

# Growth Supplement (100X)

## Description

Growth Supplement (100X) is a chemically defined recovery supplement. Use Growth Supplement to achieve efficient and optimal post-thaw recovery of cryopreserved primary cells, including human epidermal keratinocytes (HEKn and HEKa), human corneal epithelial cells (HCEC), and neurons, as well as Pluripotent Stem Cells (PSCs). Additionally use Growth Supplement for recovery from single cell passaging of PSCs using TrypLE™ Select, TrypLE™ Express, and StemPro® Accutase® dissociation reagents.

Product	Catalog no.	Amount	Storage
Growth Supplement (100X)	A24422SA	5 mL	-20°C to -5°C; Protect from light

## Product use

For Research Use Only. Not for use in diagnostic procedures. Product is a prototype and performance characteristics of this product have not been established.

## Important information

Divide thawed Growth Supplement (100X) into usage-size aliquots and store at -20°C to -5°C.

## Safety information



**CAUTION!** Skin Absorption. Contains DMSO. Components of the product may be absorbed into the body through the skin. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing and gloves.

## Recovery of cryopreserved PSCs

1. Quickly thaw cryopreserved PSCs in a 37°C waterbath until a small ice crystal remains.
2. Gently resuspend the cells and transfer the cell suspension to a 50-mL conical tube.
3. Dilute the cryopreserved cell suspension with 3 mL of growth medium by adding growth medium dropwise while gently rocking the tube back and forth to avoid osmotic shock.
4. Centrifuge the cell suspension at 200 × g for 4 minutes.
5. Aspirate the medium, being careful not to disturb the cell pellet.
6. Gently resuspend the cells in growth medium supplemented with 1X Growth Supplement.
7. Transfer the cell suspension to the appropriate growth vessel and incubate for 18–24 hours in the recommended cell culture environment.
8. Following incubation, aspirate the growth medium supplemented with 1X Growth Supplement and replace it with unsupplemented growth medium (without the addition of Growth Supplement) for the remainder of culture.

## Recovery of cryopreserved primary cells

1. Quickly thaw cryopreserved primary cells in a 37°C waterbath until a small ice crystal remains.
2. Recover the cells according to cell-type specific protocols. Perform cell counts.
3. Prepare the cells for seeding by diluting the suspension in medium containing 1X Growth Supplement.
4. Transfer the cell suspension to the appropriate growth vessel and incubate for 18–24 hours in the recommended cell culture environment.
5. Following incubation, aspirate the growth medium supplemented with 1X Growth Supplement and replace it with unsupplemented growth medium (without the addition of Growth Supplement) for the remainder of culture.

## Recovery of single cell passaged PSCs










1. Perform single cell dissociation according to recommended protocols for TrypLE™ Select, TrypLE™ Express, or StemPro® Accutase™ dissociation reagents.
2. Prepare a cell suspension containing the desired number of cells per mL in growth medium containing 1X Growth Supplement.
3. Transfer the cell suspension to the appropriate growth vessel and incubate for 18–24 hours in the recommended cell culture environment.
4. Following incubation, aspirate the growth medium supplemented with 1X Growth Supplement and replace it with unsupplemented growth medium (without the addition of Growth Supplement) for the remainder of culture.

## 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 Enzyme 1X, no phenol red	12604
CTS™ TrypLE™ Select Enzyme	A12859
Primary Rat Cortex Neurons	A10840
Neurobasal® Medium 1X, liquid	21103
B-27® Supplement 50X, liquid	17504
Human Epidermal Keratinocytes, neonatal (HEKKn)	C0015C
Human Epidermal Keratinocytes, adult (HEKa)	C0055C
Human Corneal Epithelial Cells (HCEC)	C0185C
EpiLife® Medium	MEPI500CA
Human Keratinocyte Growth Supplement (HKGS) 100X	S0015
Human Corneal Growth Supplement (HCGS) 100X	S0095

## Explanation of symbols and warnings

The symbols present on the product label are explained below:

		
Caution, consult accompanying documents	Temperature Limitation	Keep away from light
		
Sterilized using aseptic processing techniques	Consult instructions for use	Batch Code
		
Catalog number	Manufacturer	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](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

## Important licensing information

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

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

© 2014 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or their affiliate(s) or their respective owners. Accutase is a registered trademark of Innovative Cell Technologies. B-27 is a registered trademark of Board of Trustees of Southern Illinois University. Essential 8 is a registered trademark of Cellular Dynamics International, Inc.

**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](http://www.lifetechnologies.com)

*life*  
technologies™