



Biomek Tips and Labware

The Most Critical Point of Contact



BECKMAN COULTER
Biomek i-Series
• NON-STERILE
© 2016 Beckman Coulter, Inc.
Brea, CA 92821 U.S.A.

30 µl Pipette Tips
Tip Capacity: 30 µL
384 Tips/Tip Box
REF B85739

DNASE RNASE FREE
Certified
B85744-AA



ACCELERATING
answers



Table of Contents

Biomek Tips and Labware.....	3	Biomek Span-8 Fixed Probes	11
Types of Pipette Tips.....	4	Biomek Micro- and Deep-Well Plates	12
Placing Your Order.....	5	Biomek Modular Reservoir System.....	13
Size Information and Placing Your Order.....	6	Biomek Labware and Accessories.....	14
Biomek i-Series Non-Conductive Pipette Tips.....	7	Sustainability.....	16
Biomek i-Series Conductive Pipette Tips.....	10		

Biomek Tips and Labware

Excellence for the critical point of contact.

From disposable pipette tips to microplates and everything in between, we provide a full suite of accurate, precise performance you can depend on.

Biomek disposable tips from Beckman Coulter are the only tips designed, validated and approved for Biomek Liquid Handlers. Manufactured under stringent manufacturing processes, we guarantee proper fit and performance of Biomek tips on your Biomek liquid handler. Tip definitions, pipetting techniques, templates and color-coded tip racks are integrated into Biomek software to facilitate method development and make setup, method writing, and operation even more efficient.

As a crucial component to the most advanced liquid handling systems, Biomek tips from Beckman Coulter ensure high performance and quality control for the most exacting applications such as genetic analysis, protein research, cellular analysis and drug discovery.

Biomek tips are manufactured in a controlled manufacturing environment to minimize facility bio-burden. Our pipette tips, racks, and rack covers are molded out of premium-grade, virgin polypropylene to provide chemical resistance during sample preparation and to ensure only the highest quality product. Statistical process control (SPC) product sampling is conducted and samples are tested for biological parameters.

Biomek pipette tips from Beckman Coulter are certified to be*:

1

RNase- and DNase-free

Ensures high quality nucleic acid purification and reliable nucleic acid testing results.

2

DNA-free (human & mouse)/PCR inhibition

Addresses endotoxin-sensitive applications, free of contaminants that result in erroneous results and interfere with the PCR.

3

Pyrogen/endotoxin-free

Addresses endotoxin-sensitive applications, such as cell transformations and preparative protocols that use extracted biological material for in vivo experimentation.

4

Trace metal-free

Testing of the base, polypropylene tip resin reduces interferences due to chelation and denaturation.

*Based on sample results below detectable limits

Storage Conditions



It is recommended that Biomek Tips be stored in a clean and dry laboratory environment at room temperature. Exposure to high temperatures, radiation or vaporous chemicals/solvents may result in the deterioration of the design of the tip, and therefore its performance.

Pre-Sterile Info

STERILE

EO

STERILE

R

For applications requiring sterility, Beckman Coulter offers sterilized product, which is controlled under validated ethylene oxide processes, for those applications requiring sterile liquid handling. Products designated as “pre-sterile” are sterilized in accordance with ANSI/AAMI/ISO 11135 or 11137 guidelines, as appropriate. The sterilization processes certify a sterility assurance level (SAL) of 10⁻⁶. Autoclaving of Biomek Tips is not recommended, nor supported, due to the adverse effects it can have on the form and performance of the pipette tip.

Note: The use of the terms “pre-sterile” versus “sterile,” and “barrier” versus “filter,” respectively, are equivalent relative to product functionality and performance.

Types of Pipette Tips

Though small-in-size, its impact and importance in life-saving research is undeniably huge. Without pipette tips, tip-based liquid handling simply could not happen.

Pipette tips come in different types including non-sterile, pre-sterilized and filtered tips.



