

# Growth and Maintenance of BHK Cells

Catalog no. R700-01

Version B

001127  
28-0062



**U.S. Headquarters:**

Invitrogen Corporation  
1600 Faraday Avenue  
Carlsbad, CA 92008  
Toll Free Tel: (800) 955-6288  
Tel: (760) 603-7200  
Fax: (760) 602-6500  
E-mail: [tech\\_service@invitrogen.com](mailto:tech_service@invitrogen.com)  
Web: [www.invitrogen.com](http://www.invitrogen.com)

**European Headquarters:**

Invitrogen BV  
PO Box 2312, 9704 CH Groningen  
The Netherlands  
Toll Free Tel: 00800 5345 5345  
Toll Free Fax: 00800 7890 7890  
Tel: +31 (0) 50 5299 299  
Fax: +31 (0) 50 5299 281  
E-mail: [tech\\_service@invitrogen.nl](mailto:tech_service@invitrogen.nl)



# Culturing BHK Cells

---

## Shipping/Storage

Cells are shipped on dry ice. Store in liquid nitrogen upon receipt.

---

## Contents

One vial contains  $3 \times 10^6$  baby hamster kidney (BHK) cells in 1 ml of  $\alpha$ MEM, 10% fetal bovine serum, 10% DMSO.

---

## BHK Cells

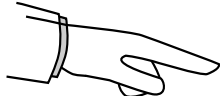
The BHK cell line was derived from baby Syrian hamster (*Mesocricetus auratus*) kidney (Macpherson and Stoker, 1962).

- The medium for BHK cells is  $\alpha$ MEM or DMEM (BioWhittaker).
  - Complete medium for BHK cells is  $\alpha$ MEM + 2 mM L-glutamine + **5% fetal bovine serum**. FBS does not need to be heat inactivated for use with BHK cells. **Note: 10% FBS makes the cells grow too fast. Infection with Sindbis viral particles will result in premature cell lysis, leading to low yields of protein expression.**
  - Cells are grown in a humidified, 37°C, 5% CO<sub>2</sub> incubator.
  - If cells are split at 1:5, it will take 1-2 days to reach 80-90% confluency.
- 

## Product Qualification

BHK cells must be 85% viable when recovered in  $\alpha$ MEM medium. The cells are thoroughly tested for the absence of mycoplasma contamination.

---



## Note

The cells have a tendency to clump in complete medium. In general, this is not a problem except when preparing the cells for electroporation. In this case, care must be taken to avoid clumps. PBS is required to keep the cells from clumping during electroporation.

---

## General Cell Handling

Use the procedures below to initiate and maintain a culture of BHK cells.

- **All solutions and equipment that come in contact with the cells must be sterile.**
  - Always use proper sterile technique and work in a laminar flow hood.
  - Use cells that are 80-90% confluent and > 90% viability for transfections and infections.
  - Before starting experiments, be sure to have cells established and also have some frozen stocks on hand.
  - For general maintenance of cells, pass BHK cells when they are 80-90% confluent (1-2 days) and split at a 1:5 dilution. For example, transfer 2 ml of a 10 ml cell suspension (without trypsin/EDTA) to a new 75 cm<sup>2</sup> flask and add 10 ml fresh medium.
  - Use trypan blue exclusion to determine cell viability. Thawed cells should be 80 to 85% viable and healthy, log phase cultures should be > 90% viable.
  - Cells may be passaged 60-70 times before re-starting a culture from frozen stocks.
- 

*continued on next page*

# Culturing BHK Cells, continued

---

## Before Starting

Be sure to have the following solutions and supplies available:

- 15 ml sterile, conical tubes
  - 5, 10, and 25 ml sterile pipettes
  - Cryovials
  - PBS (page 5)
  - 0.4% Trypan blue in PBS
  - Hemacytometer
  - $\alpha$ MEM medium (BioWhittaker)
  - Tissue culture grade 200 mM L-glutamine
  - FBS
  - Complete  $\alpha$ MEM medium ( $\alpha$ MEM + 2 mM L-glutamine + 5% FBS, page 5)
  - Freezing Medium ( $\alpha$ MEM medium + 2 mM L-glutamine + 10% FBS + 10% DMSO, page 5)
  - Table-top centrifuge
  - 75 cm<sup>2</sup> flasks, 175 cm<sup>2</sup> flasks and 35 mm plates (other flasks and plates may be used)
  - Trypsin/versene (EDTA) solution (BioWhittaker) or other trypsin solution
- 

## Initiating Cell Culture from Frozen Stock

The following protocol is designed to help you initiate a cell culture from a frozen stock. Note that the vial of BHK cells contains  $3 \times 10^6$  cells.

