

High Five™ Serum-Free Medium

Catalog No. B100-01

Information for Use

Shipping/Storage

High Five™ Serum-Free Medium is shipped at room temperature. Upon receipt, store bottled medium at +4°C in the dark. Medium can be stored up to 6 months in this manner prior to preparation for use.

Contents

Each bottle of High Five™ Serum-Free Medium contains 1 liter of sterile, complete insect medium.

Product Specifications

High Five™ Serum-Free Medium is a sterile-filtered clear liquid that contains 1.5 g/L L-glutamine, 24 µg/L Tween-80, and 910 µg/L of Pluronic Poly-all.

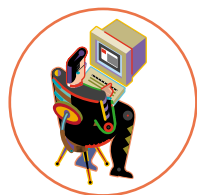
Quality Control

To assess the performance of each lot of High Five™ Serum-Free Medium, medium is prepared as described below and used to test the growth and viability of High Five™ cells. A seeding density of 1.0-2.0 x 10⁶ cells/ml is carried through three passages of suspension culture in the media. In the third passage, cells are tested for viability. Viability must be greater than 98%.

Preparation of Medium for Use

High Five™ Medium is a complete serum-free medium for growth of Sf9 cells, Sf21 cells, or High Five™ cells and requires no additional supplements for use. We recommend the addition of Gentamycin to 10 µg/ml prior to use.

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