

Transferrin, Bovine, Holo Form

Description

Transferrin has been demonstrated to be an essential requirement for most cell lines grown in serum-free or serum-reduced conditions. The iron saturated (holo) form is acknowledged as the most efficacious in promoting cell growth in a variety of applications and formulations. Bovine transferrin (MW 74 kDa and 78 kDa) provides advantages over the human product as it is readily available and alleviates concerns regarding contamination by human adventitious agents.

Product	Catalog no.*	Amount	Storage
Transferrin, Bovine, Holo Form	11107-018	100 mg	2°C to 8°C
	11107-026	1 g	
	11107-059	5 g	
	11107-067	10 g	
	11107-047	25 g	
	11107-075	100 g	

* Availability is country-specific.

Product use

Caution: For manufacturing, processing, or repacking.

Important information

- Rehydrate transferrin immediately before use.
- Concentrated stock solutions can be prepared ahead of time and stored in a sterile air tight container at 2°C to 8°C for up to 12 months.

Safety information

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

Use

Prepare a concentrated stock solution (0.5–2.0 mg/mL) by dissolving transferrin powder in Phosphate-Buffered Saline (PBS) or other suitable solvent.

1. Add 100 mg transferrin to 45 mL room temperature PBS, mix thoroughly.
2. Rinse vial with an additional 2 mL PBS and combine with previous 45 mL volume, mix thoroughly.
3. Bring total volume to 50 mL with room temperature PBS, mix thoroughly.
4. Dilute to desired concentration (if less than 2 mg/mL) with PBS.
5. Filter sterilize through a 0.2-µm membrane into a sterile container prior to use. The filtered solution may be stored at 2°C to 8°C in a sterile airtight container for up to 12 months.







Transferrin has been demonstrated to be effective in the range of 0.5–100 µg/mL medium in promoting the growth of hybridomas, adherent and suspension cell cultures. We recommend performing a titration response curve to determine the optimal concentration of transferrin for each cell line and culture system.

Related products

Product	Catalog no.
Phosphate Buffered Saline, pH 7.4 (1X)	10010
Dulbecco's Phosphate Buffered Saline (1X), liquid	14040
Transferrin, Porcine, Holo Form	PSB10045

Explanation of symbols and warnings

The symbols present on the product label are explained below:

		
Manufacturer	Catalog number	Batch code
		
Caution, consult accompanying documents	Consult instructions for use	Temperature Limitation

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

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2014 Thermo Fisher Scientific Inc. All rights reserved.

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.

lifetechnologies.com

