

# CTS™ DynaCollect™ Magnetic Separation System

Catalog Number A55867

Pub. No. 100117620 Rev. A

**Note:** For safety and biohazard guidelines, see the “Safety” appendix in the *CTS™ DynaCollect™ Magnetic Separation System User Guide* (Pub. No. MAN0026480). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Product description

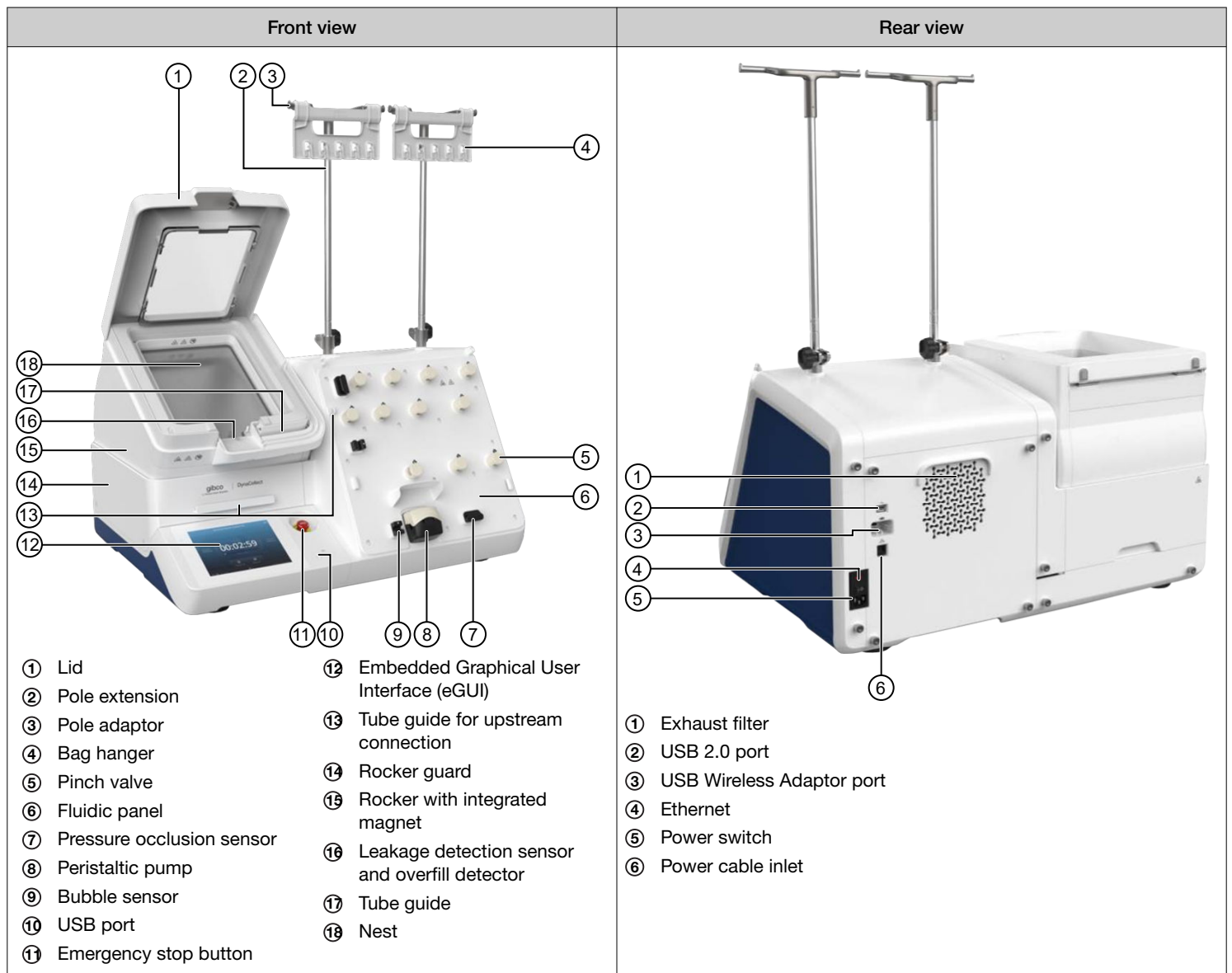
The Gibco™ CTS™ DynaCollect™ Magnetic Separation System is an instrument for automated, magnetic cell separation and bead removal. The instrument is accommodated by single-use fit-for-purpose kits for cell isolation and bead removal.

Two types of Single-Use Kits are used for performing cell isolation and bead removal.








- The CTS™ DynaCollect™ Cell Isolation Kit has 8 weldable input/output tubes for flexible setup configurations and an integrated isolation bag with maximum capacity of 1 L.
- The CTS™ DynaCollect™ Bead Removal Kit has 4 weldable input/output tubes for flexible setup configurations and an integrated bead removal bag for continuous processing.






# CTS™ DynaCollect™ Magnetic Separation System





## User interface overview

Symbol	Function
<b>Run Protocol</b>	
	Open and run protocol
<b>Create Protocol</b>	
	<ul style="list-style-type: none"> <li>• Create a new protocol</li> <li>• Create a new protocol from existing protocols</li> </ul>
<b>Manage protocol</b>	
	<ul style="list-style-type: none"> <li>• Edit existing protocols</li> <li>• Import protocols</li> <li>• Export protocols</li> <li>• Delete protocols</li> </ul>
<b>Unlock rocker lid</b>	
	Unlock rocker lid
<b>Rocker angle</b>	
	Adjust the rocker angle
<b>Open/close pinch valves</b>	
	Open and close pinch valves
<b>Manual mode</b>	
	Manual mode for executing single steps.

## Touch screen controls

Button	Function
	Returns to the previous screen
	Close the current window
	Go to menu screen

## Status indicators

Button	Function
	Indicates whether a USB device is inserted into the instrument
	Indicates whether the instrument is connected to wired network
Guest	Indicates user

## Instrument parameters

The CTS™ DynaCollect™ Magnetic Separation System is designed to operate within specific parameters. The values and limits for each parameter are listed below. If your input value exceeds the maximum value, an error is displayed.

Parameters that can be modified include:

- **Magnet:** Engaged/disengaged
- **Pinch valves:** Open/closed
- **Pump speed:** 10–150 mL/min
- **Pump:** Specified volume or duration
- **Bubble sensor:** Stop when wet/stop when dry
- **Prompt:** Active/inactive
- **Rocker:** Static, single-movement or continuous
- **Rocker angle:** -30° to 30°
- **Rocker speed:** 1–30 CPM (cycles per minute)

## Run a protocol on CTS™ DynaCollect™ Magnetic Separation System

1. Create isolation or bead removal protocol or use pre-existing template protocols.
2. Prepare cells and reagent bags according to the configuration for the protocol.
3. Unpack Single-Use Kit.
4. Weld reagent bags to Single-Use Kit.
5. To start a run, open the isolation or bead removal protocol (choose a customer created protocol or template protocol) and use the “How to install” instructions on the eGUI to install the Single-Use Kit to the instrument.
6. Select **Start run** to initiate the protocol.
7. When the protocol is complete, seal off the output bag containing the product and follow the instructions on the eGUI to remove the Single-Use Kit from the instrument.

## 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 at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

Revision history: Pub. No. 100117620

Revision	Date	Description
A	13 October 2022	New document for CTS™ DynaCollect™ Magnetic Separation System.

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.