

Instructions for Use

NutriStem[®] hPSC XF

A defined, Xeno-Free (XF), Serum-Free (SF) Media, Designed to Support the Growth of Human Embryonic Stem Cells (hESC) and Induced Pluripotent Stem Cells (iPSC)



2740227-000-02



SARTORIUS

A defined, Xeno-Free (XF), Serum-Free (SF) Media, Designed to Support the Growth of Human Embryonic Stem Cells (hESC) and Induced Pluripotent Stem Cells (iPSC)

NutriStem® hPSC XF with HSA
Xeno-free medium for feeder-free and feeder-dependent culture of hPSC and hiPSC

REF 05-100-1

 -20 to -10°C

NutriStem® hPSC XF Medium (Growth Factor-Free)

REF 06-5100-01

 -20 to -10°C

Contents

1	Product Description	4
2	Precaution and Disclaimer	4
3	Storage and Stability	5
4	Features.....	5
5	Instructions for use.....	6
6	Protocol for maintenance of undifferentiated hESC and hiPSC.....	6
7	Quality Control	14
8	Quality Assurance.....	14
8.1	Product Label Symbols	15
9	Related Products	16

1 Product Description

Traditional human Embryonic Stem Cells (hESC) and human induced Pluripotent Stem Cells (hiPSC) culture methods require the use of mouse or human fibroblast feeder layers, or feeder- conditioned medium. These culture methods are labor- intensive, hard to scale and it is difficult to maintain hPSC pluripotency due to undefined conditions.

NutriStem® hPSC XF media were developed with a leading group in stem cell research, to enable the maintenance and expansion of hESC and hiPSC in feeder-free culture or with feeder cells.

NutriStem® hPSC XF support the culture of undifferentiated hESC and hiPSC in serum-free conditions without any xeno-derived components on feeders such as Mouse Embryo Fibroblasts (MEF) and Human Foreskin Fibroblasts (HFF) as well as with Matrigel™ and recombinant laminin. The media contain recombinant human basic fibroblast growth factor (rhbFGF) and recombinant human transforming growth factor (rhTGFβ). The media have been successfully tested and proven to maintain the pluripotent nature of hESC and hiPSC.

For long-term growth of hESC and hiPSC without feeder cells it is recommended to use NutriStem® hPSC XF with HSA (Cat.# 05-100-1).

2 Precaution and Disclaimer

1. Do **not** use if a visible precipitate is observed in the medium.
2. Do **not** use beyond the expiration date indicated on the product label.
3. Do **not** use if liquid is present in the cap of the bottle.

3 Storage and Stability

- NutriStem® hPSC XF should be stored at -20 to -10°C. Upon thawing at room-temperature or at 2–8°C, the medium may be stored at 2–8°C for 2 weeks. In case smaller quantities are needed, use aseptic techniques to dispense into aliquots so repeated freezing and thawing are avoided.
- NutriStem® hPSC XF must be warmed to room temperature (15°C–25°C) before use. To ensure stability of the medium, warm only the amount needed.
- Protect the medium from light.
- Shelf Life: Refer to product label for expiration date.

4 Features

- Serum Free Media (SFM), xeno-free (XF): all components are defined and xeno-free including proteins.
- Sterile, 0.1 µm membrane filtered, under full aseptic conditions and practice.
- The proteins used: HSA (Human Serum Albumin), rh bFGF, rh TGFβ, human transferrin and recombinant human insulin.
- A complete, ready-to-use formulation (no additions are required).
- Contain Alanine glutamine. Do not contain antibiotics.
- Enable culture of hESC and hiPSC without feeder cells.
- Can be used with mouse or human feeder cells, Matrigel™, laminin and with matrix from human cells.
- Can be used with laminin 521 (e.g. LaminStem™ 521, Cat. No. 05-753-1) using single cell passage.
- Intended for use with 5% CO₂ (ordinary conditions).
- Support long-term growth of hESC and hiPSC, and maintain their ability to differentiate, without any signs of karyotype abnormalities.

5 Instructions for use

- For complete instructions on how to maintain hESC and hiPSC in NutriStem® hPSC XF, see technical manual guide.
- Specific pluripotent cell culture protocol may require optimization for best results. The following protocol is a generic guideline.
- Use aseptic techniques.

6 Protocol for maintenance of undifferentiated hESC and hiPSC

1. Feeder-dependent culture

Gelatin coating

Coat cultureware using 0.1% Gelatin solution according to the following table:

Plate dish	Volume of gelatin per well
4/24 wells	0.5 mL
6 wells	2 mL
35 mm	2 mL
10 cm ²	10 mL

Leave at room temperature or in incubator for at least two hours.

