ExpiFectamine[™] Sf Transfection Reagent

Catalog Number A38915

Pub. No. MAN0017502 **Rev.** A.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

ExpiFectamineTM Sf Transfection Reagent is a proprietary, cationic lipid formulation suitable for transfecting DNA into insect cells. This reagent is a component of the ExpiSfTM Expression System and has been optimized to transfect ExpiSf9TM cells using the Bac-to-BacTM Baculovirus Expression System.

Contents and storage

Product	Cat. No. Contents		Storage	
ExpiFectamine [™] Sf Transfection Reagent	A38915	1 mL	Store at 4°C. Do not freeze	

Required materials not supplied

Item	Source		
ExpiSf9™ Cells	A35243		
ExpiSf™ CD Medium	A3767801-A3767805		
Opti-MEM [™] I Reduced Serum Medium	31985062		
Bacmid DNA prepared using the Bac-to- Bac™ Baculovirus Expression System	10360014, A11098, A11099, A11338		
Nalgene [™] Single-Use PETG Erlenmeyer Flasks with Plain Bottom: Sterile	4115-0125		
Thomson Instrument Company 24-well deep-well plate	NC0012954		
Non-humidified, air-regulated, non-CO ₂ atmosphere incubator set at 27°C	TFS ^[1]		
Orbital shaker platform (for suspension- based transfection protocol)			
<i>Optional:</i> Sf9 or Sf21 cells			
<i>Optional:</i> Sf-900 [™] II SFM or Sf-900 [™] III SFM			
<i>Optional:</i> 6-well tissue culture-treated plate			

^[1] Available from Thermo Fisher Scientific.

Procedural guidelines

- We recommend isolating bacmid DNA using the PureLink HiPure Plasmid Prep Kits (Cat. No. K2100). To ensure sterility, you may filter your DNA preparation through a 0.22-µm filter before use.
- Allow freshly thawed cells to recover in culture for two or more passages post-thaw before transfecting.
- Gently invert the ExpiFectamine[™] Sf Transfection Reagent 5-10 times before use to ensure thorough mixing.
- ExpiFectamine[™] Sf/DNA complexes must be made in serumfree medium. We recommend using Opti-MEM[™] I Reduced Serum Medium for the complexation reaction.
- Do **not** add antibiotics to media during transfection because it may decrease transfection efficiency.
- ExpiFectamine[™] Sf Transfection Reagent can be used to transfect ExpiSf9[™], Sf9, and Sf21 cells in either adherent or suspension culture format. Two different transfection protocols are included in this manual for added flexibility.
- Make sure you have a healthy cell culture at log phase growth and ≥90% viability prior to proceeding with transfection.



Approximate experimental step times

Procedure	Time
Cell preparation	0-60 minutes
Incubation of diluted lipid reagent	5 minutes
Incubation of DNA-lipid mixture	5 minutes
Cell incubation post-transfection	72-96 hours

Suspension-based transfection of ExpiSf9[™], Sf9, and Sf21 cells with ExpiFectamine[™] Sf Transfection Reagent using the Bac-to-Bac[™] Baculovirus Expression System

Use the following procedure to transfect ExpiSf9[™], Sf9, and Sf21 suspension cell cultures in a 125-mL shake flask format. All amounts are given on a per-flask basis.

1	Prepare cells and medium	a.	At the time of transfection, dilute cells to 2.5×10^6 cells/mL ($\geq 90\%$ viability) in 25 mL of growth medium (ExpiSf TM CD Medium in a 125-mL non-baffled, vented shake flask.) Prepare the cells as follows.		
			Note: The following instructions are given on a per-shake flask basis.		
			1. Pipet 62.5×10^6 viable cells into a sterile 50-mL conical tube.		
			2. Centrifuge at $300 \times g$ for 5 minutes.		
			3. Aspirate the supernatant and gently resuspend cells in 25 mL of fresh growth medium.		
			4. Transfer the entire cell suspension to a 125-mL shake flask.		
		b.	Incubate cells for 0-30 minutes in a 27°C non-humidified, non-CO ₂ 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).		
			Note: It is important to proceed to Step 2 within 30 minutes of cell seeding. Incubating cells for longer than 30 minutes prior to transfection may result in decreased transfection efficiency.		
2	Dilute ExpiFectamine [™] Sf Transfection Reagent in Opti-MEM [™] I Medium and incubate	a.	Gently mix the ExpiFectamine [™] Sf Transfection Reagent before use by inverting 5-10 times.		
		b.	Dilute 30 µL ExpiFectamine [™] Sf Transfection Reagent in 1 mL Opti-MEM [™] I Reduced Serum Medium.		
		c.	Gently mix by inverting 5-10 times.		
		d.	Incubate the mixture at room temperature for 5 minutes.		
3	Add Bacmid DNA directly to diluted ExpiFectamine [™]	a.	Add 12.5 μg Bacmid DNA directly to the diluted ExpiFectamine [™] Sf Transfection Reagent. No pre-dilution of Bacmid DNA is required.		
	St Iranstoction Poagont				
	Sf Transfection Reagent and incubate	b.	Gently mix by inverting 5-10 times.		
			Gently mix by inverting 5-10 times. Incubate the mixture at room temperature for 5 minutes.		
4		c.			
4	and incubate Add DNA-lipid complex to	c. a.	Incubate the mixture at room temperature for 5 minutes. Slowly transfer the mixture dropwise to the prepared 125-mL shake flask from Step 1 swirling		

