

Important guidelines

- The working area must be 67.4cm (width) × 64 cm (depth including 10 cm rear clearance for ventilation) × 110 cm (front lid fully open), 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).

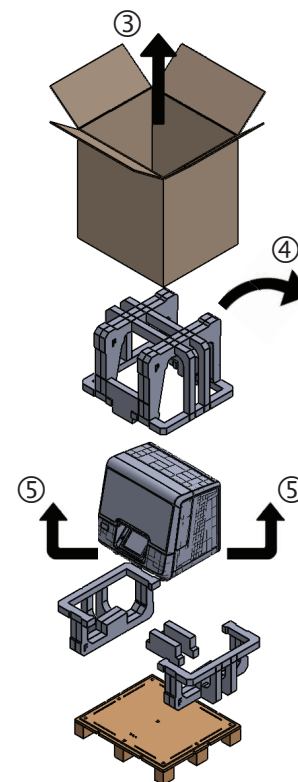
Unpack and set up the instrument

1. Remove the plastic covering from around the box.
2. Cut the packing straps to release the box from the pallet.
3. Open the top of the box and lift the box from the pallet.
4. Remove the upper packaging foam and the plastic bag covering the instrument.
5. Lift the instrument from the box using the grab rails underneath the sides of the instrument, and place it on a benchtop.

Important! The instrument weighs 70 kg (154 lbs) and requires at least two people to lift.

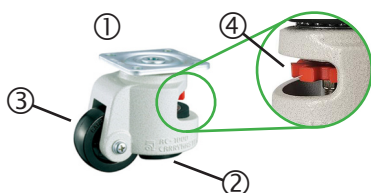


6. Place the instrument in a protected location with at least 10 cm (4 inches) to the rear for proper operation of the inlet and exhaust fans (see Important guidelines for additional details).
7. Open the door of the instrument and remove the two black foam strips from the rear edges of the door, then remove the styrofoam insert from the electroporation chamber holder.
8. Connect the power supply cable to the power inlet, and insert the plug into an electrical outlet.
9. (Optional) Connect one end of a Ethernet cable to the instrument Ethernet port, and the other end to an Ethernet port wall plug.



Level the instrument

1. Ensure that it is resting on the flat rubber foot of each castor, and not the wheels to prevent the instrument from rolling.



- ① Castor
- ② Rubber foot
- ③ Wheel
- ④ Adjustment knob

2. (Optional) If adjustment of the castors is required, use the bubble level to ensure the level placement of the instrument.

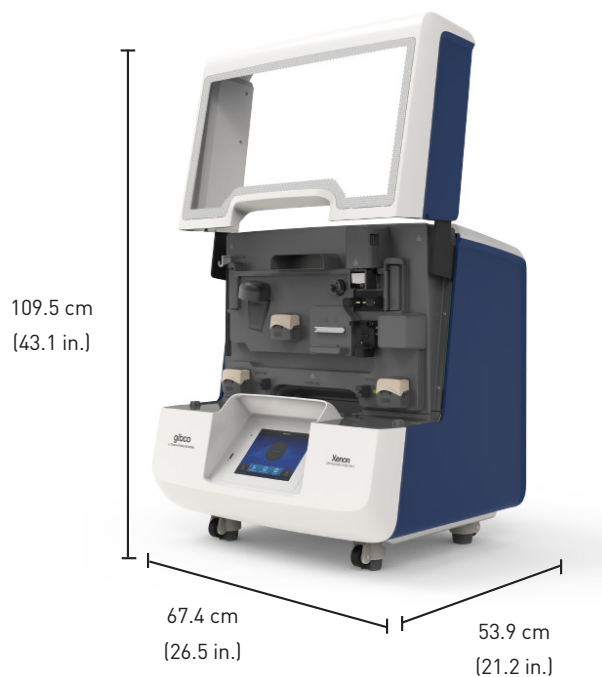
Place the bubble level in the electroporation chamber holder, then adjust the height of the castors by turning the red adjustment knob(s) by hand until the bubble is within the red circle.



- ① Electroporation chamber holder
- ② Bubble level

Contents

Components	Cat. No.
	A52727
CTS™ Xenon™ Electroporation Instrument	1
Region specific power cord (for US/Canada/Taiwan/Japan, Europe, or UK)	1
Ethernet cable	1
USB memory device (containing user guide)	1
Quick reference	1
Install guide	1
Bubble level	1



Specifications

Parameter	Value
Electrical rating	100–240 ± 10% VAC, 50/60 Hz, 1200 VA
Electroporation pulse voltage range	500–2500 V
Electroporation pulse width range	1–30 ms
Electroporation pulse interval range	500–1000 ms
Number of electroporation pulses	1–10 pulses
Cell mixer stirrer rotating speed	60 rpm
Pre-cooling technology	Peltier
Pre-cooling temperature setting range	10–30°C
Electroporation chamber volume	1 mL
Operating Temperature	15°C to 30°C
Maximum Relative Humidity	Up to 80%
Degree of Protection	IPX0
Protective Earthing	Class I (earthed)
Installation category	II
Instrument type	Benchtop unit
Instrument dimensions (door fully open)	674 mm (w) × 539 mm (l) × 1095 mm (h)
Instrument weight	70 kg (154 pounds)

For support visit thermofisher.com/support
thermofisher.com

28 September 2021

ThermoFisher
 SCIENTIFIC