NOTE It is highly recommended to prepare gelatin-coated plates 24 hours before use.

Preparation of feeder cells (Mitomycin C treated)

1. MEF feeder cells should be passaged into T75 flasks twice and used on the third passage.
2. Discard the medium and add 8 µg/mL mitomycin C (6–7 mL) into a culture flask and incubate for two hours.
3. Wash three times with D-PBS without calcium and magnesium (Cat. No. 02-023-1).
4. Add 3 mL of trypsin-EDTA solution and cover the entire culture-flask surface.
5. Incubate for 6 minutes.
6. Tap side of the flask to loosen the cells. Add 10 mL of feeder culture medium to neutralize the trypsin.
7. Aspirate cell suspension into conical tube.
8. Centrifuge for five minutes at 2000 rpm.
9. Remove supernatant, add feeder culture medium and pipette in order to re-suspend the pellet.
10. Count cells and re-suspend in desired medium volume.
11. Add cell suspension into gelatin pre-coated culture dishes.
We recommend 3.5×10^5 cells per well in six-well plate (2 mL/well).
12. Incubate the feeder cells for at least two hours before plating the hPSC (cells must be spread).
13. When plating hPSC, change the feeder medium to NutriStem® hPSC XF medium.

NOTE

- Feeder concentration can also be calculated as 3.5×10^4 cells per cm^2 .
- Do not use NutriStem® hPSC for MEF culture. Prepare MEF-covered plate using recommended culture medium and change the medium before plating the cells.
- Plates may be used within 10 days of preparation.
- It is highly recommended to prepare MEF plates 24 hours before use.

Passage of hPSC

Collagenase solution

Dissolve Collagenase type IV in DMEM/F-12 (1:1) medium (Cat. No. 01-170-1) to a concentration of 1 mg/ml and filter sterilize with a 0.22 μm pore size filter unit. Collagenase solution can be stored at 2–8°C for up to 2 weeks.

Protocol

1. Remove medium from well. Add 1.0 mL collagenase solution.
2. Incubate at 37°C or at room temperature until the edges of the cell colonies begin to loosen from the plate.

NOTE Incubation times will vary between cell lines and colony sizes. Begin checking the culture after 3 minutes. Do not over-incubate the culture, as hPSC are sensitive to enzymatic stress and may lift from the plate during the wash step.

3. Aspirate the Dissociation Solution and wash the cells twice with 2 mL of sterile DMEM/F-12 (1:1).
4. Add 1 mL of culture medium and gently wash off and detach hPSC using 5ml pipette. Most of the feeder layer cells will remain on the plate.
5. Collect cell suspension and put into conical tube.
6. Centrifuge 3 minutes at 800 rpm at room temperature.
7. Re-suspend hPSC using NutriStem[®] hPSC XF medium by gently pipetting up and down with a 5ml pipet to break hPSC clumps and plate in feeder-covered plate with 2.5–3 mL NutriStem[®] hPSC XF.
8. Change the medium daily with 2.5–3.0 mL/well fresh NutriStem[®] hPSC XF until the hPSC colonies are large enough to passage.

NOTE For effective separation of hPSC from the feeder cells, longer dissociation time is recommended (up to 1 hr.).

2. Feeder-free culture

Matrigel™ based culture

Matrigel™ preparation

Matrigel™ should be thawed on ice, aliquoted and refrozen. For full instructions and recommendations on aliquot size, please refer to the insert supplied with the product.

Preparation of Matrigel™ Aliquots

1. Thaw Matrigel™ on ice at 2–8°C for overnight to avoid the formation of a gel.
2. Dilute Matrigel™ 1:1 with cold DMEM: F12 (1:1) medium (Cat. No. 01-170-1). Mix well with a cold pipette.
3. Keep mixture on ice.
4. Aliquot into pre-chilled 15 mL tubes (for aliquot stability, please refer to the insert supplied with the product). Store at -70°C.

Coating Plates with Matrigel™

1. Slowly thaw Matrigel™ aliquot on ice at 2–8°C, to avoid the formation of a gel.
2. Dilute the Matrigel™ aliquot 1:20 in cold DMEM: F12 (1:1).
3. Add 1 mL of Matrigel™ solution to each well of 6-well plate.
4. Incubate the plate for at least 1–2 hours at room temperature or overnight at 2–8°C. Plate with Matrigel™ solution can be parafilm® sealed and stored at 2–8°C for one week. Do not use the plate if the Matrigel™ solution does not completely cover the surface of the wells.
5. Before use, warm to room temperature, remove Matrigel™ solution and immediately cover with NutriStem® hPSC XF.

