

Culture of MSCs on Nunclon™ Supra Surface Cultureware

Catalog Numbers 150476, 140680, 167013, 156376, 156380, 140682, 150478, 140681, 156372, 150470, 150472, 150474, 140683, 156374, 156378, 156382

Pub. No. MAN1000717 Rev. A



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

The Thermo Scientific™ Nunclon™ Supra Surface treatment is used for culturing a variety of fastidious, adherent cells and supports serum-free and coating-free cultures of human mesenchymal stromal cells (MSCs). MSCs cultured on the Nunclon™ Supra Surface retain their morphology and downstream differentiation potential. Nunclon™ Supra Surface cultureware used in conjunction with Gibco™ StemPro™ MSC SFM complete medium is a xeno-free solution for cell therapy research applications.

Contents and storage

| Product | Cat. No. | Storage | Shelf Life ^[1] |
|------------------------------------|--|---------|---------------------------|
| Nunclon™ Supra Surface Cultureware | 150476, 140680, 167013, 156376, 156380, 140682, 150478, 140681, 156372, 150470, 150472, 150474, 140683, 156374, 156378, 156382 | Ambient | 5 years |

^[1] Shelf life is determined from the date of manufacture.

Product use

- Nunclon™ Supra Surface cultureware: For laboratory use.
- StemPro™ MSC SFM XenoFree: For Research Use Only. Not for use in diagnostic procedures.

Important information



CAUTION! Human origin materials are non-reactive (donor level) for anti-HIV 1 & 2, anti-HCV, and HBsAg. Handle in accordance with established bio-safety practices.

Procedural guidelines

- Thaw StemPro™ MSC SFM XenoFree Supplement overnight at 2°C to 8°C prior to use (thawed supplement will have a slightly cloudy appearance). Use thawed material immediately or aliquot (i.e., 1 mL) unused material and store at -20°C to -5°C. Avoid additional freeze-thaw cycles.
- StemPro™ MSC SFM XenoFree complete medium (StemPro™ MSC SFM Basal Medium, StemPro™ MSC SFM XenoFree Supplement, and CTS™ GlutaMAX™ Supplement) is stable for 2 weeks when stored in the dark at 2°C to 8°C.

Prepare complete medium

StemPro™ MSC SFM Basal Medium requires supplementation with StemPro™ MSC SFM XenoFree supplement and CTS™ GlutaMAX™ Supplement or L-Glutamine.

- For 500 mL complete medium, aseptically add 5 mL of StemPro™ MSC SFM XenoFree supplement to StemPro™ MSC SFM Basal Medium (500 mL).
- Aseptically add 5 mL CTS™ GlutaMAX™ Supplement or 200 mM L-Glutamine to the complete medium before use.
- Optional:* add 50 µL of 50 mg/mL Gentamicin to the complete medium.

Culture conditions

Culture vessels: Nunclon™ Supra Surface cultureware

Media: StemPro™ MSC SFM XenoFree complete medium

Cell line(s): Human mesenchymal stem cells (MSCs) or adipose-derived stem cells (ADSCs)

Culture type: Adherent

Temperature range: 36°C to 38°C

Incubator atmosphere: Humidified atmosphere of 4–6% CO₂ in air. Ensure proper gas exchange and minimize exposure of cultures to light.

Note: Procedures detailed in the following sections are for cultures in T-75 culture flasks (75 cm²). Volumes should be adjusted accordingly for the desired vessel size.

Recover cryopreserved human MSCs

1. Rapidly thaw a frozen vial of human MSCs in a 37°C water bath until a small amount of ice remains.
2. Pipet the entire contents of the cryovial into a 50-mL conical tube.
3. Carefully add 5–10 mL of pre-warmed (37°C) StemPro™ MSC SFM XenoFree complete medium to the conical tube at an approximate rate of 3 to 5 drops per 5 seconds and gently swirl after every addition.
4. Centrifuge the tubes at 100–200 × *g* for 5 minutes at room temperature.
5. Resuspend the cell pellet in pre-warmed (37°C) StemPro™ MSC SFM XenoFree complete medium and add the cell suspension to an appropriate Nunclon™ Supra Surface flask at a density of $\geq 5 \times 10^3$ cells/cm²

Note: For the initial isolation of MSCs, supplementation of the complete medium with 2.5% Human Serum, Type AB facilitates cell attachment and growth. For subsequent passages, Human Serum, Type AB is not required

6. Incubate at 36°C to 38°C in a humidified atmosphere of 4–6% CO₂ in air.
7. Replace the medium in the flasks every 2 days.

Guidelines for subculturing cells in StemPro™ MSC SFM XenoFree complete medium

- StemPro™ MSC SFM XenoFree has been developed for the multi-passage expansion of human bone marrow-derived MSCs and adipose-derived stem cells (ADSCs) at greater than clonal densities ($\geq 5 \times 10^3$ cells/cm²). Reduced seeding densities may result in suboptimal cell expansion, although optimal growth conditions must be determined for each application.
- When subculturing human MSCs in StemPro™ MSC SFM XenoFree, input cell confluence should be 60–90%, cell viability should be at least 90% and the growth rate should be in mid-logarithmic phase prior to subculture. Initiating cultures under suboptimal conditions may affect product performance. Transitioning MSCs or ADSCs from serum-containing medium to StemPro™ MSC SFM XenoFree does not require an adaptation protocol.
- For optimal performance, re-feed the cultures every 2 days with StemPro™ MSC SFM XenoFree complete medium

Note: Procedures detailed on the next section are for cultures maintained in a Nunclon™ Supra Surface T-75 culture flask (75 cm²). Adjust the volumes accordingly for desired vessel size.

