

# iCycler™ Sealing Ring for iCycler 60 Well Reaction Module

## 170-8732

The iCycler Sealing Ring optimizes temperature isolation of the sample reaction block area. This BLUE sealing ring is intended for use with the iCycler 60 well 0.5 ml reaction module. Thermal isolation rings of other colors are to be used with other reaction module formats for the iCycler Thermal Cycler.

The BLUE sealing ring should be positioned directly over the fixed ridge surrounding the sample block area with the Bio-Rad logo towards the back of the instrument and the slotted flange engaged around pin, located near probe connection, and under screw head.

To properly engage sealing ring, tilt ring logo upwards and slide slotted flange under polished screw head and around pin, then tilt ring logo downwards to engage around fixed ridge surrounding sample block area. To verify installation, push down on ring and then release; ring should spring back up to original position.

Note that the thermal isolation ring also 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.

This ring is required for all 60 x 0.5 ml thermal cycling operations. The use of the sealing ring optimizes thermal uniformity and reduces condensation buildup sometimes observed after long term (24+ hours) conditions at 4 °C hold temperatures.

If you have any difficulty with this procedure or need additional information, please contact your local Bio-Rad office or in the U.S., call technical services at 1-800-BIORAD (1-800-424-6723).

