

RevitaCell™ Supplement (100X)

Description

RevitaCell™ Supplement (100X) is a chemically defined recovery supplement containing a specific ROCK inhibitor coupled with molecules that have antioxidant and free radical scavenger properties. Use RevitaCell™ Supplement to achieve efficient and optimal post-thaw recovery of cryopreserved pluripotent stem cells (PSCs) as well as primary cells, including neurons, human epidermal keratinocytes (HEKs), and human corneal epithelial cells (HCECs). RevitaCell™ Supplement can also be used for the recovery of PSCs from single-cell passaging procedures using dissociation reagents such as TrypLE™ Select, TrypLE™ Express, and StemPro® Accutase®.

| Product* | Catalog no. | Amount | Storage | Shelf life* |
|-------------------------------|-------------|--------|-----------------------------------|-------------|
| RevitaCell™ Supplement (100X) | A2644501 | 5 mL | -20°C to -5°C; Protect from Light | 12 months |

* Shelf Life duration is determined from Date of Manufacture.

Product use

For Research Use Only. Not for use in diagnostic procedures.

Safety information

Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Important information

Divide thawed RevitaCell™ Supplement (100X) into usage-size aliquots and store in a non-frost-free freezer at -20°C to -5°C.

Recovery of cryopreserved PSCs

1. Coat the culture vessels with the appropriate substrate on which to culture your PSCs. Recommended Substrates include Vitronectin (VTN-N) Recombinant Human Protein (Cat. no. A14700) and Geltrex® LDEV-Free hESC-qualified Reduced Growth Factor Basement Membrane Matrix (Cat. no. A14133).
2. Quickly thaw cryopreserved PSCs in a 37°C waterbath until only a small ice crystal remains.
3. Gently pipet the thawed cells up and down to create a cell suspension and transfer the cell suspension into a 50-mL conical tube.
4. Dilute the cell suspension with 3 mL of growth medium, adding it dropwise while gently rocking the tube back and forth to avoid osmotic shock to the cells.
5. Centrifuge the cell suspension at 200 × g for 4 minutes.
6. Aspirate the medium, being careful not to disturb the cell pellet.
7. Gently resuspend the cells in growth medium supplemented with RevitaCell™ Supplement at a 1X final concentration (i.e., 100 µL of RevitaCell™ Supplement in 10 mL of growth medium).
Note: Do not add any additional ROCK inhibitors to the growth medium.
8. Transfer the cell suspension to the culture vessel pre-coated with the appropriate substrate. Move the vessel in several quick back-and-forth and side-to-side motions to disperse the cells across its surface.
9. Incubate the cells for 18–24 hours in the recommended cell culture environment.
10. Following incubation, aspirate the growth medium supplemented with RevitaCell™ Supplement and replace it with unsupplemented growth medium (i.e., without the addition of RevitaCell™ Supplement) for the remainder of the culture.

Recovery of cryopreserved primary cells

1. Quickly thaw cryopreserved primary cells in a 37°C waterbath until only a small ice crystal remains.
2. Recover the cells according to the recommended protocol for your cell-type. Perform cell count.
3. Prepare the cells for seeding by diluting the cell suspension in medium containing RevitaCell™ Supplement at a 1X final concentration (i.e., 100 µL of RevitaCell™ Supplement in 10 mL of growth medium).
Note: Do not add any additional ROCK inhibitors to the growth medium.
4. Transfer the cell suspension to the appropriate culture vessel and incubate for 18–24 hours in the recommended cell culture environment.
5. Following incubation, aspirate the growth medium supplemented with RevitaCell™ Supplement and replace it with unsupplemented growth medium (i.e., without the addition of RevitaCell™ Supplement) for the remainder of the culture.

Recovery of single-cell passaged PSCs in Essential 8® Medium

1. Coat the culture vessels with the appropriate substrate on which to culture your PSCs. Recommended Substrates include Vitronectin (VTN-N) Recombinant Human Protein (Cat. no. A14700) and Geltrex® LDEV-Free hESC-qualified Reduced Growth Factor Basement Membrane Matrix (Cat. no. A14133).
2. Pre-warm the required volume of TrypLE™ Select, TrypLE™ Express, or StemPro® Accutase® dissociation reagent in a 37°C waterbath.
3. Aspirate the spent medium from the culture vessel.
4. Rinse the vessel once with recommended volume of Dulbecco's Phosphate Buffered Saline (DPBS) without calcium or magnesium (see Table 1).
5. Add recommended volume of prewarmed TrypLE™ Select, TrypLE™ Express, or StemPro® Accutase® dissociation reagent to the vessel (see Table 1).
6. Incubate the vessel at 37 °C, 5% CO₂ for 5 minutes, continually observing the wells for cell detachment.
Note: Avoid extended incubation of cells with dissociation reagents to minimize damage to cells.
7. Gently pipet the cells up and down 5–10 times to generate single cell suspension.
8. Transfer the cell suspension to a conical tube containing the recommended volume of Essential 8™ Medium to dilute the dissociation reagent (see Table 1).
9. Centrifuge the cells at 200 × g for 4 minutes.

