

DES[®] Serum-Free Medium

Catalog No. Q510-01

Information for Use

Shipping/Storage

DES[®] Serum-Free Medium and L-glutamine are shipped at room temperature. Upon receipt, store the bottled medium and the L-glutamine at +4° C in the dark. Medium can be stored up to 6 months in this manner prior to preparation for use as described below.

Contents

Each bottle of DES[®] Serum-Free Medium contains 1 liter of medium and is accompanied by a vial containing 10 ml of 200 mM sterile L-glutamine.

Product Specifications

Each bottle of DES[®] Serum-Free Medium contains 1 liter of sterile-filtered clear media. L-glutamine is supplied as a 200 mM sterile filtered solution.

Quality Control

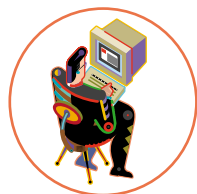
To assess the performance of DES[®] Serum-Free Medium, medium is prepared as described below and used to test the growth and viability of *Drosophila S2* cells. A seeding density of $3-8 \times 10^6$ cells/ml is carried through three passages of suspension culture in the medium. In the third passage, cells are tested for doubling time and viability. Doubling time must be between 18 and 24 hours and viability must be greater than 98%.

Preparation of Medium for Use

To prepare a complete medium for *Drosophila S2* growth, L-glutamine must be added to DES[®] Serum-Free Medium.

- 1) Pipet the 10 ml of L-glutamine provided into the DES[®] Serum-Free Medium and mix well. Final concentration of glutamine is approximately 2 mM.
 - 2) To ensure sterility, medium should be filter-sterilized with a 0.2 micron filter into a sterile container.
 - 3) After addition of L-glutamine, the medium is stable for 1 month if stored at +4 ° C in the dark.
-

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
