



*Certified RNase/DNase- and Pyrogen-Free  
for Large-Volume Pipetting up to 1 mL*

## BIOMEK® P1000 TIPS

**CONDUCTIVE, DISPOSABLE TIPS FOR  
BIOMEK NX & FX SPAN-8 SYSTEMS**

Biomek P1000 Tips are now available for the Biomek NX and FX Span-8 Systems. These new certified tips offer increased tip capacity for large-volume pipetting up to 1 mL, for such applications as:

- DNA (genomic and plasmid) purification from bacterial cell cultures, blood and related sources, body fluids, swabs, and animal and plant tissues
- Viral titer measurement from blood and related sources
- RNA isolation from bacterial cell cultures, blood and related sources, CSF, and animal and plant tissues
- Protein purification using HIS-tag chemistry
- Solid-phase extraction of target analytes using filtration plates or 1 mL SPE columns
- Broadcast dispensing of reagents for assay development and dilution studies

The P1000 conductive, disposable tips are **certified "RNase/DNase-free"** and offered in three configurations:

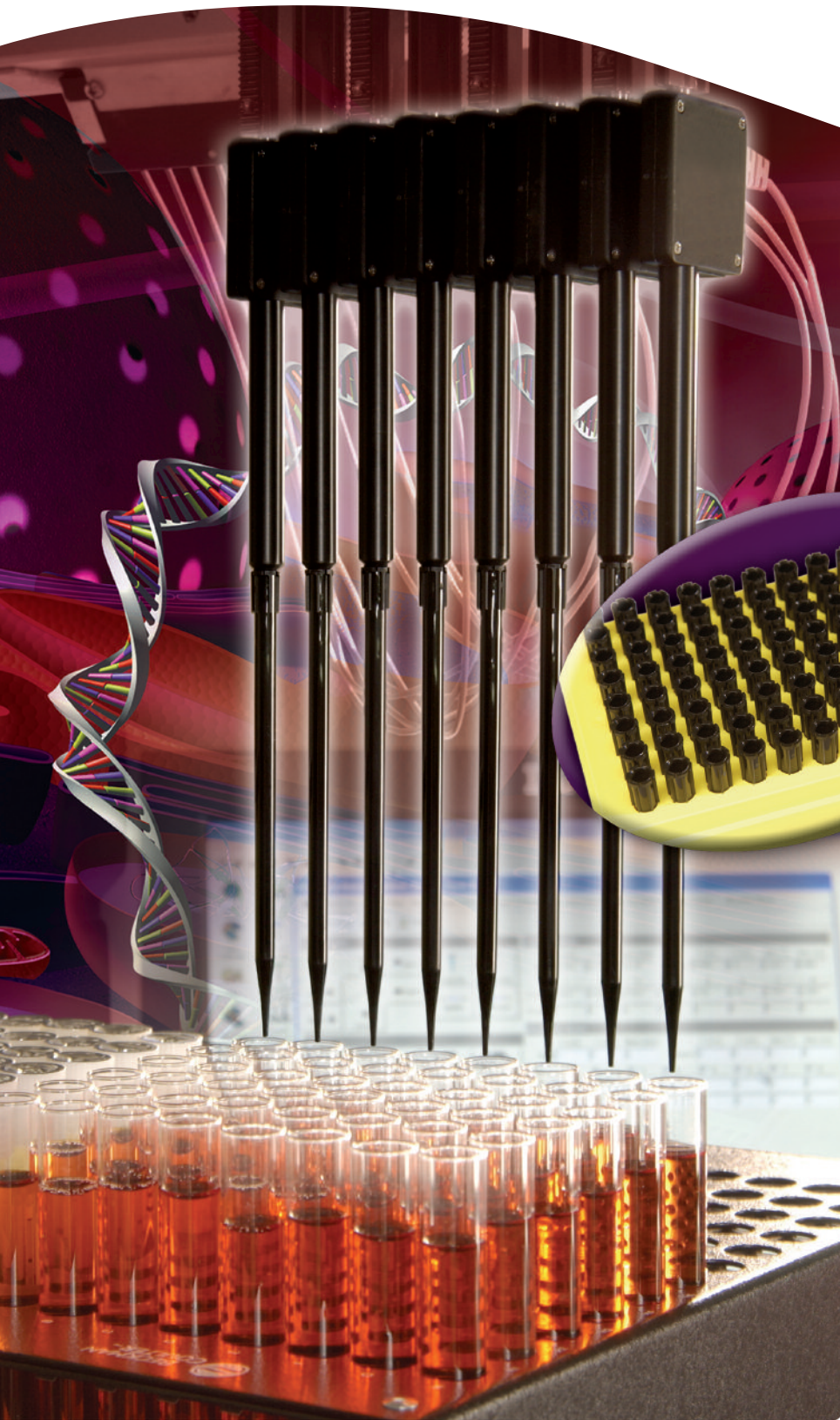
- nonsterile
- presterile
- presterile with barrier

Working capacities for the tips are:

- 1050  $\mu$ l (including air gaps) for P1000 tips without barriers
- 1025  $\mu$ L (including air gaps) for P1000 tips with barriers