1. Remove the vial of cells from the liquid nitrogen and thaw quickly at 37°C.
  2. Just before the cells are completely thawed, decontaminate the outside of the vial with 70% ethanol, and transfer the cells to a 15 ml sterile, conical tube.
  3. Add 9 ml of prewarmed (37°C), complete  $\alpha$ MEM medium dropwise to cells.
  4. Centrifuge in a table-top centrifuge at 250 x g for 5 minutes at room temperature. Decant the medium. (This removes the DMSO from the cells.)
  5. Resuspend the cells in 10 ml of complete  $\alpha$ MEM and test a small portion of the cell suspension for viability by trypan blue dye exclusion. Viability of the cells should be between 80 and 85%.
  6. Transfer the remaining cell suspension to a 75 cm<sup>2</sup> flask, and incubate at 37°C. Incubation for 1-2 days should yield an 80-90% confluent monolayer.
- 

*continued on next page*

# Culturing BHK Cells, continued

---

## Passaging the BHK Cells

1. When cells are ~80-90% confluent, remove all medium from the flask.
  2. Wash cells once with 10 ml PBS to remove medium. Serum contains inhibitors of trypsin.
  3. Add 5 ml of trypsin/versene (EDTA) solution to the monolayer and incubate 1 to 5 minutes at room temperature until cells detach. Check the cells under a microscope and confirm that most of the cells have detached. If cells are still attached, incubate a little longer until most of the cells have detached.
  4. Once the cells have detached, briefly pipet the solution up and down to break up clumps of cells.
  5. Add 5 ml of complete  $\alpha$ MEM to stop trypsinization. Centrifuge cells at 250 x g for 5 minutes.
  6. Aspirate the supernatant and resuspend the cells in 10 ml complete  $\alpha$ MEM. Check the viability of the cells. Cells should be > 90% viable.
  7. To maintain cells in 75 cm<sup>2</sup> flasks, transfer 2 ml of the 10 ml cell suspension from Step 6 to a new 75 cm<sup>2</sup> flask and add 10 ml fresh, complete  $\alpha$ MEM medium.
  8. To expand cells, transfer 3 to 4 ml of the cell suspension to a 175 cm<sup>2</sup> flask (3 flasks total) and add fresh, complete  $\alpha$ MEM medium to a final volume of 30 ml.
  9. Incubate flasks in a humidified, 37°C, 5% CO<sub>2</sub> incubator.
- Repeat Steps 1-9 as necessary to maintain or expand cells.
- 

## Freezing the BHK Cells

Before starting, label cryovials and prepare freezing medium (page 5).

1. When cells are ~80% confluent in a 175 cm<sup>2</sup> flask, remove the medium and wash the cells one time with 10 ml PBS.
  2. Add 5 ml of trypsin/versene (EDTA) solution and incubate 1 to 5 minutes until cells detach. Once cells have detached, briefly pipet solution up and down to break up clumps of cells.
  3. Add 5 ml of complete  $\alpha$ MEM to stop trypsinization. Count the cells in a hemacytometer.
  4. Pellet cells at 250 x g for 5 minutes in a table top centrifuge at +4°C and decant the medium.
  5. Resuspend the cells at a density of 3 x 10<sup>6</sup> cells/ml in freezing medium (see page 5).
  6. Aliquot 1 ml of the cell suspension per vial. Place vials at -20°C for 2-3 hours.
  7. Transfer vials to a -70 or -80°C freezer and hold overnight.
  8. Transfer vials to liquid nitrogen for long term storage.
- 

*continued on next page*

## Culturing BHK Cells, continued

---

### Complete $\alpha$ MEM Medium

To 500 ml  $\alpha$ MEM (BioWhittaker), add 5% FBS and 5 ml of 200 mM L-glutamine solution (BioWhittaker).

---

### Freezing Medium

$\alpha$ MEM containing 2 mM L-glutamine, 10% FBS, and 10% DMSO.

---

### Phosphate Buffered Saline

For washing cells only. The solution does not need to be RNase-free.

