

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



2740220-000-01



SARTORIUS

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	Features.....	4
3	Intended Use and Safety.....	4
4	Storage and Stability.....	5
5	Instructions for Use	5
6	Quality Control	6
7	Quality Assurance.....	6
8	Related Products	7

1 Product Description

A clear, colorless 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 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)

3 Intended Use and Safety

- For research use or for use as ancillary material in manufacturing cell, gene or tissue-based products
- Not intended for in vitro diagnostics use or use as a medical device
- Not intended for human in vivo applications
- Do not use if visible particles and | or precipitate are observed.
- Do not use beyond the expiration date indicated on the product label.
- Maintain aseptic work conditions.
- Do not use if there is any package leakage or exposure to environmental conditions, as the sterility of the product might be compromised.
- Refer to the Material Safety Data Sheet (MSDS) for hazard information.

4 Storage and Stability

- Store at 15°C to 30°C.
- Protect from direct light.
- Shelf life: Refer to product label for expiration date.

5 Instructions for Use

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 μ L 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 cm ²	Volume of the 1:500 diluted solution
96-well	0.34	0.1 mL
24-well	1.9	0.5 mL

Cultureware	Surface area cm ²	Volume of the 1:500 diluted solution
12-well	3.9	1 mL
6-well	9.6	2.5 mL
T25 Flask	25	6.5 mL
T75 Flask	75	19 mL

6 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). For full specifications, please check the lot specific Certificate of Analysis (CoA).

7 Quality Assurance







- Composed of cGMP medical grade raw materials
- Manufactured under ISO 13485 QMS and ISO 9001 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 – Cleanrooms and associated controlled environments

Manufacturer



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

Product Label Symbols

	Indicates the manufacturer's catalogue number so that the product can be identified.
	Indicates the manufacturer's batch code so that the batch or lot can be identified. NOTE Synonyms for batch code are lot number and batch number.
	Indicates the date after which the product is not to be used.
	Indicates the temperature limits to which the product can be safely exposed.
	Indicates a product that has been manufactured using accepted aseptic techniques.
	Indicates the need for the user to consult the instructions for use.

8 Related Products

Product	Cat. No.
MSC NutriStem® XF Basal Medium and Supplement Mix	05-200-1 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

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 all genders.

Copyright notice:

These instructions, including all components, are 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.

Last updated:

03 | 2024

© 2024
Biological Industries Israel Beit Haemek Ltd.
2511500 Kibbutz Beit Haemek
Israel

LH | Publication No.: SCM6036-e240302
DIR No.: 2740220-000-01