

CTS™ DynaCollect™ Magnetic Separation System

Catalog Number A55867

Pub. No. 100117621 Rev. A

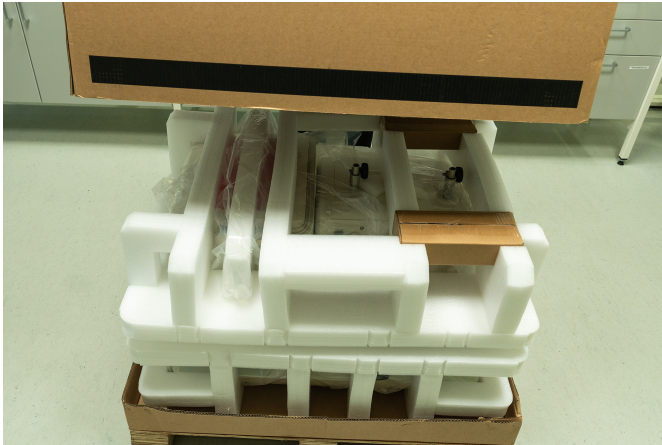
Note: For safety and biohazard guidelines, see the “Safety” appendix in the *CTS™ DynaCollect™ Magnetic Separation System User Guide* (Pub. No. MAN0026480). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Important guidelines

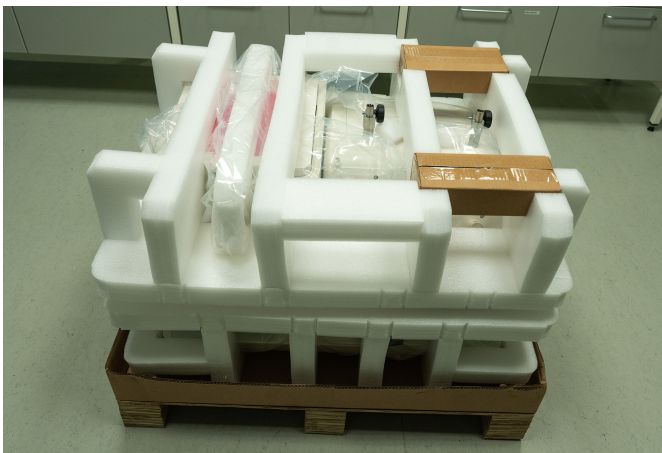
- The working area must be at least 749.5 mm (29.5") (width) × 562 mm (22.12") (depth) (include 10 cm rear clearance for ventilation) × 495 mm (19") (with pole base) (height), flat, dry, clean, and vibration-proof.
- Ambient humidity should be no more than 80% at temperatures up to 30°C (86°F) to avoid condensation.
- Ambient temperature range should be between +15°C (59°F) and +30°C (86°F).

Unpack and set up the instrument

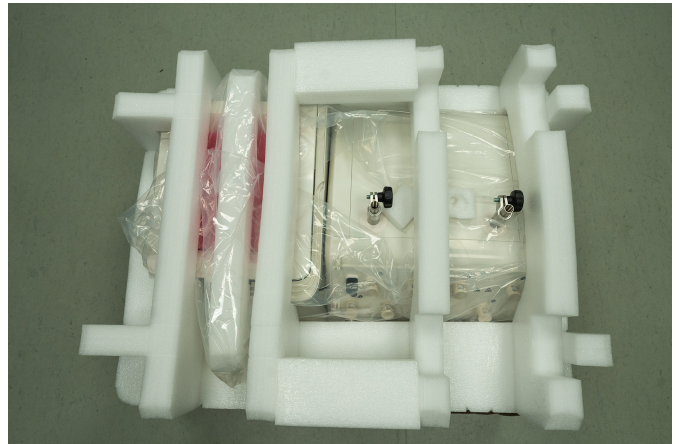
1. Lift the cardboard box off the pallet.



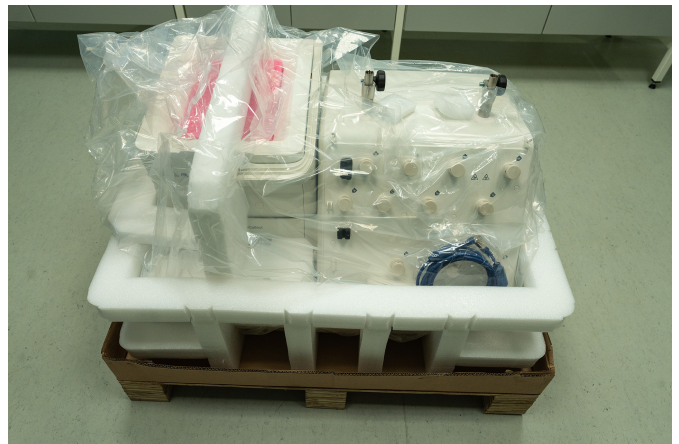
2. Remove the two small cardboard boxes containing the bag hangers, the pole extensions and the pole adaptors.



3. Remove the top foam.



4. Unwrap packaging bag and remove the ethernet cable and power cable.

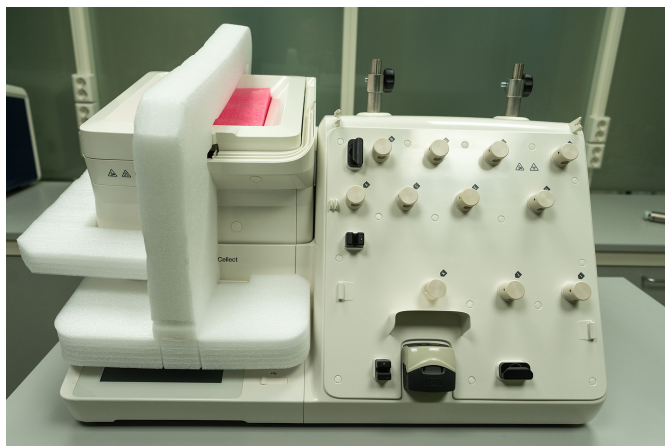


- Lift the instrument to bench or cart. The operation requires two people. Use the four holding recesses (front left side, back right side, and on both sides) to lift in a diagonal position.