Enzymatic passaging of hPSC grown in NutriStem® hPSC XF on Matrigel™

We recommend using Collagenase IV at 1 mg/mL. If using other enzyme-based dissociation solutions (Dispase or recombinant trypsin), optimal conditions should be determined by the user.

Protocol

1. Aspirate the culture medium from the well.
2. Wash the well once with warm DMEM: F12 (1:1) medium (Cat. No. 01-170-1).
3. Add 1 mL per well of warmed Collagenase IV (1 mg/mL).
4. Return the plate to the 37°C 5% CO₂ incubator. Check the culture every 5–10 minutes. Do not over-expose the cells to Collagenase IV. Observe the colonies under the microscope. Incubate until the edges of the colonies begin to curl up but the colonies should remain attached to the plate surface.
5. Aspirate the Collagenase solution and wash very gently at least once with DMEM: F12 (1:1) medium (Cat. No. 01-170-1).
6. Add 2 mL of pre-warmed NutriStem[®] hPSC XF medium to the well.
7. Gently scrape and wash the colonies off with 5ml glass pipette.
8. Repeat the scraping and pipetting action 3–4 times until all the colonies have been removed from the well surface. Pipette gently to avoid breaking up the colonies into too small clumps.
9. Transfer the detached colonies into a sterile conical tube.
10. Use another 2 mL of medium to wash the well. Transfer the cell clumps into the same tube.

NOTE The split ratio of NutriStem[®] hPSC XF medium-based culture is usually 1:6 – 1:8 every 3 – 5 days.

11. To obtain the desirable split ratio, increase the volume of the clumps suspension in the tube with NutriStem[®] hPSC XF medium accordingly.
12. Prepare a new Matrigel[™]-coated 6-well plate by washing once with DMEM: F12 (1:1). Gently add 3 mL NutriStem[®] hPSC XF medium to each well to be seeded.
13. Mix the hPSC clumps suspension in the tube to break cell clumps by gently pipetting up and down with a 5 mL pipet.
14. Gently and evenly add 1ml of the suspended hPSC clumps to each new well to be seeded.
15. Place the plate in the 37°C 5% CO₂ incubator and carefully move the plate back and forth and side to side to evenly distribute the clumps throughout the well.
16. Allow the colonies to attach in 37°C 5% CO₂ incubator.

17. After 48 hours, change the medium daily with 2.5–3.0 mL/well fresh NutriStem® hPSC XF until the hPSC colonies are large enough to passage.
18. It is possible to perform higher volume feed (4–5 mL) on the weekend and the next medium change two days later.
19. If the culture is at optimal density, the cells can be split every 3–4 days using a 1:6–1:8 splitting ratio (colonies from 1 well of a 6–well plate can be plated in 6–8 new wells of a 6–well plate). If the colonies are too dense or too sparse, adjust the splitting ratio accordingly.

Enzyme-free passaging of hPSC grown in NutriStem® hPSC XF on Matrigel™

Dissociation solution

0.5mM EDTA in DPBS: prepare 0.5mM EDTA by combining 50 µL of 0.5M EDTA (Cat. No. 01-862-1), pH 8.0 with 50 mL of DPBS without calcium and magnesium (Cat. No. 02-023-1). Filter sterilize the solution and store at room temperature.

Protocol

This procedure describes the passage of the colonies as very small clumps. Volumes are for 1 well of 6–well plate.

1. Wash cells twice with 2 mL DPBS without calcium and magnesium (Cat. No. 02-023-1). Add 1 mL 0.5 mM EDTA solution, swirl the vessel to coat the entire cell surface and quickly discard.

NOTE Do not expose cells to the EDTA solution for more than needed for a quick wash at this point.

2. Add 1 mL of 0.5mM EDTA solution and incubate for 3–4 minutes at room temperature. Do not move the plate during EDTA exposure.
3. Gently remove the EDTA solution and add carefully 1 mL NutriStem® hPSC XF medium.
4. Detach and break colonies by gently pipetting up and down 3–4 times with a 1 mL tip. Make sure the pipetting washes the entire well.
5. Plate the cell aggregates at the desired density in Matrigel™ coated wells with 4 mL pre-warmed NutriStem® hPSC XF medium. Usually a splitting ratio of 1:8–1:10 every 4 days is required.

6. Place the plate in a 37°C 5% CO₂ incubator. Move the plate several times back and forth and side to side motions to distribute the aggregates evenly in the well.
7. After 48 hours, change the medium daily with 2.5–3.0 mL/well fresh NutriStem® hPSC XF until the hPSC colonies are large enough to passage.
8. It is possible to perform higher volume feed (4–5 mL) on the weekend and the next medium change two days later.

NOTE Do not move the plate during the first 48 hours post-split (this may increase differentiation of hPSC).

