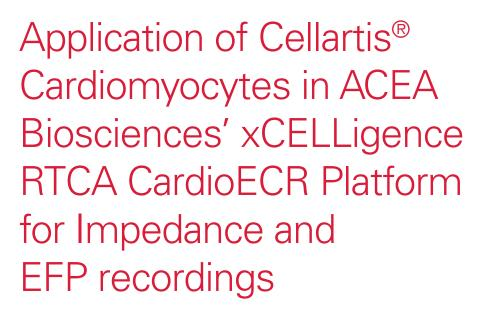


Stem Cell Application Protocol





I. Introduction

Cellartis Cardiomyocytes are derived from human induced pluripotent stem cells and provide a promising physiologically-relevant, human model for pre-clinical safety evaluation and drug screening. The hybrid system xCELLigence RTCA CardioECR allows for both impedance readout and extracellular field potential (EFP) recordings in high throughput format. Cellartis Cardiomyocytes used in combination with this technique form an excellent platform to accurately predict cardiotoxic responses and to screen compound efficacy.

II. Materials Required

- Cellartis Cardiomyocytes (from ChiPSC22) Kit (Takara Bio, Cat. No Y10075)
 - Cellartis Cardiomyocytes (from ChiPSC22)
 - Cellartis CM Thawing Base
 - Cellartis CM Culture Base
- Fetal Bovine Serum (FBS) (Life Technologies, Cat. No. 16140)
- Y-27632
- Electronic 48-well microtiter plate (E-Plate® CardioECR 48, ACEA Biosciences)
- Fibronectin (Sigma-Aldrich, Cat. No. F0895)
- PBS Dulbecco's with Ca²⁺ & Mg²⁺ (D-PBS +/+)
- xCELLigence RTCA CardioECR instrument (ACEA Biosciences)
- General cell culture equipment used in cell culture laboratory

III. Protocol

NOTE: Avoid contact with the electrodes in all of the following procedures as they are fragile. These procedures should be performed under aseptic conditions as much as possible.

A. Coating of the E-Plate® CardioECR 48



- 1. Dilute the required volume of Fibronectin in D-PBS +/+ to a final concentration of 10 µg/ml.
- 2. Add the diluted Fibronectin solution into each well to be used. Use 50 μl/well.
- 3. Incubate at 37°C for 3 hrs.
- 4. Aspirate the Fibronectin solution from the cell culture plate just before use.

B. Medium Preparation

Preparing Cellartis CM Thawing Medium

- Thaw Cellartis CM Thawing Base.
- 2. Decontaminate the external surface of all bottles with an appropriate disinfectant and place into the biological safety cabine.
- 3. Add 8 ml FBS per 32 ml Cellartis CM Thawing Base to achieve Cellartis CM Thawing Medium.
- 4. Cellartis CM Thawing Medium should be stored at 4°C and expires one month after the date of preparation.
- 5. Always discard any leftover warmed Cellartis CM Thawing Medium.

Preparing Cellartis CM Thawing Medium with Y-27632

- On the day of use, prepare Cellartis CM Thawing Medium with Y-27632 by adding Y-27632 to a final concentration of 10
 μM to Cellartis CM Thawing Medium.
- 2. Cellartis CM Thawing Medium with Y-27632 should be used on the day of preparation.

Preparing Cellartis CM Culture Medium

- Thaw Cellartis CM Culture Base.
- Decontaminate the external surface of supplement and medium bottle with appropriate disinfectant and place into the biological safety cabinet.
- 3. Add 10 ml FBS per 90 ml Cellartis CM Culture Base to achieve Cellartis CM Culture Medium.
- 4. Cellartis CM Culture Medium should be stored at 4°C and expires one month after the date of preparation.
- 5. Always discard any leftover warmed Cellartis CM Culture Medium.

Preparing Cellartis CM Culture Medium with Y-27632

- On the day of use, prepare Cellartis CM Culture Medium with Y-27632 by adding Y-27632 to a final concentration of 10 μM
 to Cellartis CM Culture Medium.
- 2. Cellartis CM Culture Medium with Y-27632 should be used on the day of preparation.

C. Thawing and Plating of Cellartis Cardiomyocytes

NOTE: It is recommended that not more than of two to three vials are thawed at once.