IMPORTANT! The instrument weighs 50 kg (109.7 lbs) and requires two people to lift.



- Place the instrument in a protected location with at least 10 cm (4 inches) to the rear for proper operation of the exhaust fans. See "Important guidelines" on page 1.
- Remove the rocker homing position foam.

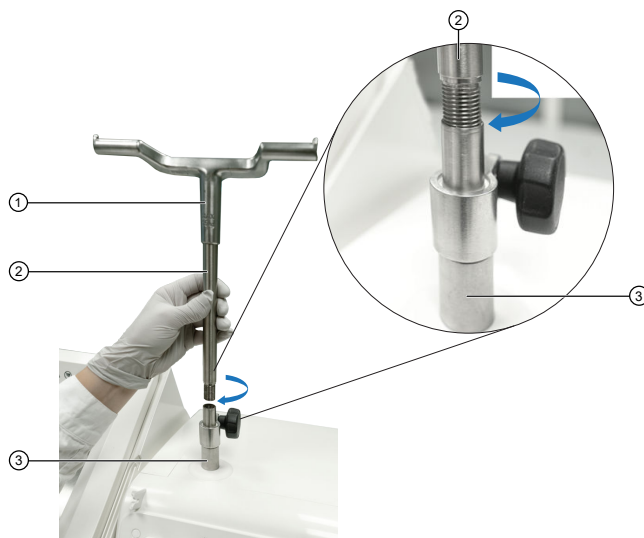


Note: It is acceptable for the rocker to freefall by gravity after removing the foam. Do not manually move the rocker once it has settled into its locked resting position.

Contents

Component	Quantity
CTS™ DynaCollect™ Magnetic Separation System (Cat. No. A55867)	1
Region specific power cord (for US/Canada/Taiwan/Japan, Europe, or UK)	1
Ethernet cable	1
Pole extension	2
Pole adaptor	2
Bag hanger	2
Quick Reference Guide	1
Installation Guide	1

- Install the two pole adaptors to the two pole extensions, and install both on the pole base on the instrument.



- Pole adaptor
- Pole extension
- Pole base

- Place the bag hangers on the pole adaptors.
- Connect the power cable to the power inlet, and insert the plug into an electrical outlet.
- (Optional) Connect one end of a ethernet cable to the instrument ethernet port, and the other end to an ethernet port wall plug.

Working space

The CTS™ DynaCollect™ Magnetic Separation System is designed for benchtop use and requires a working space of approximately height 495 mm (19") with pole base and 434 mm (17") without pole base, width 749.5 mm (29.5"), depth 562 mm (22.12"). The system can also be placed on a cart.



Product specifications

Category	Property	Value
Electrical		
	Supply voltage	100–240 VAC ±10%, 50/60 Hz, 600 W
	Phases	Single
	Maximum rated input current	6 A
	Fuses	2 × 10 A
Instrument		
	Regulatory requirements	HEPA filter meets ISO 14644-14
	Installation category	II
	Instrument type	Benchtop/cart unit
	Instrument dimensions	495 mm (19") with pole base/ 434 mm (17") without pole base (H) × 749.5 mm (29.5") (W) × 562 mm (22.12") (D)
	Instrument weight	49.75 kg (109.7 lbs) (50.6 kg/111.6 lbs with poles and bag hangers)
	Built-in features	8" touch screen display (800 × 600 px)
Sound level		
	Operating sound level	73 dBA (measured 1 m from instrument)
	Non-operating sound level	63 dBA (measured 1 m from instrument)
Environmental ranges		
	Ambient temp	15°C to 30°C
	Transport temperature	-30°C to 60°C (extreme end temperatures at maximum 72 hours)
	Storage temperature	15°C to 30°C
	Relative humidity	15%–80% (non-condensing)
Operating limits		
	Magnet	Engaged/Disengaged
	Pinch valves	Open/Closed
	Pump speed	10–150 mL/min
	Rocker angle	-30° to 30°
	Rocker speed	1-30 CPM (cycles per minute)
	Pressure occlusion sensor	Max 2 Bar. Only detect liquid pressure (not pressure build up by air). Following lines are not covered: <ul style="list-style-type: none"> • CTS™ DynaCollect™ Cell Isolation Kit – Ports B and K • CTS™ DynaCollect™ Bead Removal Kit – Ports B, D and E
	Height rocker lid (isolation)	35 mm (1.4")
	Height rocker lid (bead removal)	5 mm (0.2")
	Overfilled detection	1120 mL (static and rocking)
Magnet		
	Magnet dimensions	265 × 205 mm (10.4" × 8.07")
	Magnetic strength	≥ 6000 G
Pole		
	Max height	1070 mm (42.13") from the bottom of the rubber foot of the instrument to the tip of the bag hanger hook point

Category	Property	Value
Pole adaptor		
	Max weight	3 kg (6.6 lbs) per hook on bag hanger. Total 15 kg (33 lbs) per bag hanger
Use		
	Indoor use only	
	Not intended for use in a wet environment	
	Intended for use in Pollution degree 2 environment	

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. If you have any questions, please contact Life Technologies at 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.

Revision history: Pub. No. 100117621A

Revision	Date	Description
A	17 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.