


# FreeStyle™ 293-F Cells

Catalog Number R79007

Pub. No. MAN0007834 Rev. C.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

## Product description

Gibco™ FreeStyle™ 293-F Cells are derived from the 293 cell line and are intended for use with the FreeStyle™ MAX 293 Expression System or FreeStyle™ 293 Expression System.

FreeStyle™ 293-F Cells can be thawed, grown, maintained, and transfected in FreeStyle™ 293 Expression System.

## Contents and storage

Contents	Cat. No.	Amount	Storage
FreeStyle™ 293-F Cells	R79007	1 vial (1 × 10 <sup>7</sup> cells)	Liquid nitrogen. Protect cultures from light.

## Required materials not supplied

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Item	Source
FreeStyle™ 293 Expression Medium	12338018
125-mL polycarbonate, disposable, sterile, vent-cap Erlenmeyer shaker flask or other appropriate vessel for culturing suspension cells	MLS
Orbital shaker in temperature and CO <sub>2</sub> controlled incubator	MLS
Reagents and equipment to determine cell viability (e.g., hemocytometer with trypan blue or cell counter)	MLS

## Procedural guidelines

- Subculture the FreeStyle™ 293-F Cells a minimum of three times to allow them to recover from thawing before using them in transfection experiments.
- Keep cell densities between  $1 \times 10^6$ – $3 \times 10^6$  cells/mL of culture for best performance.
- We recommend maintaining cells in a 125-mL or a 250-mL polycarbonate, disposable, sterile Erlenmeyer flask containing 25–40 mL or 50–80 mL total working volume of cell suspension, respectively.
- Glass flasks without baffles may be used, but clean them thoroughly after each use to avoid potential toxicity.

## Guidelines for thawing and storing cells

- On receipt, either thaw the cells immediately or immediately place the frozen cells into vapor phase liquid nitrogen storage until ready to use. Do not store the cells at  $-80^\circ\text{C}$ .
- Avoid short-term, extreme temperature changes. When storing cells in liquid nitrogen following receipt on dry ice, allow the cells to remain in liquid nitrogen for 3–4 days before thaw.
- Before starting experiments, ensure to have cells that are established and have frozen stocks on hand. On receipt, grow and freeze multiple vials of cells to ensure that you have an adequate supply of early-passage cells.

## FreeStyle™ 293-F Cells characteristics

**Growth properties:** Suspension

**Doubling time:** 25 hours. Doubling times may vary based on cell health, handling, and passage number.

**Viability during log phase culture:** >90%

**Subculture conditions:** Grow to  $1 \times 10^6$ – $3 \times 10^6$  cells/mL, and split cells to  $0.2 \times 10^6$ – $0.5 \times 10^6$  cells/mL every 2–3 days. Do not grow above  $3 \times 10^6$  cells/mL for best performance. Discard cells when they reach passage number 30.

## Scale up FreeStyle™ 293-F cell culture


You can scale up the FreeStyle™ 293-F cultures in spinner flasks or bioreactors. Determine the optimal spinner or impeller speed and seeding density for your culture system. We recommend that the cells be seeded at  $0.5 \times 10^6$  viable cells/mL. Optimum spinner speed is approximately 100–130 rpm, and optimum impeller speed in Celligen™ stirred tank bioreactors is 70–100 rpm. If the split ratio of cells to fresh media is less than 1:2, centrifuge the cell suspension and re-suspend the cell pellet in fresh medium before inoculating the culture.

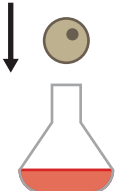
## Cryopreserve FreeStyle™ 293-F Cells

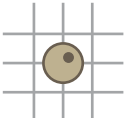
FreeStyle™ 293-F Cells can be frozen directly in FreeStyle™ 293 Expression Medium.


- Freeze FreeStyle™ 293-F Cells at a final density of  $1 \times 10^7$  viable cells/mL.
- Use a freezing medium composed of 90% fresh FreeStyle™ 293 Expression Medium and 10% DMSO.
- Freeze cells in an automated or manual, controlled-rate, freezing apparatus following standard procedures. For ideal cryopreservation, the rate of temperature decrease should be  $1^\circ\text{C}$  per minute.
- Transfer frozen vials to liquid nitrogen for storage.
- Check the viability and recovery of frozen cells 24 hours after storing cryovials in liquid nitrogen by following the procedures outlined in this protocol.


# Thaw and passage FreeStyle™ 293-F Cells in FreeStyle™ 293 Expression Medium

- 1 Day 1: Thaw cells**  


Rapidly thaw the cells in a water bath, decontaminate the vial using 70% ethanol, and open the cryovial in a class II biological cabinet.
- 2 Day 1: Add cells to medium**  


Add cells to 29 mL of pre-warmed medium in a 125-mL shake flask.
- 3 Day 1: Count cells and determine viability**  


Within 1–2 hours post-thaw, count cells and determine viability. Use hemocytometer and trypan blue exclusion method or automated cell counter. Cell density should be approximately  $0.3 \times 10^6$  cells/mL and cell viability >90%.
- 4 Day 1: Incubate**  


**Temperature:** 37°C  
**Humidified atmosphere:** 8% CO<sub>2</sub> in air  
**Orbital shaker platform:** 125 rpm
- 5 Day 3–4: Subculture cells**  


**First passage:** When cell density reaches  $>1 \times 10^6$  cells/mL at  $\geq 90\%$  viability (typically 2–3 days post-thaw), split the culture to  $0.2 \times 10^6$ – $0.5 \times 10^6$  cells/mL in FreeStyle™ 293 Expression Medium.  
**Subsequent passages:** Every 2–3 days, cells should reach  $1 \times 10^6$ – $3 \times 10^6$  cells/mL. Split to  $0.2 \times 10^6$ – $0.5 \times 10^6$  cells/mL. Do not grow above  $3 \times 10^6$  cells/mL. We recommend using a 125-mL or a 250-mL flask containing 25–40 mL or 50–80 mL of medium, respectively. When cell density reaches  $>1 \times 10^6$  cells/mL at  $\geq 90\%$  viability (typically 2–3 days post-thaw), split the culture to  $0.2 \times 10^6$ – $0.5 \times 10^6$  cells/mL in FreeStyle™ 293 Expression Medium.

## 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 at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, CA 92008

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**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.

©2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.