

Instructions for Use

# 4Cell<sup>®</sup> XtraCHO Media System

A Chemically Defined and Animal Component Free  
Cell Culture Media for any CHO Cell Lines

2557820-000-04



**SARTORIUS**



# Contents

<b>1</b>	<b>Introduction</b>	<b>4</b>
1.1	Safety Statements	4
1.2	Stability	5
1.3	Unpacking and Storage Instructions	5
1.4	Recommended Materials	5
<b>2</b>	<b>Instructions for Use</b>	<b>6</b>
2.1	Adapting Cell Lines to 4Cell® XtraCHO Medium	6
2.2	Cell Cultivation	7
2.3	Thawing of Cells   Initiation of Culture Process	7
2.4	Freezing of Cells   Storage	8
2.5	Protein Production in Fed-batch Mode	8
2.6	Fed-batch Protocol	9
<b>3</b>	<b>Instructions for Reconstitution of 4Cell® XtraCHO Media Powder (Optional)</b>	<b>12</b>
3.1	4Cell® XtraCHO Stock & Adaptation Medium	12
3.2	4Cell® XtraCHO Production Medium	13
3.3	4Cell® XtraCHO Feed Medium A	13
3.4	4Cell® XtraCHO Feed Medium B	14
<b>4</b>	<b>References   Contacts</b>	<b>15</b>

# 1 Introduction

4Cell® XtraCHO Media System consists of four different media:

- 4Cell® XtraCHO Stock & Adaptation Medium (SAM)
- 4Cell® XtraCHO Production Medium (PM)
- 4Cell® XtraCHO Feed Medium A (FMA)
- 4Cell® XtraCHO Feed Medium B (FMB)

These media are chemically defined non-animal origin growth media formulated to maximize the product titer in CHO cell lines, optimized for CHO DG44 cells. It provide robust performance in production systems of small and large scale, e.g. flask, Ambr® 15/250 and bioreactors.

## Product Characteristics

- Complete medium
- Animal Component Free (ACF) medium
- Serum-free (SF) medium
- Does **not** require the addition of serum.
- Does **not** contain antibiotics.
- The medium requires the addition of Glutamine.
- Available in as non-sterile powder and as sterile liquid.

## 1.1 Safety Statements

**For the intended use, please refer to Certificate of Analysis.**

**Not** approved for human or veterinary use. **Not** for application in humans or animals, or for use in vitro diagnostic or clinical procedures.

## 1.2 Stability

4Cell® XtraCHO Stock & Adaptation Medium (SAM), and 4Cell® XtraCHO Production Medium (PM) needs to be supplemented with L-Glutamine prior to use following the instructions for use, usually at 4 – 6 mM final concentration. After addition of L-Glutamine, the media are stable for 4 weeks when stored at 2 – 8°C protected from light.

Both feeds, 4Cell® XtraCHO Feed Medium A (FMA) and 4Cell® XtraCHO Feed Medium B (FMB), are ready for use and do **not** require L-Glutamine supplementation.

## 1.3 Unpacking and Storage Instructions

1. Check all containers for leakage or breakage.
2. When **not** in use store 4Cell® XtraCHO Media System components at 2°C to 8°C protected from light.

## 1.4 Recommended Materials

- DMSO for cell preservation
- 125 mL Erlenmeyer flask, Ambr® 15/250, 2 L – 50 L bioreactors
- 100 – 400 g/L sterile filtered Glucose
- L-Glutamine, 200 mM
- PES membrane filter with 0.2 µm pore size, e.g. Sartopore® 2

## 2 Instructions for Use

### 2.1 Adapting Cell Lines to 4Cell® XtraCHO Medium

In general, there are two approaches to adapt a cell line to a new culture medium formulation.

- **Option 1:** Passage directly the culture directly from the initial medium into 4Cell® XtraCHO Stock & Adaptation Medium (SAM). Choose a high seeding cell density at each passage (e.g.  $5 \times 10^5$  cells/mL) for a minimum of two weeks. When the cells achieve a stable growth rate and viability >90% for 2 passages, the adaptation is considered complete.
- **Option 2:** Passage the culture into a mixture of original culture medium and 4Cell® XtraCHO SAM and gradually increase the content of 4Cell® XtraCHO SAM. An example for a step-wise adaptation protocol is given below.

