

iCycler™ Plate Removal Tool for iCycler 384 Well Reaction Module 170-8735

The iCycler Plate Removal Tool facilitates removal of 384 well reaction plates from the Reaction Module.

The small reaction volumes typical of 384 well formats require optimal contact between reaction plate and reaction module. The use of the iCycler Amber Pad is strongly recommended to attain a tight fit of the sample plate in the Reaction Module.

During the reaction, the heating and cooling cycles further secure a tight fit of the sample plate in the Reaction Module so that at the end of the reaction, it can be difficult to directly remove the plate from the iCycler.

The Plate Removal tool should be positioned with the hook shaped edge under the sample plate. Lift the inserted tool to release the sample plate from the Reaction Module.

