Gentamicin/Amphotericin Solution

Gentamicin & Amphotericin B

Catalog Number: R-015-10 10 x 1 ml vials per package

Product Description

Gentamicin/Amphotericin Solution is a sterile, concentrated (500x) antibiotic/antimycotic solution. When added to medium, the final mixture will contain 10 μ g/ml gentamicin and 0.25 μ g/ml amphotericin B.

Intended Use

Gentamicin/Amphotericin Solution solution may be used as an antibiotic/antimycotic solution in any of Cascade Biologics' cell culture media.

This product is for research use only. Not for use in animals, humans, or diagnostic procedures.

Caution: If handled improperly, some components of this product may present a health hazard. Take appropriate precautions when handling this product, including the wearing of protective clothing and eyewear. Dispose of properly.

Storage and Stability

Gentamicin/Amphotericin Solution is stored at $-20\,^{\circ}$ C at our facility and is shipped on dry ice. Upon receipt, Gentamicin/Amphotericin Solution should be stored at $-20\,^{\circ}$ C in a freezer that is <u>not</u> self-defrosting. When stored at $-20\,^{\circ}$ C, the product is stable until the expiration date shown on the label.

Preparation of Gentamicin/Amphotericin-Supplemented Medium

- 1. Obtain one bottle of Cascade Biologics' medium (500 ml) and one vial of Gentamicin/Amphotericin Solution.
- 2. Wipe the outside of the containers with a disinfecting solution such as 70% ethanol or isopropanol.
- 3. Using sterile technique in a laminar flow culture hood, transfer the entire contents of the vial of Gentamicin/Amphotericin Solution to the bottle of medium.
- 4. Tightly cap the bottle of supplemented medium and swirl to ensure a homogeneous solution. Avoid causing the medium to foam.

Terms and Conditions of Sale

Cascade Biologics, Inc. (hereinafter, CBI) warrants that its products will perform according to the information provided in various publications that it distributes, and as described herein for the intended shelf life of the product when stored under the conditions prescribed by CBI. If any product does not perform according to the published information provided by us, CBI will replace the product free of charge to the original customer. The remedy of product replacement shall be the customers' sole and exclusive remedy for defective product, unless CBI is unable to deliver replacement product, in which case CBI shall reimburse the customer for the purchase price of the defective product. Customer understands that the foregoing limited warranty is in lieu of all other warranties and CBI hereby disclaims all other warranties including, but not limited to, implied warranty of merchantability or fitness for adequacy for any particular purpose or use. CBI shall not be liable to customer or to any party claiming through customer for any incidental or consequential damages, including, but not limited to any lost profits, lost savings, or lost business, whether arising out of contract, tort, or otherwise. Customer acknowledges that products are intended for research use only and are not to be used in animals, humans, or diagnostic procedures.