Adaptation step	Ratio of original medium to 4Cell® XtraCHO SAM	Acceptance criterion to proceed to next adaptation step
1	75:25	Viability $\geq$ 90% of original medium normal doubling time for 2 passages
2	50:50	Viability $\geq$ 90% of original medium normal doubling time for 2 passages
3	25:75	Viability $\geq$ 90% of original medium normal doubling time for 2 passages
4	0:100	Adaptation complete if viability >90% in 4Cell® XtraCHO SAM medium normal doubling time for 2 passages

## 2.2 Cell Cultivation

- Cultivate the cells in an incubator with a shaking platform and humidified atmosphere, containing  $7.5\% \pm 0.5\% \text{CO}_2$ .
- Other cultivation parameters may be adapted to each cell line's individual requirements.
- A recommended starting point is a temperature of  $36.8^\circ\text{C} \pm 0,2^\circ\text{C}$  and 103 rpm\* on an orbital shaking platform.
- By regular passaging of the cells, ensure that the culture remains in midexponential growth phase at all times. Determine cell density and viability of the culture every 1 – 2 days and dilute the culture to a suitable seeding density with fresh pre-warmed medium (e.g.  $3 \times 10^5$  viable cells/mL).

## 2.3 Thawing of Cells | Initiation of Culture Process

1. Pre-warm 4Cell® XtraCHO Stock & Adaptation Medium (SAM) to cultivation temperature before use. The required medium volume depends on the cell density in frozen cryovials. The cell density after thawing should be  $3 - 5 \times 10^5$  viable cells/mL.
2. After removing cryovial from storage, wipe the cryovial with 70% v/v ethanol or isopropanol before opening. In a Biological Safety Cabinet (BSC), briefly twist the cap a quarter turn to relieve pressure, and then retighten.
3. Quickly thaw the cryovial in a  $37^\circ\text{C}$  water bath (do **not** submerge the cryovial completely) or heating block at  $37^\circ\text{C}$  until only a small grain of ice remains. Thawing the cells for longer than 3 minutes may result in reduced cell viability.
4. Dry the cryovial with a lint-free wipe, spray with 70% v/v ethanol or isopropanol, and then wipe to remove excess liquid.
5. Immediately transfer the thawed cell suspension with a pipette into 10 mL of 4Cell® XtraCHO SAM and centrifuge at  $180 - 200 \times g$  for 3 minutes. Remove the supernatant carefully.
6. Carefully reconstitute the cell pellet in fresh pre-warmed 4Cell® XtraCHO SAM by gently mixing by pipetting up and down.
7. Transfer the suspension as inoculum into the culture vessel. Proceed with cell cultivation as described above.

\* Shaking rate for an Infors Multitron cell incubator with 50 mm orbital diameter.  
For shakers with other orbital diameters: shaking rate in rpm =  $103 \times (50/\text{orbital diameter})^{1/2}$

## 2.4 Freezing of Cells | Storage

The cell culture should be in mid-logarithmic growth phase and >90% viable at the point of freezing.

1. Prepare the necessary volume of freezing medium by supplementing 4Cell® XtraCHO Stock & Adaptation Medium (SAM) with 7.5% Dimethylsulfoxide (DMSO). L-Glutamine can be added optionally (4 – 6 mM final concentration) to the freezing medium, but is **not** necessary. Store the freezing medium at 2 – 8°C until use.
2. Transfer the required volume of cell suspension into centrifugation vessels and spin down the cells at 180 – 200 × g for 5 minutes. Gently remove the supernatant.
3. Reconstitute the cell pellet in the required volume of freezing medium to achieve a cell density of at least  $1 \times 10^7$  viable cells/mL. Dispense the suspension into cryovials, taking care that the suspension remains homogeneous at all times.
4. Use a suitable controlled cooling method to freeze the vials, ideally a controlled-rate freezer. Alternatively, place them in a cell freezing container overnight at -80°C. For storage keep the vials at a temperature below -130°C preferably in vapor phase LN2 for frequent access and in liquid nitrogen freezer for long-term storage.

## 2.5 Protein Production in Fed-batch Mode

The 4Cell® XtraCHO Media System Starting Kit includes all necessary medium components for adaptation of CHO cell lines, cultivation from stock culture of adapted cells up to 3 L of fed-batch cultivation in 4Cell® XtraCHO Production Medium (PM) plus Feed Medium A (FMA) and Feed Medium B (FMB). Stirred bioreactors and Sartorius Ambr® 15/250 are ideally suited to carry out comprehensive evaluation of fed-batch parameter in the course of process development.

The Table 1 below outlines volumes needed for each individual medium component dependent on the target cultivation volume from seed train to the end of fed-batch culture.

## NOTE

The actual volume needed for adaptation to 4Cell® XtraCHO Stock & Adaptation Medium is dependent on the cell type and adaptation method used.

