


ExpiSf9™ Cells

Catalog Number A35243

Pub. No. MAN0017552 Rev. B.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.

Product description

The Gibco™ ExpiSf9™ cell line is a derivative of the Sf9 (*Spodoptera frugiperda*) cell line. ExpiSf9™ cells are adapted to high-density, serum-free suspension culture in ExpiSf™ CD Medium. The cells can be thawed directly into ExpiSf™ CD Medium. Transfection, baculovirus production, and protein expression experiments can be performed in ExpiSf™ CD Medium.

Contents and storage

Contents	Cat. No.	Amount	Storage
ExpiSf9™ Cells (1×10^7 cells/mL)	A35243	1.5 mL	Liquid nitrogen, vapor phase

Required materials not supplied

Unless otherwise indicated, all materials are available through thermofisher.com.

Item	Source
ExpiSf™ CD Medium	A3767802
Nalgene™ Single-Use PETG Erlenmeyer Flasks with Plain Bottom: Sterile	4115-0125

Procedural guidelines

- Store the frozen ExpiSf9™ cells in vapor phase liquid nitrogen until ready to use.

IMPORTANT! Do not store the cells at -80°C .

Culture conditions

Media: ExpiSf™ CD Medium

Cell line: ExpiSf9™ cells

Culture type: Suspension

Shake flask type: It is recommended to use PETG, non-baffled, vented Erlenmeyer flasks; however, baffled Erlenmeyer flasks may also be used.

Temperature range: $27.5^{\circ}\text{C} \pm 0.5^{\circ}\text{C}$

Shaker speed: For shakers with a 19-mm or 25-mm shaking diameter, set the shake speed to 125 ± 5 rpm. For shakers with a 50-mm shaking diameter, set the shake speed to 95 ± 5 rpm.

Incubator type: Non-humidified, air regulated, non-CO₂ atmosphere. Ensure proper gas exchange and minimize exposure of culture to light.

Guidelines for ExpiSf9™ cell culture

- ExpiSf9™ is a robust cell line adapted to high density growth in ExpiSf™ CD Medium with a doubling time of approximately 24 hours during log-phase growth.
- The cells have a broad log-phase growth window spanning approximately from 4×10^6 – 12×10^6 cells/mL with a maximum density of approximately 20×10^6 cells/mL in standard shake flask culture.
- For general cell maintenance, passage ExpiSf9™ cells when they reach a density of 5×10^6 – 10×10^6 viable cells/mL (i.e., early to mid-log phase growth), typically every 3–4 days.

Note: Cells that are subcultured at densities outside of this log-phase growth window may show longer doubling times and lower baculovirus and/or protein titers over time. If necessary, modify the initial seeding density to attain the target cell density of 5×10^6 – 10×10^6 viable cells/mL at the time of subculturing.
- Use a hemocytometer with the trypan blue exclusion method or an automated cell counter to determine cell viability. Log-phase cultures should be $\geq 90\%$.
- When thawing or subculturing cells, transfer cells into pre-warmed (i.e., room temperature) ExpiSf™ CD Medium.
- We recommend thawing a fresh low-passage vial of cells every 2.5–3 months or 20–30 passages.

Thaw ExpiSf9™ cells

1. Remove one vial of cells from liquid nitrogen and swirl in a 37°C water bath for 1 to 2 minutes to thaw the cells rapidly until only a small amount of ice remains.
Note: Do not submerge the vial in the water.
2. Just before the cells are completely thawed, decontaminate the vial by wiping with 70% ethanol before opening it in a laminar flow hood.
3. Transfer the entire contents of the cryovial with a 2-mL or 5-mL pipette into a 125-mL PETG, sterile, non-baffled, vented shake flask containing 25 mL of pre-warmed (i.e., room temperature) ExpiSf9™ CD Medium.

IMPORTANT! ExpiSf9™ CD Medium comes in a ready-to-use format. For suspension growth, baculovirus production, and protein expression applications, use the ExpiSf9™ CD Medium without any supplementation.