NOTE: For your protection, wear a protective face mask and protective gloves. Use forceps when handling a frozen vial. Never hold the vial in your hand as it may explode due to rapid temperature changes.

- 1. Prepare the appropriate volume of Cellartis CM Thawing Medium with Y-27632 (see Section B) and warm to room temperature (RT, 15–25°C).
- 2. Transfer, as guickly as possible, the frozen vial from liquid nitrogen to a 37°C ± 1°C water bath using forceps.
- 3. Thaw the cells by gently pushing the vial under the surface of the water, without swirling the vial. Do not submerge the cap of the vial in the water bath as this could contaminate the cells.
- 4. Take the vial out of the water bath as soon as the thawing is completed (approximately 3 min; the vial should still be cold on the outside).



- 5. Wipe the vial with an appropriate disinfectant and place into the biological safety cabinet.
- 6. As soon as possible, gently transfer the cell suspension into a sterile 50 ml tube by using a pipette.
- 7. Rinse the vial with 1 ml of Cellartis CM Thawing Medium with Y-27632 and carefully add it to the cell suspension dropwise.
- 8. Add 8 ml of Cellartis CM Thawing Medium with Y-27632 dropwise. Gently swirl the tube a few times in between.
- 9. Centrifuge the tube at 200*g* for 5 min at RT and remove the supernatant.
- 10. Carefully re-suspend the cell pellet with Cellartis CM Thawing Medium with Y-27632, using 4 ml medium per thawed vial.
- 11. Count the cells and measure viability.
- 12. Adjust the number of viable cells to 5 x 10⁵ cells/ml with Cellartis CM Thawing Medium with Y-27632.

NOTE: Aspirate the Fibronectin solution just before adding the cell suspension. Prepare one column at a time since drying of the surface might result in crystallization of the Fibronectin and subsequent damaging of the cells.

- 13. Aspirate the Fibronectin solution just before adding the cell suspension, making sure the wells do not run dry.
- 14. Carefully mix your cell suspension and pipet 50 µl into each well (corresponding to 2.5 x 10⁴ cells/well).
- 15. Proceed rapidly with the remaining columns.
- 16. Place the plate in the incubator (37°C \pm 1°C, 5% CO2, and >90% humidity).
- 17. After 3 hrs, carefully add 130 µl Cellartis CM Thawing Medium with Y-27632 to each well to reach a final volume of 180 µl.

D. Medium change

It is recommended to do the first medium change 48 hours after thawing and plating, and further on every 2–3 days (e.g. Monday, Wednesday, Friday) until analysis.

Medium preparation

1. Prepare the appropriate volume of Cellartis CM Culture Medium with 10 μ M Y-27632 as described in section B and warm to 37°C \pm 1°C before use.

Medium change

NOTE: Work very gently in order not to detach the cells.

- Replace the medium with 180 µl of fresh Cellartis CM Culture Medium with Y-27632.
- 2. Place the plate in the incubator (37°C \pm 1°C, 5% CO2, and >90% humidity).

NOTE: Impedance studies are best conducted after 7 days of post-thaw culture, depending on the stabilization of the signals.



	P R O D U C T S		
Cat.	#	Product	Size
Y100)75	Cellartis Cardiomyocytes (from ChiPSC22)	1 kit

Notice to Purchaser

Your use of these products and technologies is subject to compliance with any applicable licensing requirements described on the product's web page at http://www.clontech.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, in vitro diagnostic purposes, therapeutics, or in humans. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Takara Bio USA, Inc.

Partnering for Your Success

We are committed to developing and providing optimal tools that are tailored to your projects. To ensure your success, our Cellartis iPS Cell Services team delivers professional expertise through close collaboration.



Generating clinical-grade human pluripotent stem cell lines? We can help!

Derivation of proprietary, clinical-grade hES cell lines from blastocysts, or hiPS cell lines from blood samples, includes:

- Project report including all relevant documents, QA and QC package
- Seed Bank(s) of up to three customer-proprietary donor hPS cell lines
- Master Cell Bank of one customer-proprietary hPS cell line with a project-specific QC

Interested in learning more about how our services can support the goals of your project? Visit www.clontech.com/stem-cell-services