Propagate cells on Nunclon™ Supra Surface cultureware

1. Observe the stock culture flask under the microscope and confirm that the cells are ready to be sub-passaged (~60–90% confluent).
2. Pre-warm CTS™ TrypLE™ Select Enzyme and StemPro™ MSC SFM XenoFree complete medium to 37°C before use.
3. Add 10 mL of pre-warmed StemPro™ MSC SFM XenoFree complete medium to a 50-mL conical tube for each flask being harvested.
4. Remove the spent medium from the T-75 flask and discard.
5. Wash the cell surface with 10 mL of CTS™ DPBS without calcium chloride, without magnesium chloride, remove and discard.
6. Add 3–5 mL of CTS™ TrypLE™ Select Enzyme to the T-75 flask and tilt the flask in all directions to evenly distribute. Incubate the cells in CTS™ TrypLE™ Select Enzyme for 2–10 minutes in the incubator.

Note: Cells coming out of serum-containing medium may require a longer incubation time (5–10 minutes), while cells growing under serum-free conditions should detach more readily (2–3 minutes).

7. After incubation, check the flask under the microscope for cell detachment. Firmly tap the flask as necessary to facilitate complete cell detachment.
8. Add 7 mL of pre-warmed StemPro™ MSC SFM XenoFree to each flask and collect the cell suspension in the 50-mL conical tube containing complete medium from step 3. Firmly tap the flask, re-wash with 10 mL StemPro™ MSC SFM XenoFree and collect.
9. Centrifuge the tubes at 100–200 × *g* for 5 minutes at room temperature.
10. Resuspend the cells in a minimal volume of pre-warmed StemPro™ MSC SFM XenoFree complete medium for cell counting, using a preferred counting method (e.g., Countess™ 3 FL Automated Cell Counter).
11. Add 15 mL of StemPro™ MSC SFM XenoFree complete medium to as many new Nunclon™ Supra Surface flasks as desired.

12. Add enough cell suspension to each flask to provide $\geq 5 \times 10^3$ cells/cm² (i.e. 3.75×10^5 cells per T-75 flask). Mix or swirl the cell suspension to ensure even distribution.
13. Place the culture flask(s) in the incubator at 36°C to 38°C with a humidified atmosphere of 4–6% CO₂.
14. Replace the spent culture medium every 2 days with pre-warmed StemPro™ MSC SFM XenoFree complete medium.

Cryopreservation

1. Prepare cryopreservation solution (2X) by supplementing StemPro™ MSC SFM XenoFree complete medium with 10–20% Dimethyl Sulfoxide (DMSO). Store at 4°C until use; make cryopreservation medium on day of intended use.
2. Reconstitute the harvested cell pellet to twice the desired final concentration (i.e., 2×10^6 cells/mL) in pre-warmed StemPro™ MSC SFM XenoFree complete medium.
3. Slowly add cryopreservation solution to the cell suspension, and gently mix to ensure even cell distribution.
4. Immediately add the desired volume of cell suspension (i.e., 1 mL) to pre-chilled (2°C to 8°C) cryovials.
5. Place the cryovials at -70°C in a cryogenic freezing container (e.g., “Mr. Frosty™ Freezing Container (1°C)”).
6. After 24 hours, transfer the frozen cells to liquid nitrogen (vapor phase); storage at -200°C to -125°C is recommended.

Related products

Unless otherwise indicated, all materials are available through thermofisher.com.

| Product | Catalog No. |
|--|---|
| StemPro™ MSC SFM XenoFree | A1067501 |
| CTS™ GlutaMAX™ Supplement | A1286001 |
| L-Glutamine (200 mM) | A2916801 |
| Gentamicin | 15710064 |
| CTS™ TrypLE™ Select Enzyme | A1285901 |
| CTS™ DPBS without calcium chloride, without magnesium chloride | A1285601 |
| Human Serum, Type AB, Fisher BioReagents | BP2525100 |
| Thermo Scientific™ Nunc™ EasYFlask™ Flasks with Nunclon™ Supra Surface: <ul style="list-style-type: none">• Filter cap: T-25, T-75, T-175• Solid cap; T-25, T-75, T-175 | 156372 , 156376 , 156380 , 156374 , 156378 , 156382 |
| Thermo Scientific™ Nunc™ Multidishes with Nunclon™ Supra Surface 12-well, 6-well, 24-well, 48-well | 140681 , 140680 , 140682 , 140683 |
| Thermo Scientific™ Nunc™ EasYDish™ Dishes with Nunclon™ Supra Surface, 35-mm, 60-mm, 100-mm, 100-mm, 150-mm | 150470 , 150472 , 150474 , 150476 , 150478 |
| Countess™ 3 FL Automated Cell Counter | AMQAF2001 |

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Revision history: Pub. No. A MAN1000717

| Revision | Date | Description |
|----------|------------------|------------------|
| A | 21 November 2024 | Initial release. |

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.