

Life Technologies recommends that the following 0.5 mL microcentrifuge tube be used to mate with the Arcturus[®] CapSure[®] and ExtracSure[™] Device assembly during incubation:

GeneAmp[®] Tube (#N8010611)
Applied Biosystems