

Instructions for Use

MSC Attachment Solution

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



2740201-000



SARTORIUS

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

REF

05-752-1



2-8°C

Contents

1	Product Description	4
2	Features.....	4
3	Precaution and Disclaimer	4
4	Storage and Stability	4
5	Instructions for use.....	5
6	Quality Control	6
7	Quality Assurance.....	6
	7.1 Product Label Symbols	7
8	Auxiliary Products.....	8

1 Product Description

MSC Attachment Solution is a clear sterile solution containing human fibronectin (hFN) obtained by affinity purification on gelatine-sepharose from human plasma. This solution is particularly useful for the culture of cells that are not capable of synthesizing their own biomatrix or when culturing cells in serum-free medium. The solution was qualified for the attachment of human mesenchymal stem cells (hMSC) when culturing in serum-free media.

2 Features

- Xeno-free
- Easy-to-use
- Supplied as a ready-to-use stock solution
- Design for use with serum-free culture system

3 Precaution and Disclaimer

1. Do **not** use if a visible precipitate is observed in the MSC Attachment Solution.
2. Do **not** use beyond the expiration date indicated on the product label.

4 Storage and Stability

Stable at 2–8°C until the expiry date stated on the label.

5 Instructions for use

1. Dilute the MSC Attachment Solution 1:100 using sterile DPBS (without Ca⁺⁺ and Mg⁺⁺ (Cat. No. 02-023-1) and gently mix using a pipette.
2. Add the diluted MSC Attachment Solution to the cultureware. Gently agitate the coated cultureware and verify complete covering of the surface. Use Table 1 for recommended volumes.
3. Incubate the coated cultureware for at least 30 minutes in a humidified CO₂ incubator (37°C).
4. Following 30 minutes incubation:

For immediate use:

- Gently wash the cultureware with DPBS (For T-25 use at least 5 mL).
- Seed cells immediately.

It is critical that the coating does not dry out.

For later use:

- Wrap the coated cultureware with Parafilm® and incubate at 2–8°C. Coated cultureware stored under sterile conditions at 2–8°C are stable for 1 week.
- Gently wash the cultureware with DPBS.
- Seed cells immediately.

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 1:100 MSC Attachment Solution
96-well	0.34	0.1 mL
24-well	1.9	0.4 mL
12-well	3.9	0.8 mL
6-well/35 mm plate	9.6	2 mL
6 cm/T25 flask	25	5 mL
T75 flask	75	15 mL

6 Quality Control

Each lot is tested for sterility as well as for biological performance (the ability to promote the attachment and spreading of MSC in serum-free medium). For full specifications please check the lot specific Certificate of Analysis (CoA).

7 Quality Assurance

- For in vitro diagnostic use, research use or for use as ancillary material in manufacturing cell, gene or tissue based products.
- Listed in Europe under CE IVD class I, thus comply with European In-Vitro Diagnostic Devices Directive (98/79/EC) requirements.
- 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 – Cleanrooms and associated controlled environments



Authorized Representative in the European Community:

MedNet GmbH

Borkstraße 10, 48163 Münster, Germany



Manufacturer

Biological Industries Israel Beit Haemek Ltd.

Kibbutz Beit Haemek 2511500, Israel

7.1 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.

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.

STERILE A

Indicates a product that has been manufactured using accepted aseptic techniques.



Indicates that the product meets the requirements of the applicable EC directives

IVD

Indicates a product that is intended to be used as an in vitro diagnostic medical device.



Indicates the need for the user to consult the instructions for use.

8 Auxiliary Products

Product	Cat. No.
MSC NutriStem [®] XF Basal Medium	05-200-1
MSC NutriStem [®] XF Supplement Mix	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 (w/o 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 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: 2740201-000-00
Revision 01