

Anti-Clumping Agent

Description

Anti-Clumping Agent is a concentrated liquid reagent supplement designed to reduce cell aggregation or “clumping” in suspension cell culture without the use of enzymes. This reagent is an animal origin-free, chemically-defined formulation containing no proteins or hydrolysates. Cell lines adapted to suspension culture often can revert to a more adherent or aggregated clumping state. Anti-Clumping Agent has been evaluated for use in a variety of serum-free culture media for mammalian cell lines, including Chinese Hamster Ovary (CHO) and 293 cells. The addition of Anti-Clumping Agent has been shown to significantly reduce cell clumping (see Figure 1).

Product	Catalog no.	Amount	Storage	Shelf life*
Anti-Clumping Agent	01-0057AE	20 mL	2°C to 8°C	24 months
	01-0057DG	100 mL		

*Shelf life duration is determined from Date of Manufacture.

Product use

Caution: For manufacturing, processing, or repacking.

Important information

- Anti-Clumping Agent interferes with cationic lipid transfection and other transient transfection reagents, such as polyethylenimine (PEI). Do not use Anti-Clumping Agent until after transfection.
- Anti-Clumping Agent may also interfere with endotoxin testing using Limulus amoebocyte lysate (LAL).
- Not intended for use to dissociate adherent cells. If cell dissociation is desired, consider using TrypLE™ Select.

Safety information

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

Use

Anti-Clumping Agent is a reagent supplement designed to improve suspension cell culture condition by significantly reducing cell aggregation or “clumping”. Anti-Clumping Agent has been tested in a variety of cell lines grown in suspension culture.

The amount necessary to reduce cell clumping has typically been titrated at a dilution ranging from 1:100 to 1:1000. These dilutions are guidelines; the appropriate dose response for this reagent needs to be established for each application.

Anti-Clumping Agent should be added aseptically to cell culture media at the determined dilution prior to cell inoculation.

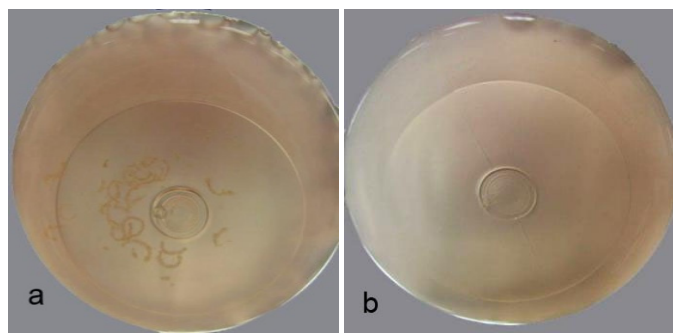


Figure 1









Suspension cell culture shake flasks with: (a) significant cell clumping, (b) very low level of cell clumping using Anti-Clumping Agent.

Related products

Product	Catalog no.
Freedom™ CHO-S™ Kit	A13696
CD FortiCHO™ Medium	A11483
CD CHO Medium	10743
TrypLE™ Select	12563

Explanation of symbols and warnings

The symbols present on the product label are explained below:

			
Use By:	Manufacturer	Batch code	Temperature Limitation
			
Catalog number	Consult instructions for use	Caution, consult accompanying documents	Sterilized using aseptic processing techniques

Limited Use Label License: Internal Research and Bioproduction Use

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product, under intellectual property rights that are owned and/or controlled by Life Technologies and relate specifically to the product, to perform (a) internal research for the sole benefit of the purchaser; and (b) research or manufacturing services conducted by the purchaser on a fee for service or contract basis for or on behalf of third parties. The purchase of this product does not grant the purchaser any additional rights, including (without limitation) the right to transfer or resell the product in any form, the right to use the product as a therapeutic agent or diagnostics test component, or to use the product to perform tests other than what is indicated in this Limited Use Label License on a contract or fee per test basis for or on behalf of third parties. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

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.

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.

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

