### Instructions for Use

## NutriCoat™ Attachment Solution

A xeno-free solution that facilitates the attachment and spreading of human mesenchymal stem cells cultured in serum-free medium



2740220-000





A xeno-free solution that facilitates the attachment and spreading of human mesenchymal stem cells cultured in serum-free medium

REF

05-760-1-15 (1.5 mL/vial)



15-25°C (Room temperature)

# Contents

| 1 | Product Description                              | . 4 |
|---|--|-----|
| 2 | Precaution and Disclaimer                        | . 4 |
| 3 | Features   | . 4 |
|   | Instructions for Use                             |     |
| 5 | Quality Control                                  | . 6 |
|   | Quality Assurance.     6.1 Product Label Symbols |     |
| 7 | Related Products                                 | . 7 |

# 1 Product Description

A clear, colourless to yellowish, sterile solution, based on human Fibrinogen. NutriCoat™ Attachment Solution was qualified for the clinical grade of mesenchymal stem cells from various sources when culturing in MSC NutriStem® XF Medium.

## 2 Precaution and Disclaimer

- Not for human use; for research or further manufacturing use only.
- Do not use solution that is cloudy or has deposits.
- Do not use beyond the expiration date indicated on the product label.
- In order to maintain the sterility of the product, it is necessary to use a strict aseptic technique and use sterile materials and tools only.

## 3 Features

- Xeno-free
- Sterile by membrane filtration
- Easy-to-use
- Supplied at room temperature as a ready-to-use stock solution
- Design for use with serum-free, xeno-free culture system
- Enables cells dissociation using recombinant trypsin solutions (e.g. 03-078-1)

## 4 Instructions for Use

### 4.1 Pre-coating Procedure

- 1. Dilute the solution 1:500 in saline and gently mix using a pipette.
- 2. Coat the culture dish with the 1:500 solution. The volume should be adequate for covering the surface of the cultureware. Use table 1 for the recommended volumes (e.g. per 1 well of 12 w/p add 2 µL of NutriCoat™ in 1 mL saline).
- 3. Gently agitate and incubate for 1 hour in a humidified CO₂ incubator (37°C).

**NOTE** Longer incubation of coated culture dish (up to 24 hours) will not affect cell attachment and spreading.

4. Wash step is not required.

NOTE One wash using Dulbecco's PBS (without Ca & Mg) is an option.

5. Prior to cells seeding, carefully discard the solution and add a culture medium.

NOTE It is critical that the coating does not dry out.

Table 1: Recommended volume for the coating procedure

| Cultureware | Surface area cm² | Volume of the 1:500 diluted solution |
|-------------|------------------|--------------------------------------|
| 96-well     | 0.34             | 0.1 mL                               |
| 24-well     | 1.9              | 0.5 mL                               |
| 12-well     | 3.9              | 1 mL                                 |
| 6-well      | 9.6              | 2.5 mL                               |
| T25 Flask   | 25               | 6.5 mL                               |
| T75 Flask   | 75               | 19 mL                                |

# 5 Quality Control

Each lot is tested for sterility as well as its biological performance (the ability to promote the attachment and spreading of cells in serum-free medium).

# 6 Quality Assurance

- For research use or for use as ancillary material in manufacturing cell, gene or tissue-based products.
- Composed of cGMP medical grade raw materials.
- Manufactured under ISO 13485 QMS and in compliance with applicable cGMP guidelines.
- Manufactured under controlled environments and processes in accordance with:
  - 1. ISO 13408 Aseptic Processing of Health Care Products;
  - 2. ISO 14644 Airborne Particulate Cleanliness Classes in Clean Rooms and Clean Zones.

### Manufacturer



Biological Industries Israel Beit Haemek Ltd. Kibbutz Beit Haemek 2511500, Israel

### 6.1 Product Label Symbols

| REF         | Indicates the manufacturer's catalogue number so that the product can be identified. |
|-------------|--|
| LOT         | Indicates the manufacturer's batch code so that the batch or lot can be identified.  |
|             | <b>NOTE</b> Synonyms for batch code are lot number and batch number.                 |
| $\subseteq$ | Indicates the date after which the product is not to be used.                        |
| 1           | Indicates the temperature limits to which the product can be safely exposed.         |
| STERILE A   | Indicates a product that has been manufactured using accepted aseptic techniques.    |
| $\bigcap$ i | Indicates the need for the user to consult the instructions for use.                 |

# 7 Related Products

| Product   | Cat. No.             |
|---|----------------------|
| MSC NutriStem® XF Basal Medium and Supplement Mix | 05-200-1<br>05-201-1 |
| Recombinant Trypsin Solution                      | 03-078-1             |
| Recombinant Trypsin EDTA Solution                 | 03-079-1             |
| NutriFreez® D10 Cryopreservation Medium           | 05-713-1             |
| Dulbecco's PBS (without Ca & Mg)                  | 02-023-1             |
| Saline – Sodium Chloride 0.9% Solution            | 03-043-1             |

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen, Germany

Phone: +49 551 308 0 www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote the other gender as well.

### Copyright notice:

This instruction manual, including all of its components, is protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated: 02 | 2021

Printed in the EU on paper bleached without chlorine. | MB DIR: 2740220-000-00 Revision 01