml rinse from the well to the 15 ml centrifuge tube containing the iCell Mesenchymal Stem Cells suspension to quench the TrypLE.
- f. Resuspend cells by gently pipetting. Remove a sample of cells to perform a cell count using a hemocytometer or an automated cell counter.
- g. Centrifuge the cell suspension at 400 x g for 5 minutes.
- h. Carefully aspirate the supernatant, taking care not to disturb the cell pellet.
- i. Resuspend cells in the appropriate volume of Maintenance Medium to plate iCell Mesenchymal Stem Cells at a density of 35,000 cells/cm<sup>2</sup> (i.e. 336,000 cells/well in 2 ml of Maintenance Medium for a 6-well cell culture plate). Suggested cell numbers for different vessel formats are provided in Table 1 on page 9.

## Notes

6. Add iCell Mesenchymal Stem Cells into the culture vessel.

**Note:** *iCell Mesenchymal Stem Cells can be maintained by routine splitting for up to 5 passages. To avoid a poor quality cell culture, do not allow cells to become over-confluent at any time. It is recommended to passage at 80% - 90% confluency.*

---

### 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 were made.

**C. DATA.** Customer agrees that if described on Customer's product quotation from Cellular Dynamics it will provide data and information as described therein to Cellular Dynamics regarding Customer's use of the Products.

---

### Appendix B. Product Provided "AS IS"

A. The Products are sold or provided "AS IS."

B. TO THE FULLEST EXTENT PERMITTED BY APPLICABLE LAW, CELLULAR DYNAMICS DISCLAIMS, ALL REPRESENTATIONS, AND WARRANTIES, EXPRESS OR IMPLIED (INCLUDING BUT NOT LIMITED TO, ANY IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE OR NON-INFRINGEMENT), AND ALL LIABILITY, WITH RESPECT TO THE PRODUCTS; AND BY ORDERING AND ACCEPTING THE PRODUCTS CUSTOMER WAIVES ALL RIGHTS AND REMEDIES OTHERWISE AVAILABLE WITH RESPECT TO THE PRODUCTS.

C. Customer will be solely responsible for (i) Customer's use of the Products for a purpose or in a manner other than that for which they were designed or that is permitted or in breach of the Use Restrictions above; (ii) Customer's failure to follow this User's Guide for the use,



## Notes

storage, and handling of the Products however such failure is caused; (iii) Customer's failure to comply with any of the provisions of Appendix A above; and (iv) any abuse, other misuse or neglect of the Products by Customer or 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).

D. 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. Because the number of cells in a unit 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.

---

### 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.

© 2016 Cellular Dynamics International, Inc. All rights reserved.