Laminin based culture

LaminStem™ 521 (Cat. No. 05-753-1) coating procedure using 0.5–1 µg/cm² in a 6-well plate. For a detailed protocol, follow laminin manufacturer instructions.

General coating protocol: (0.5 µg/cm²)

1. Slowly thaw recombinant laminin (LaminStem™ 521) at 2–8°C.
2. Thawed laminin stock may be stored at 2–8°C under aseptic conditions.
3. Dilute 300 µL LN521 (0.1 mg/mL) with 12 mL DPBS with calcium and magnesium, (50 µL/2 mL/well).
4. Add 2 mL/well of the diluted LaminStem™ 521.
5. Seal the plate with plastic film (e.g., Parafilm®) to prevent evaporation, and incubate overnight at 2–8°C. Make sure the laminin solution is evenly spreaded across the surface. Note that the laminin matrix will be inactivated if dried.

NOTE

- Optimal coating concentration is cell-dependent and should be calibrated. 1 µg/cm² should work well for most hPSC.
- Coated plates may be stored aseptically at 2–8°C.
- Rapid coating may be done at 37°C for 2 hours.

Single cell passage of hPSC cultured on LaminStem™ 521 (enzymatic procedure, volumes are for 1 well of 6-well plate)

1. Wash cells twice with 2 mL/well DPBS without calcium and magnesium
2. Add 1 mL of recombinant Trypsin-EDTA Solution (Cat. No. 03-079-1) and incubate at 37°C for 2–4 minutes. Exposure time may vary and should be adjusted.

NOTE Do not over-expose the cells to the recombinant Trypsin-EDTA Solution.

3. Detach the cells by gently pipetting up and down. Add 4 mL 1xSBT1 (Cat. No. 03-048-1) to the recombinant Trypsin-EDTA Solution.
4. Detach and break colonies by pipetting 5–6 times up and down with a 1 mL pipette. Make sure the pipetting washes the entire well.
5. Collect cells into new sterile tube.
6. Centrifuge at 200 g for 5 minutes.
7. Suspend the pellet with 1 mL of fresh medium.
8. Count cells.
9. Plate cells at the desired density in LaminStem™ 521 pre-coated wells with 3–4 mL pre-equilibrated medium. Usually, 10–20,000/cm² for splitting every 4–5 days.
10. Place the plate in a 37°C 5% CO₂ incubator. Move the plate several times in back and forth and side to side motions to evenly distribute the cells in the well.
11. Perform daily medium change after 48 hours with 2.5–3.0 mL/well fresh NutriStem® hPSC XF until the hPSC colonies are large enough to passage. Do not move the plate during the first 48 hours post-split.
12. It is possible to perform higher volume feed (4–5 mL) for the weekend and skip 2 days without changing the medium. In this case, the cells must be passaged at the end of the week.

NOTE

- It is critical not to move the seeded plates during the first 48 hours after splitting as this may increase differentiation of hPSC.
- Cells are ready to be passaged when cell culture is ≥60% confluent. Optimal seeding densities will vary from one cell line to another and must be determined empirically for your system. With optimal culture conditions and seeding density, most cell lines will reach confluency within 4–6 days and expand 10–25 folds.

7 Quality Control

NutriStem® hPSC XF Medium performance is tested for optimal maintenance of hESC (H1) for NLT 3 passages and IF Staining for pluripotency markers. Additional parameters tested are: pH, osmolality, endotoxins, mycoplasma and sterility tests. For full specifications please check the lot specific Certificate of Analysis (CoA).

8 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:
 - ISO 13408 – Aseptic Processing of Health Care Products
 - ISO 14644 – Airborne Particulate Cleanliness Classes in Clean Rooms and Clean Zones.
- A DMF (Drug Master File) has been acknowledged by the US-FDA.



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

8.1 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.
CE	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.

9 Related Products

Product	Cat. No.
Dulbecco's PBS (without Ca & Mg)	02-023-1
Dulbecco's PBS (with Ca & Mg)	02-020-1
Trypsin-EDTA Solution B (EDTA 0.05%, Trypsin 0.25%) with Phenol Red	03-052-1
Bio-Pure Human Serum Albumin (HSA), 10% solution	05-720-1
LaminStem™-521	05-753-1
Recombinatn Trypsin-EDTA Solution	03-079-1
0.1% Gelatin Solution	01-944-1
SBTI Solution (50X)	03-048-1
EDTA 0.5M Solution	01-862-1
NutriFreez® D10 Cryopreservation Medium	05-713-1

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:

01 | 2022

© 2022
Biological Industries Israel Beit Haemek Ltd.
2511500 Kibbutz Beit Haemek
Israel

AM | DIR: 2740227-000-02
Revision 03