

Choosing the right water bath for your requirements

I want to keep my samples at a temperature of 37°C. **Which type of water bath is ideal for this requirement: general purpose, circulating, or shaking?**

General purpose, circulating, and shaking heated water baths are all designed to keep samples at your desired temperature. So, the ideal type of bath depends on whether your samples need to be thawed, need shaking to keep them emulsified, or need circulation to meet a uniformity requirement.

Water baths are used to process a large variety of samples in a number of applications, including chemical, biological, and microbiological research, as well as in production laboratories of the food and pharmaceutical industries. Certain applications have specific requirements, whether it is temperature uniformity or sample uniformity. It is important to understand your unique requirements in order to select your ideal type of bath.



What are the differences in the various types of water baths?

Precision redefined for your lab

General purpose water baths

Thermo Scientific™ Precision™ general purpose water baths range between 2 and 28 liters to meet a variety of volume requirements. Operational components such as heaters and sensors are attached to the underside of the bath chamber, allowing full use and easy cleaning of the entire bath area. Precision general purpose water baths are designed to minimize noise and maintenance.



Circulating water baths

Thermo Scientific™ Precision™ circulating water baths are designed to provide greater temperature uniformity compared to general purpose water baths, making them more suited for samples that require precise temperature control. Precision circulating water baths use an efficient heating element combined with a magnetically coupled circulating pump, resulting in outstanding temperature uniformity.

Shaking water baths

Thermo Scientific™ Precision™ shaking water baths are designed to prevent separation through continual mixing for outstanding sample uniformity. Precision shaking water baths feature a next-generation shaker motor and pump, as well as presets for the oscillation speed for your convenience.

Summary

The portfolio of Precision water baths is designed to simplify workflows and maximize productivity, with features like intuitive graphical displays, audible timers, and temperature presets. Precision baths also have global voltage for easy ordering.

Learn more about Precision water baths at
thermofisher.com/precisionbaths

ThermoFisher
SCIENTIFIC