4. Incubate the cells at 27.5°C ±0.5°C in a non-humidified, non-CO₂ atmosphere incubator on an orbital shaker platform set at 125 ±5 rpm (for shakers with a 19-mm or 25-mm shaking diameter) or 95 ±5 rpm (for shakers with a 50-mm shaking diameter).
5. Three days post-thaw, determine viable cell density and percent viability.
Note: Cell viability should be ≥80% by three days post-thaw.
6. Continue to monitor cell density and viability and subculture the cells once the culture has reached 5 × 10⁶ – 10 × 10⁶ viable cells/mL (typically 4–5 days post-thaw).

Guidelines to subculture ExpiSf9™ cells

- Passage ExpiSf9™ cells directly in ExpiSf9™ CD Medium.
- It is recommended to use a 125-mL or 250-mL PETG, sterile, non-baffled, vented shake flask containing 20–32% total working volume of cell suspension. When using larger flasks, the total working volume should be between 25–33%.
- Subculture ExpiSf9™ cells when they attain a density of 5 × 10⁶–10 × 10⁶ viable cells/mL.
- Cells should exhibit only minimal clumping during routine cell culture maintenance.

Subculture ExpiSf9™ cells

1. Using the viable cell density, calculate the volume of cell suspension required to seed a new shake flask according to the recommended seeding densities in Table 1 and the recommended culture volumes in Table 2.
2. Transfer the calculated volume of cells to fresh, pre-warmed (i.e., room temperature) ExpiSf9™ CD Medium in a shake flask.

3. Incubate flasks at 27.5°C ±0.5°C in a non-humidified, non-CO₂ atmosphere incubator on a shaker platform set at 125 ±5 rpm (for shakers with a 19-mm or 25-mm shaking diameter) or 95 ±5 rpm (for shakers with a 50-mm shaking diameter) until cultures reach a density of 5 × 10⁶–10 × 10⁶ viable cells/mL.

Note: If necessary, modify the initial seeding density to attain the target cell density of 5 × 10⁶–10 × 10⁶ viable cells/mL at the time of subculturing.

4. Repeat Steps 1–3 to maintain or expand the cells for baculovirus production or protein expression.

Table 1 Recommended seeding densities for routine cell culture

Subculture timing	Seeding density
Cells ready 3 days post-passaging	0.7 × 10 ⁶ –1.0 × 10 ⁶ viable cells/mL
Cells ready 4 days post-passaging	0.4 × 10 ⁶ –0.6 × 10 ⁶ viable cells/mL

Table 2 Recommended culture volume using different non-baffled shake flask sizes

Shake flask size	Recommended culture volume
125-mL	25–30 mL
250-mL	50–60 mL
500-mL	100–120 mL
1-L	200–240 mL
2-L	400–480 mL
3-L	600–800 mL

Guidelines to cryopreserve ExpiSf9™ cells

- ExpiSf9™ cells can be frozen directly in ExpiSf9™ CD Medium.
- Freeze ExpiSf9™ cells at a final density of 1 × 10⁷ viable cells/mL in 1.5 mL total volume of 92.5% conditioned ExpiSf9™ CD Medium and 7.5% DMSO.
- Allow cells to attain a viable cell density of 3 × 10⁶–4.5 × 10⁶ cells/mL and ≥95% viability before harvest.

Note: For cryopreservation, the viable cell density at time of harvest is critical for optimal cell health. Therefore, make sure to only harvest cells when they are within the recommended 3 × 10⁶–4.5 × 10⁶ viable cells/mL range. If viable cell density is too low at the time of harvest, return cells to the incubator until the cells reach the recommended density. If viable cell density is too high at the time of harvest, subculture the cells at 0.5 × 10⁶–0.6 × 10⁶ cells/mL and prepare for harvest again after 3–4 days.

Cryopreserve ExpiSf9™ cells

1. Centrifuge the cells at 300 × g for 5 minutes.
2. Decant the spent “conditioned” medium into a sterile conical tube or bottle.

IMPORTANT! Do not discard the supernatant.

3. Resuspend the cell pellet using the appropriate volume of conditioned medium collected above to achieve a final cell density of 1×10^7 viable cells/mL, and gently resuspend the cell pellet by pipetting up/down.
4. Add the required volume of DMSO (7.5% final) to the cell suspension and gently mix.
5. Immediately aliquot 1.5 mL cell suspension volume per cryovial.
6. Freeze the cells in an automated or manual controlled-rate freezing apparatus following standard procedures.
Note: For ideal cryopreservation, the freezing rate should be a decrease of 1°C per minute.
7. Transfer frozen vials to liquid nitrogen (vapor phase) for long-term storage.