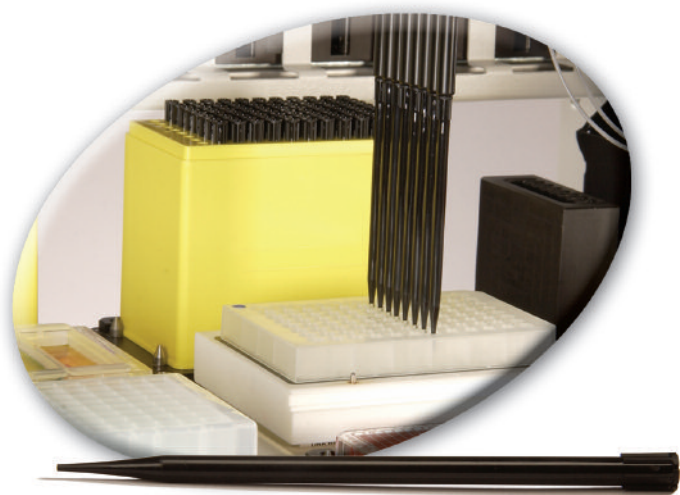
Compatible labware commonly used with P1000 tips includes:

- 1 and 2 ml, 96-well deepwell plates
- 96-well micro- and amplification plates
- 10, 12, and 13 mm diameter test tubes
- Microfuge® tubes
- 1 and 2 ml SPE filtration plates
- Reagent reservoirs



Use of the Biomek P1000 Tips requires the appropriate Biomek Software and Span-8 Disposal ALP (Automated Labware Positioner) hardware. Contact your local Beckman Coulter representative for upgrading Biomek NX and FX Systems.

With Beckman Coulter, you are assured that extensive verification and validation testing has been performed on the P1000 tip to assess both product and system level performance.



## CERTIFIED RNase/DNase-FREE

Biomek P1000 Tips are certified to be free of RNase, DNase, DNA (human and mouse)/PCR Inhibition, Pyrogen/ Endotoxin, and Trace Metals.

RNase- and DNase-free certification ensures high quality nucleic acid purification and reliable nucleic acid testing results.

DNA-free (human and mouse)/PCR® Inhibition certification ensures that amplification applications are free of contaminants that result in erroneous results and interfere with the PCR.

Pyrogen/Endotoxin-free certification addresses endotoxin-sensitive applications, such as cell transformations and preparative protocols, that use extracted biological material for in vivo experimentation.

Trace metal-free testing of the base, polypropylene tip resin reduces possible interferences due to chelation and denaturation.

## STORAGE CONDITIONS

Biomek Tips are precision plastic accessories intended for use on Biomek Laboratory Automation Workstations. Validation and performance of the systems are dependent on the proper storage of these precision components.

It is recommended that Biomek Tips are stored in a clean and dry laboratory environment at room temperature for no longer than two years. Exposure to high temperatures, radiation, or vaporous chemicals/solvents may result in the deterioration of the design of the tip, and therefore its performance.

## PRESTERILE PRODUCT

Presterile Biomek Tips with packaging intact and undamaged, and stored under recommended conditions is considered to retain its sterile nature indefinitely.

Due to the precise geometries and tolerances required of Biomek Tips in automated systems, autoclaving should not be used to provide sterility.

## ORDERING INFORMATION

Part No.	Description	Conductive	Working Capacity	No. of Tips/ Case	Package Quantity
987935	<b>Span-8 P1000 Tips, nonsterile</b>	Yes	1050 µL	480	96 tips/rack, 5 racks/case
987936	<b>Span-8 P1000 Tips, presterile</b>	Yes	1050 µL	480	96 tips/rack, 5 racks/case
987925	<b>Span-8 P1000 Tips, presterile with barrier</b>	Yes	1025 µL	480	96 tips/rack, 5 racks/case

**For the best in system performance, use only Biomek validated tips on Biomek instruments.**



*Developing innovative solutions in Systems Biology.*

**Innovate Automate**  
SIMPLIFY

Beckman Coulter, Inc. • 4300 N. Harbor Boulevard, Box 3100 • Fullerton, California 92834-3100  
Sales & Service: 1-800-742-2345 • Telex: 678413 • Fax: 1-800-643-4366 • www.beckmancoulter.com

### Worldwide Biomedical Research Division Offices:

**Australia** (61) 2 9844-6000 **Canada** (905) 819-1234 **China** (86) 10 6515 6028 **Eastern Europe, Middle East, Africa** (41) 22 994 07 07  
**France** 01 49 90 90 00 **Germany** 49 21 513335 **Hong Kong** (852) 2814 7431/2814 0481 **Italy** 02-953921 **Japan** 03-5404-8359  
**Mexico** 525-605-77-70 **Netherlands** 0297-230630 **Singapore** (65) 6339 3633 **South Africa** (27) 11-805-2014/5 **Spain** 91 3836080  
**Sweden** 08-564 85 900 **Switzerland** 0800 850 810 **Taiwan** (886) 2 2378 3456 **Turkey** 90 216 309 1900 **U.K.** 01494 441181 **U.S.A.** 1-800-742-2345