

iCycler™ Sealing Ring for iCycler 96 Well Reaction Module

170-8731

The iCycler Sealing Ring optimizes temperature isolation of the sample reaction block area. This sealing ring is intended for use with the iCycler 96 well 0.2ml reaction module. This sealing ring is green, thermal isolation rings of other colors are to be used with other reaction module formats for the iCycler Thermal Cycler.

The Thermal Sealing Ring also offers structural support minimizing the number of tubes required in any experiment. Note that the thermal isolation ring has a slot for use with the sample probe. When using the sample probe for temperature control, position the sample probe to feed through this slot.

The green ring should be positioned directly over the ridge surrounding the sample block area with the Bio-Rad logo towards the back of the instrument.

Although the Sealing Ring is not required for operation, use of the sealing ring reduces condensation buildup sometimes observed after long time (24+ hours) at 4 °C hold temperatures.