5. Repeat Step 4, increasing stepwise the ratio of ExpiSf™ CD Medium to original medium (75:25 followed by 90:10) until the cells are transferred into 100% ExpiSf™ CD Medium.

Note: Multiple passages at each step may be needed. Subculture cells when viable cell density is $\geq 5 \times 10^6$ cells/mL at each step.

6. Once fully adapted to ExpiSf™ CD Medium, the viable cell density should exceed 5×10^6 cells/mL with a viability $\geq 90\%$ within 3–4 days of subculture (when using a seeding density of 1×10^6 cells/mL).

At this stage the seeding density may be reduced to 0.5×10^6 – 1.0×10^6 viable cells/mL for subsequent passaging.

7. At this point, it is recommended to create a cell bank of the adapted cells. As with any cell line adaptation, it is best practice to continue to culture and monitor the cells as they gain more passages in ExpiSf™ CD Medium and generate additional cell banks, as necessary.

Guidelines to adapt Sf9 and Sf21 cells to ExpiSf™ CD Medium

- You may adapt your Sf9 and Sf21 cells for growth in ExpiSf™ CD Medium.
- We recommend using a sequential adaptation protocol.
- It is critical that cell viability be at least 90% and the growth rate be in mid-logarithmic phase prior to initiating adaptation procedures.

Note: This procedure is meant to serve as a guideline for your adaptation process. Different growth kinetics and maximum cell density may be observed during the adaptation of your Sf9 or Sf21 cell line.

Adapt Sf9 and Sf21 cells to ExpiSf™ CD Medium

1. Subculture Sf9 or Sf21 cells into a 25:75 ratio of ExpiSf™ CD Medium to the original media.
Note: In the event that cells lag during this first transition, it is possible to reduce the ratio to 10:90 as a first step and/or to use conditioned media for this first step.
Note: During the adaptation procedure use a seeding density of 1.0×10^6 viable cells/mL at each sub-culturing step.
2. Incubate at $27.5^\circ\text{C} \pm 0.5^\circ\text{C}$ in a non-humidified, non- CO_2 atmosphere incubator on an orbital shaker platform set at 125 ± 5 rpm (19-mm or 25-mm shaking diameter) or 95 ± 5 rpm (50-mm shaking diameter).
3. Three to four days after subculture, remove a small amount of the cell suspension and perform a cell count to determine viable cell density and viability.
Note: If cells are $>90\%$ viable and $\geq 5 \times 10^6$ viable cells/mL, they are ready to be passaged. If cells are at $\geq 90\%$ viability, but have not reached the desired density, return cells to the incubator for 1–2 days, until the viable cell density is $\geq 5 \times 10^6$ cells/mL.
4. Subculture when the viable cell density is $\geq 5 \times 10^6$ cells/mL by passaging cells into a 50:50 ratio of ExpiSf™ CD Medium to original medium.

Related products

Unless otherwise indicated, all materials are available through thermofisher.com.

Item	Source
ExpiSf™ Expression System Starter Kit	A38841 ^[1] A39112 ^[2] A39111 ^[3]
ExpiFectamine™ Sf Transfection Reagent	A38915
ExpiSf™ Protein Production Kit, 1 L	A3767806
ExpiSf™ Protein Production Kit, 10 L	A3767807
ExpiSf™ Protein Production Kit, 5 × 10 L	A3767808
Bac-to-Bac™ Baculovirus Expression System	10359-016
Bac-to-Bac™ C-His TOPO™ Expression System	A11100
Bac-to-Bac™ N-His TOPO™ Expression System	A11101
Bac-to-Bac™ HBM TOPO™ Secreted Expression System	A11339
Nalgene™ Single-Use PETG Erlenmeyer Flasks with Plain Bottom: Sterile	4115-0125
Trypan Blue Stain	15250

^[1] North America, Europe

^[2] Latin America, Asia Pacific, Japan

^[3] China

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. If you have any questions, please contact Life Technologies at 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.

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.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.