

Orbit mini. Temperature Control.

- Active heating and cooling of whole bilayer chamber
- Wide temperature range of 5 - 50 °C applicable
- Automated temperature control throughout whole experiment
- Experiments at physiological temperature

The **Orbit mini** temperature control add-on

The optional temperature control add-on for Nanion's Orbit mini bilayer workstation allows for the convenient application and automated control of the temperature while performing bilayer experiments. The desired temperature can simply be defined via the Orbit mini's control software and is automatically kept constant thanks to various sensors inside the device. Active liquid cooling ensures precisely constant temperatures even during long-term experiments.

Key features of the **Orbit mini** temperature control

- Temperature freely definable between 5 - 50 °C
- Convenient control via Orbit mini's recording software
- Small footprint (35 x 22 x 11 cm) to be usable with the Orbit mini device anywhere
- Powerful 10 W Peltier element to actively heat up and cool down the whole bilayer chamber
- Liquid cooling of Peltier element to ensure stable temperature controlled operation for hours



Precise control of the temperature during a bilayer experiment is not only advantageous for the investigation of temperature-sensitive species like TRP channels. Conducting experiments at lowered temperatures enables the active alteration of kinetics and therefore e.g. the assessment of events otherwise too short to be detected. Heating of the chamber allows for experiments at physiological temperatures and the consideration of lipid phase transition temperatures.

Orbit mini

nanjion