

Xuri T Cell Expansion Medium

Instructions for use

Introduction

Product codes

29185230

29185231

Description

Xuri™ T Cell Expansion Medium is a chemically defined medium for *ex vivo* expansion of human T lymphocytes. It promotes proliferation of CD3/CD28 activated and Interleukin-2 stimulated human T lymphocytes in static culture as well as high density cultures in Xuri Cellbag™ bioreactors with the use of human serum. The product is for single use only.

Important

Read these instructions carefully before using the products.

Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

Meets USP<1043>. For further manufacturing only. Single use only. This product meets USP<1043> 'ancillary materials for cell, gene, and tissue-engineered products', within the responsibilities applicable to a supplier. Cytiva cannot fulfil USP<1043> in regards of application and therapy specific aspects (e.g. use in finished therapeutic, assessment of removal from a finished therapeutic and possibly biocompatibility, cytotoxicity or adventitious agent testing).

Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

Shipping

The product is shipped with ice packs.

Storage and stability

Do not freeze. Upon receipt, store the product at 2°C to 8°C. Protect from light. Please contact your Cytiva representative for stability information.

Expiry

See product label and Certificate of Analysis for lot specific information.

Volume size

Product code	Unit size ¹
29185230	500 mL
29185231	1000 mL

¹ variation in filling volume for 500 mL (494 mL to 546 mL) and 1000 mL (999 mL to 1041 mL) do apply.

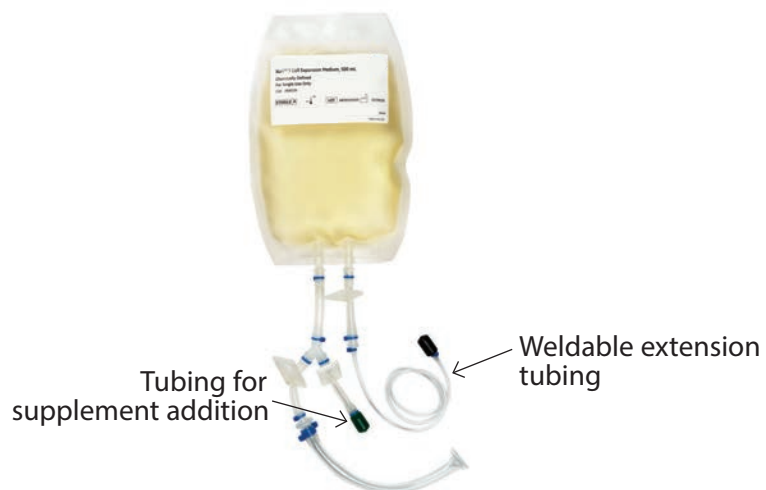
Components

- Chemically defined, human-and animal-derived component free (ADCF) medium.
- Free from antibiotics and phenol-red.
- Contains purified plant extracts that are characterized, recombinant human serum albumin (rec HSA) and L-Alanyl-L-Glutamine.

Preparation

Under aseptic conditions, use one of the tubings in the Y-connection to add supplements (human serum and growth factors) to the medium bag by attaching a needleless syringe to the Luer lock port.

Use the weldable extension tubing to connect Cell Expansion System Cellbag bioreactor by sterile welding using an appropriate tube welder.



Culture conditions

Xuri T Cell Expansion Medium is established and recommended to be used with human serum and growth promoting cytokines (e.g. Xuri IL-2). The use and amount of human serum and cytokines should be determined according to each application. Culture media should be equilibrated to room temperature for at least 2 hours prior to use.

Complete Media:

Xuri T Cell Expansion Medium + human serum + Xuri IL-2

Cells:

Peripheral blood mononuclear cells (PBMC)

Culture:

Static tissue culture vessel in a humidified incubator at 37°C, 5% CO₂.

Xuri Cellbag bioreactor in a Xuri Cell Expansion System W25 or W5 at 37°C, 5% CO₂.

Related products

Product	Product code
Xuri IL-2	29062790
Xuri IL-15	29112116
Xuri IL-21	29112119
Xuri Cell Expansion System W5	29036837
Xuri Cell Expansion System W25	29064568
Xuri Cell Expansion System Cellbag 2L Perfusion	29108442
Xuri Cell Expansion System Cellbag 2L Perfusion pH and DO	29105498
Xuri Cell Expansion System Cellbag 10L Perfusion	29108443
Xuri Cell Expansion System Cellbag 10L Perfusion pH and DO	29105499
Xuri Cell Expansion System Cellbag Basic 2L	29105492
Xuri Cell Expansion System Cellbag Basic 10L	29105493
Xuri Cell Expansion System Cellbag 2L pH and DO	29105494
Xuri Cell Expansion System Cellbag 10L pH and DO	29105495

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Cellbag and Xuri are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

All other third-party trademarks are the property of their respective owners.

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

29221749 AC V:6 03/2021

