

T Cell Serum Replacement Media Supplement

Description

T Cell Serum Replacement (SR) Media Supplement is designed to support the growth and expansion of human T cells in a cell culture medium containing only human-derived or human recombinant proteins. T Cell SR Media Supplement is xenofree and does not contain bovine or other non-human, animal-derived components.

Product	Catalog No.	Amount	Storage*	Shelf Life†
T Cell Serum Replacement Media Supplement	A24906DG	100 mL	Store at -5°C to -20°C. Protect from light.	N/A

* Store T Cell Serum Replacement Media Supplement in a **non-frost-free** freezer at -5°C to -20°C.

† Product is under development and the shelf life has not been validated.

Intended Use

For Investigational Use Only. The performance characteristics of this product have not been established.

Important Information

For best results, pre-equilibrate complete medium to temperature (36°C to 38°C) and gases (4–6% CO₂ in humidified air) before use. Ensure that proper gas-exchange is achieved in culture vessels. Avoid overexposure of cultures to light.

Safety Information

Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

T Cell Culture Conditions

Media: Complete medium comprised of a basal T cell media or T cell media kit (AIM V[®] Medium CTS™, OpTmizer™ CTS™, or other supportive media) containing T Cell SR Media Supplement.

Culture Type: Suspension

Culture Vessels: T-flasks or cell culture bag

Temperature Range: 36°C to 38°C

Incubator Atmosphere: Humidified atmosphere of 4–6% CO₂. Ensure that proper gas exchange is achieved in culture vessels.

Thaw T Cell SR Media Supplement

- To thaw T Cell SR Media Supplement, place it at 2°C to 8°C overnight. The next day, place the bottle in a 37°C water bath and continue thawing until the ice is gone and the supplement is clear. Gently swirl the bottle occasionally while in the water bath.

Alternatively, frozen T Cell SR Media Supplement can be thawed in a 37°C water bath, as long as the bottle is swirled frequently throughout the thawing process.

- Store thawed T Cell SR Media Supplement at 2°C to 8°C in the dark for up to 4 weeks.

Note: It is normal to see some flocculent material in the T Cell SR Media Supplement while thawing, but this material will go into solution with gentle swirling at 37°C. Minimize dwell time in water bath.

Use

- T cell cultures may be grown in Complete Medium comprised of a basal T cell medium or T cell media kit (AIM V[®] Medium CTS™, OpTmizer™ CTS™ T-cell Expansion SFM, or other supportive media) containing T Cell SR Media Supplement. As all T cell culture methods and cell sources may vary, it is recommended that the user determine what concentration of T Cell SR Media Supplement is optimal for their culture system. It is recommended that testing should be carried out with media containing a range of T Cell SR Media Supplement content, with 2%, 5%, and 10% spanning the likely functional ranges.
- Some media (such as OpTmizer™ CTS™ T-Cell Expansion Basal Medium) require the addition of glutamine and growth factors (e.g., IL-2) to support growth. Please follow standard media protocols provided by the media vendor, or your own internal protocols developed for your specific process. It is recommended that T Cell SR Media Supplement be tested against users' standard medium, where the user may normally add a serum supplement to the medium (e.g., 5% heat-inactivated pooled human Type AB serum).
- T Cell SR Media Supplement is supplied as a sterile-filtered solution. However, in order to reduce the risk of contamination during formulation of complete medium, filtration of complete medium containing T Cell SR Media Supplement may be performed (e.g., 0.2 µm). Users should test to ensure filtered complete medium continues to support T cell culture and expansion.
- T Cell SR Media Supplement in complete medium is stable for at least 10 days when stored in the dark at 2°C to 8°C. Avoid repeated warming and chilling of the complete medium. Warm only the volume required for that day's use.
- T Cell SR Media Supplement has been shown to work well in standard rapid expansion protocols using feeder cells, as well as Dynabeads[®] CD3/CD28 CTS™ bead-based expansion systems. T Cell SR Media Supplement may also work in other T cell culture protocols, although it has not been tested for such applications (e.g., NK T cells, regulatory T cells). T Cell SR Media Supplement has also been shown to support T cell gene-transduction protocols using lentiviral vectors.

Technical Support

For technical questions, contact us at the numbers below.

Sandy Kuligowski: sandra.kuligowski@lifetech.com











Karoline Schjetne: karoline.schjetne@lifetech.com

Related Products

Product	Cat. No.
AIM V [®] Medium CTS [™]	0870112
OpTmizer [™] CTS [™] T-cell Expansion SFM	A10485
AB-Human Serum	34005-100
Dynabeads [®] CD3/CD28 CTS [™]	40203D
L-Glutamine, 200 mM (100X), liquid	25030

Explanation of Symbols and Warnings

The symbols present on the product label are explained below:

				
Caution, consult accompanying documents	Temperature Limitation	Keep away from light	Use By:	Consult instructions for use
				
Batch Code	Catalog number	Manufacturer	Sterilized using aseptic processing techniques	Read Safety Data Sheet

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Important Licensing Information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

For additional technical information such as Safety Data Sheets (SDS), Certificates of Analysis, visit www.lifetechnologies.com/support. For further assistance, email techsupport@lifetech.com

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

DISCLAIMER: LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

www.lifetechnologies.com

