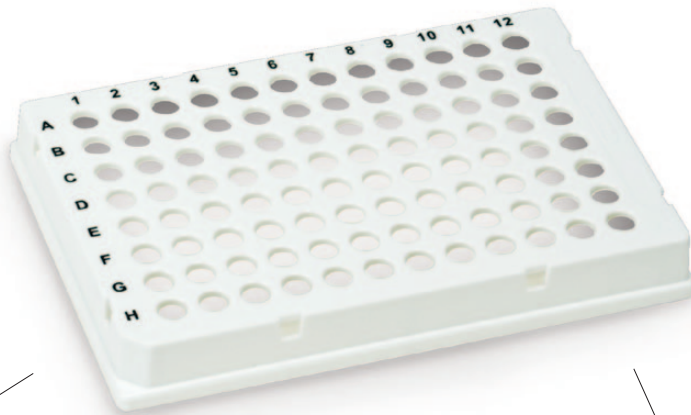


AMPLIFICATION PCR Tube Rack

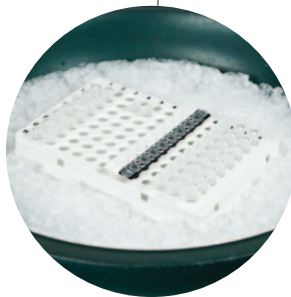
- Holds 0.2 ml and 0.1 ml PCR tubes and unskirted and semi-skirted PCR plates
- Has ANSI/SBS standard footprint for robotic handling
- Prevents tubes from sticking to rack
- Allows organization of PCR tubes on ice
- Supports tubes in centrifuges and compatible thermal cyclers
- Is autoclavable and affordable

PCR Tube Rack — New Design!

Secure your PCR tubes and plates on ice, robots, centrifuges, and thermal cyclers.



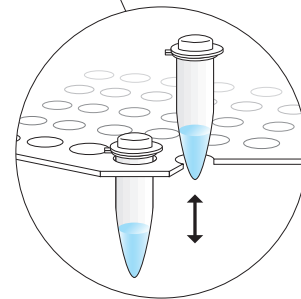
Smooth robotic handling
(ANSI/SBS standard)



Easy organization on ice



Support in centrifuges and thermal cyclers



No more sticking



Other Capping Tools and Racks

96-Place PCR Tube Rack and Cover

These stackable storage units for tubes and PCR plates provide a stable platform for preparing or centrifuging reactions.

Easy Cap™ Tool

The Easy Cap tool provides the pressure necessary to achieve a tight seal, one tube at a time, when capping individual or strip tubes. The narrow end securely fastens domed caps to 0.2 ml tubes or 96-well PCR plates. The wide end firmly holds thin-wall 0.5 ml tubes to prevent accidental crushing when opening or closing. The side slot allows easy opening of tight-fitting caps without generating aerosols.

Strip Cap Tool

This tool quickly and easily seats 8- and 12-cap strips on PCR plates or tubes. A grooved channel on one side is designed to seal domed caps, while the flat edge on the opposite side seats flat caps. For best results, seal tube strips while they are in a thermal cycler block or in a 96-place rack.

Ordering Information

Catalog #	Description
TRC-9601	PCR Tube Rack , ANSI/SBS standard, white, 10
TRC-0501	96-Place Racks , with covers, for PCR tubes and unskirted and semi-skirted microplates, assorted colors, 5
ECT-1000	Easy Cap Tool , ensures tight seal for 0.2 ml PCR tubes or 96-well microplates
ECT-2000	Strip Cap Tool , for seating 8- and 12-cap strips on PCR plates or tubes

Practice of the polymerase chain reaction (PCR) may require a license.



Strip Cap Tool



96-Place Racks and Covers

Easy Cap Tool

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600
Canada 905 364 3435 China 86 20 8732 2339 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000