

iCycler™ Amber Pad for iCycler 384 Well Reaction Module

170-8739

The iCycler Amber Pad optimizes both lid pressure and temperature isolation of samples for the 384 well Reaction Module.

The small reaction volumes typical of 384 well formats require optimal contact between reaction plate and reaction module. The use of the Amber Pad ensures that a tight fit of plate to reaction module can be attained.

The Amber Pad also serves to isolate the liquid sample from the heated lid so that accurate sample temperature may be achieved. Small reaction volumes may be adversely affected by the heat generated at the lid. The Amber Pad prevents any temperature effect due to the heated lid.

The Amber Pad should be positioned over the sealed 384 well reaction plate just prior to sliding the reaction module lid to the closed position. This pad does not serve as a replacement for plate sealing. The Amber Pad is strongly recommended for all reactions in the 384 well Reaction Module.

Position Amber Pad
directly over sealed
reaction plate

