

# Smart Notes



## How can the new design of Thermo Scientific General Purpose Pro Centrifuges improve its ergonomics and ease-of-use?

After careful consideration of the voice of the customer, Thermo Scientific™ General Purpose Pro Centrifuges feature a new, modern design that has been extensively tested by customers and provides improved ease-of-use and safety during operation.

### Safe operation

Improved reach-in during rotor installation for safer loading and unloading

- Loading height decrease 1.2 cm (0.5in)
- Lid grapping height decreased 5.1 cm (2 in)
- Reduced closing force of the lid
- Hand grip cue on the lid indicates the easiest and best one-handed closure location, which can also extend the life of your lid gas spring.
- Improved finger squeeze situation: New design fully adheres to finger squeeze specifications and requirements.

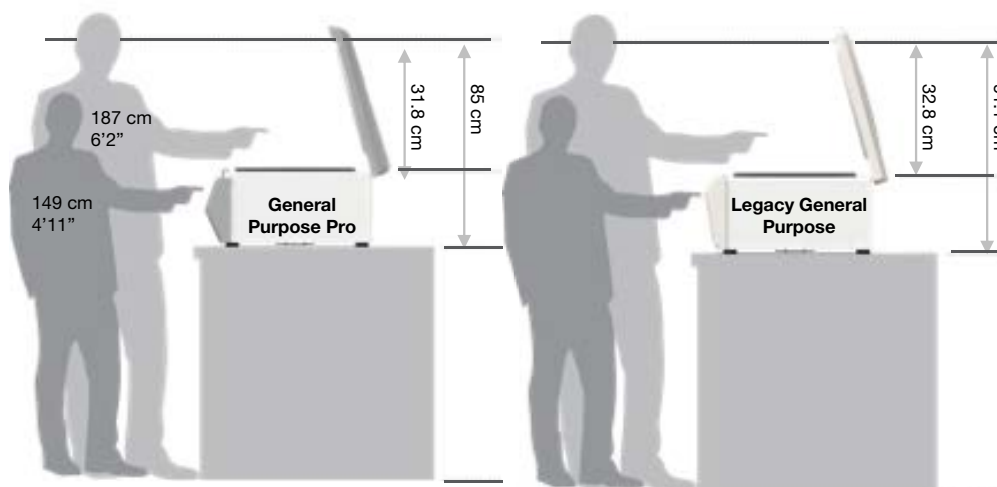
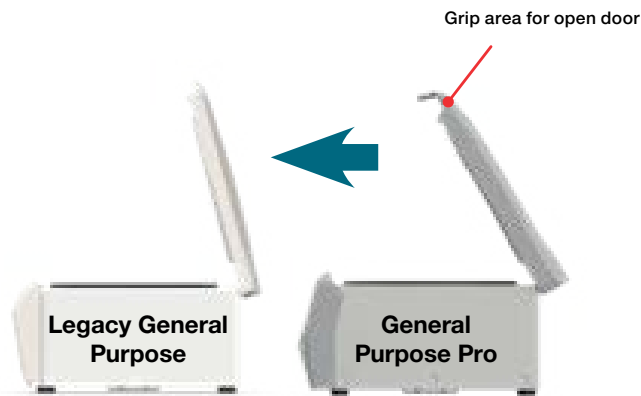


Figure 1: Improved reach-in and loading height decrease



**Figure 2:** New door shape and open angle provide a reach area closer to the user than the previous design. Grip area for open door position is lower and closer to the user.

## General Purpose Pro Centrifuges feature multiple ergonomic and safety features that make their operation simpler and safer.

Enhancements to centrifuge loading and unloading provides safer use and accessibility during reach-in.

Additional enhancements and new benefits include:

- Accessory tray that holds tools or samples as you prepare for your run
- Front USB port for easier service technician plug in
- Compliance with most-recent applicable regulatory and safety standards
- Available energy-savings settings
- Over 19 rotor options to choose from
- Better wear-and-tear of the instrument by changing the centrifuge body surface to one that attracts less dirt

### Summary

Following usability studies, multiple ergonomics features have been added to the NEW! Thermo Scientific General Purpose Pro Centrifuges to simplify their use.



Small recess to hold items while using centrifuge



Experience General Purpose Pro Centrifuges at [thermofisher.com/gp-pro-centrifuge](https://thermofisher.com/gp-pro-centrifuge)

**ThermoFisher**  
SCIENTIFIC