137 mM NaCl  
2.7 mM KCl  
10 mM Na<sub>2</sub>HPO<sub>4</sub>  
1.8 mM KH<sub>2</sub>PO<sub>4</sub>

1. Dissolve:
  - 8 g NaCl
  - 0.2 g KCl
  - 1.44 g Na<sub>2</sub>HPO<sub>4</sub>
  - 0.24 g KH<sub>2</sub>PO<sub>4</sub>

in 800 ml deionized water.

2. Adjust pH to 7.4 with concentrated HCl.
  3. Bring the volume to 1 liter and autoclave for 20 minutes on liquid cycle.
  4. Store at +4°C or room temperature.
- 

### Reference

Macpherson, I. A. and Stoker, M. G. P. (1962) *Virology*, **16**: 147.

---

## Technical Service

---

### World Wide Web



Visit the [Invitrogen Web Resource](#) using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- Subscribe to InvitroWire™, our individually-tailored electronic news service
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Post a question at one of our many user forums
- Request catalog and product literature

Once connected to the Internet, launch your web browser (Netscape 3.0 or newer), then enter the following location (or URL):

**<http://www.invitrogen.com>**

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

---

### Technical Service

For Technical Service, please call, write, fax, or E-mail:

#### **U.S. Headquarters:**

Invitrogen Corporation  
1600 Faraday Avenue  
Carlsbad, CA 92008  
Toll Free Tel: (800) 955-6288  
Tel: (760) 603-7200  
Fax: (760) 602-6500  
E-mail: [tech\\_service@invitrogen.com](mailto:tech_service@invitrogen.com)  
Web site: [www.invitrogen.com](http://www.invitrogen.com)

#### **European Headquarters:**

Invitrogen BV  
PO Box 2312, 9704 CH Groningen  
The Netherlands  
Toll Free Tel: 00800 5345 5345  
Toll Free Fax: 00800 7890 7890  
Tel: +31 (0) 50 5299 299  
Fax: +31 (0) 50 5299 281  
E-mail: [tech\\_service@invitrogen.nl](mailto:tech_service@invitrogen.nl)

---

### Limited Warranty

Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, please contact our Technical Service Representatives at:

1-800-955-6288, extension 2 (U.S. and Canada)

00800 5345 5345 (toll free) or +31 (0) 50 5299 299 (Europe, Middle East, and Africa)

1-760-603-7200 extension 2 (all others)

Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives.

**Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.**

---

# MSDS Information

---

## Introduction

In an effort to provide our customers with the best possible information any time they need it, Material Data Safety Sheets (MSDSs) are now available 24 hours a day, 7 days a week from either our web site or the 3E Company (see below). Please note that MSDSs will no longer be included with product shipments.

---

## MSDS Information Online

Chemicals that require MSDS information are listed on our web site:

<http://www.invitrogen.com/tech/index.html>

Select the "Safety & MSDS" link in the right-hand column. The table lists the compound name, CAS number, product catalog number, and MSDS part number. Click the compound name to view the MSDS information.

---

## MSDS Information by Phone

Customers of Invitrogen can call the 3E Company, 24 hours a day, 7 days a week for MSDS information. This information can be obtained directly over the phone, faxed, or mailed to the customer. Use the compound name, product catalog number, or compound part number listed in the upper right-hand corner of the label to obtain MSDS information.

3E Company

1905 Aston Avenue

Carlsbad, CA 92008

Voice: 1-800-451-8346 (U.S., Canada, and Guam)

Voice: 1-760-602-8700 (See below for other toll-free numbers)

Fax: 1-760-602-8888

---

## Emergency Information

In the event of an emergency, the 3E Company can help with disposal or spill information. They can also connect the customer with poison control or the University of California at San Diego Medical Center doctors.

---

## Toll-Free Numbers

If you reside in one of the countries listed below, you can contact the 3E Company using the toll-free number listed.

Country	Toll-Free Number	Country	Toll-Free Number
Belgium	008071178	Mexico	958008342735
China	108001100017		958004518346
France	0800903046		958003603220
Germany	0130829154		958003466737
Indonesia	00180316570620	Netherlands	08000220091
Italy	167870990	Singapore	8001100987
Japan	006633800185	Spain	900931263
Korea	0308116570820	Turkey	00800136570900
Malaysia	1800808714	United Kingdom	0800967491

---

©1995-2000 Invitrogen Corporation. Reproduction forbidden without permission.