Cultivation Volume	Liquid Volumes				Cultivation Volume	Liquid Volumes			
	SAM	PM	FMA	FMB		SAM	PM	FMA	FMB
L	L	L	L	L	L	L	L	L	L
2	0.52	1.20	0.60	0.08	5	1.3	3	1.5	0.2
2.5	0.65	1.50	0.75	0.10	10	2.6	6	3	0.4
3	0.78	1.80	0.90	0.12	15	3.9	9	4.5	0.6
4	1.04	2.40	1.20	0.16	20	5.2	12	6	0.8

Table 1

## 2.6 Fed-batch Protocol

The following protocol is best carried out in a stirred tank bioreactor. Alternatively, Ambr® 15/250 systems or shake flasks can be used.

After successful adaptation, seed the cells in the desired volume of completed 4Cell® XtraCHO Production Medium (PM) at a suitable viable cell density (recommended:  $3 \times 10^5$  cells/mL). If the volume of inoculum is less than 20% of the final volume, cells can be transferred directly from 4Cell® XtraCHO SAM to 4Cell® XtraCHO PM; otherwise it is recommended to centrifuge the cells to remove the used 4Cell® XtraCHO SAM and reconstitute them in fresh 4Cell® XtraCHO PM.

From day 3 after inoculation, perform daily sampling and monitoring of cell density, viability, product titer and key metabolites.

Begin the addition of 4Cell® XtraCHO Feed Medium A (FMA) and 4Cell® XtraCHO Feed Medium B (FMB) on day 3 after inoculation, or when the viable cell density reaches  $2 \times 10^6$  cells/mL. Three suggested feeding options for a fed-batch are outlined in the Tables 2a-c below. Other cultivation conditions are the same as outlined above in paragraph “Cell cultivation”. Add 4Cell® XtraCHO FMA at 2 – 6% of original culture volume and 4Cell® XtraCHO FMB at 10% of Feed Medium A volume. Add both feed media slowly and ensure that the feed media are quickly dispersed within the cultured cells. FMB has a pH of 10 – 11, so its addition may lead to a short spike in culture pH.

4Cell® XtraCHO FMA contains sufficient glucose to supply the fed-batch culture until day 5 after inoculation. Glucose concentration should be measured daily and kept at a value of 2 – 6 g/L by addition of a separate concentrated solution (recommended concentration: 400 g/L) when required.

End the cultivation when a pre-determined end criterion is reached, e.g. a specific time point after inoculation or when viability falls below 50 – 70%.

Table 2: Three suggested options for a fed batch procedure to determine the optimal daily feeding rate. The ratio of Feed Medium A to Feed Medium B should always be 10:1. Values are calculated as percentages of starting values at day 0.

Day	0	1	2	3	4	5	6	7	8	9	10	11	12	13
FMA (%)	-	-	-	2	2	2	2	2	2	2	2	2	2	2
FMB (%)	-	-	-	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2	0.2
Glucose	-	-	-	-	-	To target concentration, 4 – 6 g/L								

Table 2a) Option 1

Day	0	1	2	3	4	5	6	7	8	9	10	11	12	13
FMA (%)	-	-	-	4	4	4	4	4	4	4	4	4	4	4
FMB (%)	-	-	-	0.4	0.4	0.4	0.4	0.4	0.4	0.4	0.4	0.4	0.4	0.4
Glucose	-	-	-	-	-	To target concentration, 4 – 6 g/L								

Table 2b) Option 2

Day	0	1	2	3	4	5	6	7	8	9	10	11	12	13
FMA (%)	-	-	-	6	6	6	6	6	6	6	6	6	6	6
FMB (%)	-	-	-	0.6	0.6	0.6	0.6	0.6	0.6	0.6	0.6	0.6	0.6	0.6
Glucose	-	-	-	-	-	To target concentration, 4 – 6 g/L								

Table 2c) Option 3

## IMPORTANT

- For cells depending on continuous supply with L-Glutamine the cultures may need to be supplemented with additional L-Glutamine to prevent its depletion.
- Store 4Cell® XtraCHO FMB in an airtight container when in use. If solid crystals become visible in a container of FMB, discard and replace it with fresh 4Cell® XtraCHO FMB. If FMB is exposed to ambient oxygen and temperature during culture replace every three days with fresh 4Cell® XtraCHO FMB from storage. e.g. in Ambr® 15 & 250 systems.

### 3 Instructions for Reconstitution of 4Cell® XtraCHO Media Powder (Optional)

The media is packaged and intended to be hydrated completely from one container.

