

Grace's Antheraea Insect Media

Catalog No. B823-01, B823-02

Information for Use

Shipping/Storage

Grace's Antheraea Insect Media and the supplements are shipped at room temperature. Upon receipt, store the bottled medium and the supplements at +4°C in the dark. Medium and supplements can be stored up to 6 months in this manner prior to preparation for use (described below).

Contents

Each bottle of Grace's Antheraea Insect Media contains 500 ml of sterile-filtered medium and is accompanied by a pouch of supplements containing yeastolate, lactalbumin and L-glutamine.

Product Specifications

Grace's Antheraea Insect Medium is a clear liquid, free of particulate matter with the following characteristics:

- pH of 6.2 +/- 0.2
- Osmolality of 350-365 mOSM/kg
- Endotoxin levels are less than 1 ng/ml.

Grace's Insect supplements are a non-sterile powder containing 1.66 g of TC yeastolate, 1.66 g lactalbumin hydrolysate, and 300 mg of L-glutamine. Medium formulation is available from Technical Services (see next page).

Quality Control

To assess the performance of each lot of Grace's Antheraea Insect Medium, medium is prepared as described below and used to test the growth and viability of Sf9 cells. A seeding density of 0.5×10^6 cells/ml is carried through three passages of suspension culture in the medium. In the third passage, cells are tested for doubling time and viability. Doubling time must be between 18 and 24 hours and viability must be greater than 99%.

continued on next page

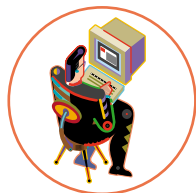
Information for Use, continued

Preparation of Medium for Use

To prepare a complete medium for insect cell growth, the provided supplements and fetal bovine serum (FBS) must be added to Grace's Antheraea Insect Media. With supplements and serum, the medium is called Complete TNM-FH (*Trichoplusia ni* Medium-Formulation Hink). The following procedure should be used to make Complete TNM-FH.

- 1) Dissolve the contents of one supplements pouch into one 500 ml bottle of Grace's Antheraea Insect Medium.
 - 2) Add 55 ml of tissue culture grade FBS and 500 µl of a 10 mg/ml stock of Gentamycin to the mixture of Grace's Medium and supplements. Final concentration of the FBS is 10% and final concentration of Gentamycin is 10 µg/ml.
 - 3) Filter the resulting solution through a 0.2 micron filter into a sterile container.
 - 4) Reconstituted medium should be stored at +4° C in the dark and is stable for one month.
-

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) 603-7201
E-mail: tech_service@invitrogen.com
Web: 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