Adherent-based transfection of ExpiSf9[™], Sf9, and Sf21 cells with ExpiFectamine[™] Sf Transfection Reagent using the Bac-to-Bac[™] Baculovirus Expression System

Use the following procedure to transfect ExpiSf9[™], Sf9, and Sf21 adherent cell cultures in a 6-well plate format. All amounts are given on a per-well basis.

1	Prepare cells and medium	a.	At the time of transfection, seed cells at 1 × 10 ⁶ cells/well (≥90% viability) in complete growth medium (ExpiSf [™] CD Medium in a 6-well plate (3 mL total volume per well)).	
		b.	Allow cells to attach for 30-60 minutes in a 27°C non-humidified, non-CO ₂ incubator.	
2	Dilute ExpiFectamine [™] Sf	a.	Gently mix the ExpiFectamine [™] Sf Transfection Reagent before use by inverting 5-10 times.	
Transfection Reagent in Opti-MEM [™] I Medium and incubate		b.	Dilute 10 µL ExpiFectamine [™] Sf Transfection Reagent in 250 µL Opti-MEM [™] I Reduced Serum Medium.	
		c.	Gently mix by inverting 5-10 times.	
		d.	Incubate the mixture at room temperature for 5 minutes.	
3	Add Bacmid DNA to diluted ExpiFectamine [™] Sf	a.	Add 1 µg Bacmid DNA directly to the diluted ExpiFectamine [™] Sf Transfection Reagent. No pre- dilution of Bacmid DNA is required.	
Transfection Reagent and incubate		b.	Gently mix by inverting 5-10 times.	
		c.	Incubate the mixture at room temperature for 5 minutes.	
4	Add DNA-lipid complex to cells	a.	Slowly transfer the mixture from Step 3 dropwise to the appropriate well of the prepared 6-well plate from Step 1.	
		b.	Incubate cells in a 27°C non-humidified, non-CO ₂ incubator until you see signs of viral infection (typically 72-96 hours).	
			Note: ExpiFectamine [™] Sf Transfection Reagent exhibits low cytotoxicity; therefore, no media change is required post-transfection.	

Scaling-up or -down transfections

Refer to the table below for the volumes of ExpiFectamine[™] Sf and other reagents required to transfect suspension insect cultures in different formats.

Note: The volumes given are on a per-well or per-shake flask basis.

Denemeter	Culture Type Suspension					
Parameter						
Vessel type	Deep-well plate	Vented, non-baffled shaker flask				
Vessel size	24 DWP	125 mL	250 mL	500 mL		
Number of cells required	10 × 10 ⁶ cells	62.5 × 10 ⁶ cells	125 × 10 ⁶ cells	250 × 10 ⁶ cells		
Culture volume to transfect	4 mL/well	25 mL	50 mL	100 mL		
	250 ± 5 rpm (19-mm shaking diameter)	125 ± 5 rpm (19-mm shaking diameter)				
Shake speed ^[1]		125 ± 5 rpm (25-mm shaking diameter)				
		95 ± 5 rpm (50-mm shaking diameter)				
Amount of plasmid DNA	1 µg	12.5 µg	25 µg	50 µg		
ExpiFectamine [™] Sf Transfection Reagent volume	5 µL	30 µL	60 µL	120 µL		
Opti-MEM [™] I Reduced Serum Medium volume	250 μL	1 mL	2 mL	4 mL		
Bacmid DNA volume ^[2]	2-4 µL	25-50 μL	50-100 μL	100-200 µL		

^[1] Optimal shake speed should be determined empirically based on the specific laboratory equipment used.

 $^{[2]}$ Bacmid DNA volume provided assumes a stock concentration of 250-500 ng/µL

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

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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: 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.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

