

microarray systems

The VersArray™ Colony Picker and Arrayer Systems

Revolutionary Robotic Systems With Versatile Capabilities

VersArray colony arrayer and VersArray colony picker and arrayer systems are revolutionary robotic systems that let you easily and efficiently perform colony picking, arraying, gridding, or replicating of high-density plates of colonies or cultures. Features include:

- 96-, 384-, and 768-pin replicator heads
- Optional 4-level elevator for high-throughput capacity
- Ability to transfer samples from:
 - Gel to gel (replication)
 - Plate to plate (liquid transfer)
 - Plate to gel (arraying)
 - Plate to membrane (gridding or microarraying)
 - Gel to plate (colony picking)



BIO-RAD



The VersArray Colony Picker and Arrayer Systems

Colony Arrayer: Unique Colony Arraying and Gridding Capability

The VersArray colony arrayer forms an ordered matrix of colonies on a gel plate that can then be screened for the effect of virtually any genetic or chemical interaction.

Colony Picker and Arrayer: Added Versatility With Colony-Picking Functions

The colony picker and arrayer uses a sophisticated image acquisition system to prioritize, pick, and inoculate over 9,000 colonies per hour based on colony size or blue/white selection with unattended operation. It can also perform colony arraying and gridding.

Liquid Handling Module

Either system can be easily transformed to provide liquid handling capabilities. The four-channel liquid handling module uses disposable-tip pipets to replicate or rearrange 0.2–250 µl samples from and into 96- or 384-well plates. In addition, these are the only instruments of their kind to provide open source code for the robotics, allowing you unparalleled flexibility for your experiments. And like the VersArray™ ChipWriter Pro system, they are unmatched in reliability and are completely contained within an environmental chamber that protects your samples and materials.

Specifications

Superior resolution	1.22 µm (x-, y-axes), 0.24 µm (z-axis) with 3 µm repeatability
Arraying throughput	>23,000 colonies/hr (1 min wash cycle — variable)
Pin cleaning	Heat sterilizer and brush station for ultimate pin cleaning
Wash station	Pump (optional), sonicator, and air-drying unit
Replicator head types	96, 384, or 768 pins, 0.5 or 1.0 mm
Colony-picking head	High-speed head with 16 indexable pins
Gridding tools	96- or 384-pin head with 0.5 or 1 mm pins
Liquid handling tools	Two 4-channel disposable-tip pipets (0.2–10 µl and 10–250 µl)
Stacker module	Holds 32 shallow plates in 8-plate stacks
Environmental chamber	Fully enclosed, HEPA filtered with positive air pressure, humidity control
Robotic gripper arm	For plate or lid transfer
Dimensions (W x D x H)	127 x 102 x 143 cm (51 x 41 x 57")

Ordering Information

Catalog #	Description
169-0010	VersArray Colony Arrayer System, 110/120 V
169-0012	VersArray Colony Arrayer System, 220/240 V
169-0011	VersArray Colony Picker and Arrayer System, 110/120 V
169-0013	VersArray Colony Picker and Arrayer System, 220/240 V

For complete information on the full line of VersArray systems, contact your local Bio-Rad sales representative or visit us on the Web at www.bio-rad.com/versarray/

Certain arrays and/or the method of preparation, analysis, or use may be covered by patents or other intellectual property rights owned by third parties. Use of this product is authorized only for applications for which the end user has a license under proprietary rights of third parties or for applications for which a license is not required.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 2527 3454 Canada (905) 712-2771 China (86-21) 63052255 Czech Republic + 420 2 41 43 05 32 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84-0 Hong Kong 852-2789-3300 Hungary 36 1 455 8800 India (91-124)-6398112/113/114, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20 The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 331 99 99 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-6415 3188 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000