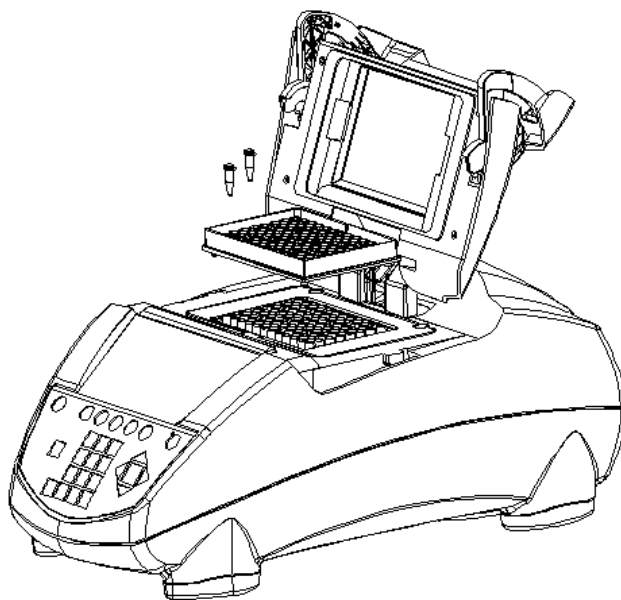


# MyCycler™ Sample Loading Tray for the MyCycler Thermal Cycler

Catalog Number 170-9709

The MyCycler Sample Loading Tray is designed to function as an optimization tool for the MyCycler Thermal Cycler. When correctly positioned in the MyCycler reaction block, the tray evenly distributes lid pressure across samples loaded into the MyCycler instrument. The array of openings in the MyCycler Sample Loading Tray should be positioned directly over the 96-wells of the sample block area, allowing the sample tubes to be placed directly into wells of the reaction block.



The MyCycler Sample Loading Tray is compatible with individual 0.2 ml thin wall tubes with domed caps (such as Bio-Rad 0.2 ml Thin-Wall Tubes, catalog number 223-9473). Once correctly positioned in the MyCycler instrument, the Sample Loading Tray can be left in the reaction block at all times. The multifunctional tray can also be used to transport PCR tubes between the workbench and the reaction block, or stacked upon an empty tray to serve as a post-PCR storage system for 0.2 ml tubes.

Use of the MyCycler Sample Loading Tray is strongly recommended with individual 0.2 ml thin wall tubes. The tray ensures uniform lid pressure across the sample block area, and therefore provides optimum contact between the heated lid and each PCR sample.