#### 3.1 4Cell® XtraCHO Stock & Adaptation Medium

1. Fill WFI (Water For Injection) into the appropriate mixing vessel. The WFI should be at room temperature. To allow pH adjustment later on, the volume should be 97% of the final volume.
2. Add 20.04 g/L of the media powder 4Cell® XtraCHO Stock & Adaptation Medium and stir rapidly for 60 min at room temperature.
3. Without suspending stirring, stepwise add 3 mL of 2 M NaOH solution per liter medium and continue to stir for 30 min at room temperature.
4. Add 1.80 g/L NaHCO<sub>3</sub> and stir until completely dissolved (~15 min).
5. If required, adjust the pH to 6.90 – 7.30 by adding 2 M NaOH.
6. Add WFI to the final volume. The osmolality value of the liquefied 4Cell XtraCHO Stock & Adaptation Medium is expected to stand within 270 – 330 mOsmol/kg H<sub>2</sub>O.
7. Sterile filter the medium using a PES membrane filter with 0.2 µm pore size (Sartopore® 2).

Store at 2 – 8°C. Protect from light.

## 3.2 4Cell® XtraCHO Production Medium

1. Fill WFI (Water For Injection) into the appropriate mixing vessel. The WFI should be at room temperature. To allow pH adjustment later on, the volume should be 97% of the final volume.
2. Add 22.31 g/L of the media powder 4Cell® XtraCHO Production Medium and stir rapidly for 60 min at room temperature.
3. Without suspending stirring, stepwise add 6.2 mL of 5 M NaOH solution per liter medium and continue to stir for 30 min at room temperature.
4. Add 1.80 g/L NaHCO<sub>3</sub> and stir until completely dissolved (~15 min).
5. If required, adjust the pH to 6.90 – 7.35 by adding 2 M NaOH.
6. Add WFI to the final volume. The osmolality value of the liquefied 4Cell® XtraCHO Production Medium is expected to stand within 275 – 330 mOsmol/kg H<sub>2</sub>O.
7. Sterile filter the medium using a PES membrane filter with 0.2 µm pore size (Sartopore® 2).

Store at 2 – 8°C. Protect from light.

## 3.3 4Cell® XtraCHO Feed Medium A

1. Fill WFI (Water For Injection) into the appropriate mixing vessel. The WFI should be at room temperature. To allow pH adjustment later on, the volume should be 85% of the final volume.
2. Add 154.12 g/L of the media powder 4Cell® XtraCHO Feed Medium A and stir rapidly for 60 min at room temperature.
3. Without suspending stirring, stepwise add 15.7 mL of 5 M NaOH solution per liter medium and continue to stir for 30 min at room temperature.
4. If required, adjust the pH to 6.50 – 6.80 by adding 2 M NaOH.
5. Add WFI to the final volume. The osmolality value range of the liquefied 4Cell® XtraCHO Feed Medium A is expected to stand within 208 – 268 mOsmol/kg H<sub>2</sub>O, measured at a 1:5 dilution.
6. Sterile filter the medium using a PES membrane filter with 0.2 µm pore size (Sartopore® 2).

Store at 2 – 8°C. Protect from light.

## 3.4 4Cell® XtraCHO Feed Medium B

1. Fill WFI (Water For Injection) into the appropriate mixing vessel. The WFI should be at room temperature. To allow pH adjustment later on, the volume should be 75% of the final volume.
2. Add 94.60 g/L of the media powder 4Cell® XtraCHO Feed Medium B and stir rapidly for 60 min at room temperature.
3. Without suspending stirring, stepwise add 160 mL of 5 M NaOH solution per liter medium and continue to stir for 30 min at room temperature.
4. If required, adjust the pH to 11.00 – 11.40 by adding 5 M NaOH.
5. Add WFI to the final volume. The osmolality value range of the liquefied 4Cell® XtraCHO Feed Medium B is expected to stand within 218 – 266 mOsmol/kg H<sub>2</sub>O, measured at a 1:5 dilution.
6. Sterile filter the medium using a PES membrane filter with 0.2 µm pore size (Sartopore® 2).

Store at 2 – 8°C. Protect from light.

# 4 References | Contacts

Europe | Asia: [CellCultureMedia.EU@sartorius.com](mailto:CellCultureMedia.EU@sartorius.com)

Americas: [CellCultureMedia.NA@sartorius.com](mailto:CellCultureMedia.NA@sartorius.com)

Or at: [Info\\_CellCultureMedia@sartorius.com](mailto:Info_CellCultureMedia@sartorius.com)

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

Phone: +49 551 308 0  
[www.sartorius.com](http://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:

12 | 2022

© 2022

Sartorius Stedim France S.A.S.  
Zone Industrielle les Paluds, Avenue de Jouques CS 71058  
13781 Aubagne Cedex, France

LM | DIR: 2557820-000-04