- Discard the supernatant, flick the tube 3–5 times to loosen the pellet, and then resuspend the cells by pipetting them up and down 5–10 times in the recommended volume of Essential 8[®] Medium supplemented with RevitaCell[™] Supplement at a 1X final concentration (see Table 1).
 - Determine viable cell density and percent viability using a Countess[®] Automated Cell Counter or a similar automated or manual method.
 - Adjust the concentration of the cell suspension using Essential 8[®] Medium supplemented with RevitaCell[™] Supplement at a 1X final concentration to achieve the cell seeding density recommended for your culture vessel (see Table 2).
- Note:** Do not add any additional ROCK inhibitors to the growth medium.
- Transfer the cell suspension to the culture vessel pre-coated with the appropriate substrate. Move the vessel in several quick back-and-forth and side-to-side motions to disperse the cells across its surface.
 - Incubate the cells for 18–24 hours in the recommended cell culture environment.
 - Following incubation, aspirate the Essential 8[®] Medium supplemented with RevitaCell[™] Supplement and replace it with unsupplemented Essential 8[®] Medium (i.e., without the addition of RevitaCell[™] Supplement) for the remainder of the culture.

Table 1 Reagent volumes (per well or per dish)

| Culture vessel (surface area) | DPBS for wash | Dissociation reagent* | Essential 8 [®] Medium** | Essential 8 [®] Medium + 1X RevitaCell [™] Supplement*** |
|-------------------------------|---------------|-----------------------|-----------------------------------|--|
| 6-well (10 cm ²) | 2 mL | 1 mL | 3 mL | 2 mL |
| 12-well (4 cm ²) | 1 mL | 0.4 mL | 1.2 mL | 1 mL |
| 24-well (2 cm ²) | 0.5 mL | 0.2 mL | 0.6 mL | 0.5 mL |
| 35-mm (10 cm ²) | 2 mL | 1 mL | 3 mL | 2 mL |
| 60-mm (20 cm ²) | 4 mL | 2 mL | 6 mL | 4 mL |
| 100-mm (60 cm ²) | 12 mL | 6 mL | 18 mL | 12 mL |

* TrypLE[™] Express, TrypLE[™] Select, or StemPro[®] Accutase[®] reagent

** For dilution *** For resuspension

Table 2 Recommended cell seeding densities and volumes of medium for plating (per well or per dish)

| Culture vessel (surface area) | Number of viable cells added* | | Essential 8 [®] Medium + 1X RevitaCell [™] Supplement** |
|-------------------------------|-------------------------------|------------------------------|---|
| | 12,500 cells/cm ² | 25,000 cells/cm ² | |
| 6-well (10 cm ²) | 125,000 | 250,000 | 2 mL |
| 12-well (4 cm ²) | 50,000 | 100,000 | 1 mL |
| 24-well (2 cm ²) | 25,000 | 50,000 | 0.5 mL |
| 35-mm (10 cm ²) | 125,000 | 250,000 | 2 mL |
| 60-mm (20 cm ²) | 250,000 | 500,000 | 4 mL |
| 100-mm (60 cm ²) | 750,000 | 1,500,000 | 12 mL |

* Time to confluency is 4–5 days for a seeding density of 12,500 cells/cm² and 3–4 days for a seeding density of 25,000 cells/cm²

** For resuspension

For additional technical information such as Safety Data Sheets (SDS), Certificates of Analysis, visit www.lifetechnologies.com/support. For further assistance, email techsupport@lifetech.com.

© 2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Essential 8 is a registered trademark of Cellular Dynamics International, Inc. Accutase is a registered trademark of Innovative Cell Technologies.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

www.lifetechnologies.com

Related Products

| Product | Cat. no. |
|--|-----------|
| Essential 8 [®] Medium | A15170 |
| Vitronectin, truncated human recombinant (VTN-N) | A14700 |
| Geltrex [®] LDEV-Free, hESC-Qualified, Reduced Growth Factor Basement Membrane Matrix | A14133 |
| UltraPure [™] 0.5 M EDTA, pH 8.0 | 15575 |
| StemPro [®] Accutase [®] Cell Dissociation Reagent | A11105 |
| TrypLE [™] Express, no phenol red | 12604 |
| TrypLE [™] Select, no phenol red | 12563 |
| DPBS, no calcium, no magnesium | 14190 |
| Primary Rat Cortex Neurons | A10840 |
| Neurobasal [®] Medium 1X, liquid | 21103 |
| B-27 [®] Supplement 50X, liquid | 17504 |
| Human Epidermal Keratinocytes, neonatal (HEKn) | C0015C |
| Human Epidermal Keratinocytes, adult (HEKa) | C0055C |
| Human Corneal Epithelial Cells (HCEC) | C0185C |
| EpiLife [®] Medium, with 60 µM calcium | MEPI500CA |
| Human Keratinocyte Growth Supplement (HKGS) 100X | S0015 |
| Human Corneal Growth Supplement (HCGS) 100X | S0095 |

Explanation of Symbols and Warnings

| | | | | |
|---|---|---|---|---|
|  |  |  |  |  |
| Caution, consult accompanying documents | Temperature Limitation | Keep away from light | Use By: | Consult instructions for use |
|  |  |  |  |  |
| Batch Code | Catalog number | Manufacturer | Sterilized using aseptic processing techniques | Read Safety Data Sheet |

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Important licensing information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

 life technologies