Non-Sterile

The most commonly used pipette tip. Non-sterile tips are often used in laboratory applications where sterility is not important to the experiment or test being performed.



Pre-sterilized/Sterile

Pre-sterilized tips are designed to prevent contamination and are free from living organisms. Our pre-sterilized tips are sterilized through an ethylene oxide process.



Filtered/Barrier

Filter tips are used to prevent potential contamination or damage of the pipetting system by aerosols. They are often used for pipetting nucleic acids or volatile liquids.



Multichannel

Multichannel tips are used on 96 - or 384 - channel pipetting heads in a variety of high-throughput applications which require multiple wells to be filled simultaneously.



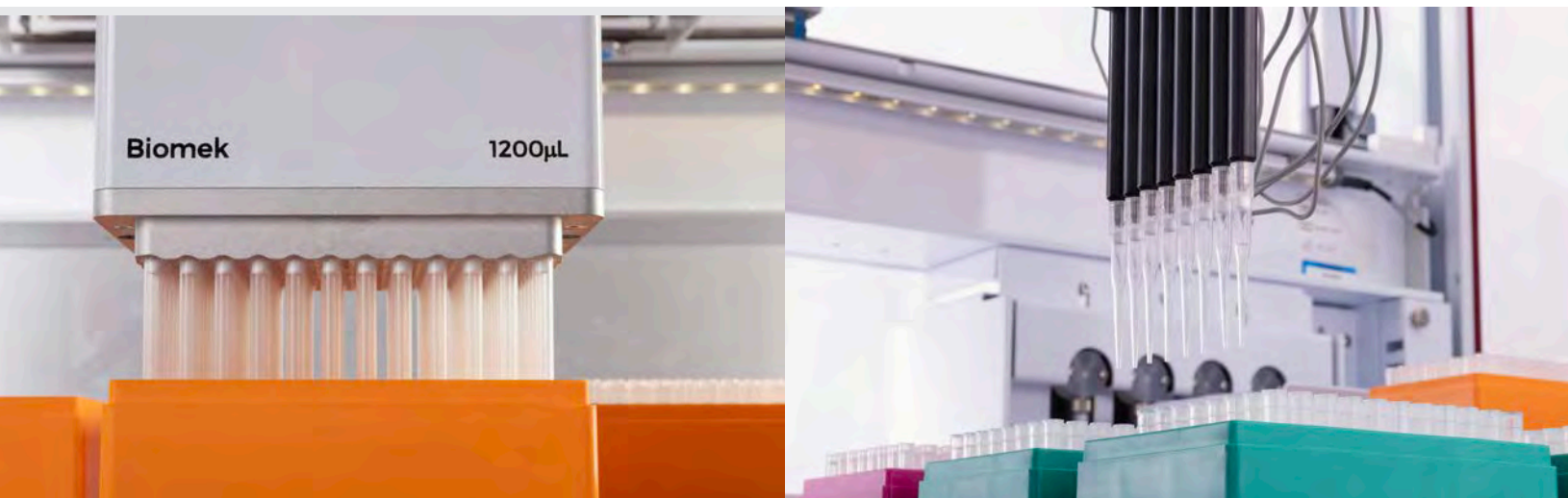
Span-8

Tips used on a Span-8 pipettor and which enable processes such as hit-picking, sample normalization and reformatting to and from a variety of labware types.



Fixed Probes

Fixed probes are used on the Span-8 pipettor and are great for piercing blood tubes or transfers through other tube septa and are used for pipetting chemicals that are incompatible with plastics. They are capable of pipetting larger volumes and are available in stainless steel or teflon-coated.



Types of Pipette Tips

Non-Conductive vs Conductive

Biomek tips are made from 100% virgin polypropylene in two materials.



Non-Conductive Tip

Clear/non-conductive. Made from 100% clear virgin polypropylene. These tips do not contain carbon.



Conductive Tip

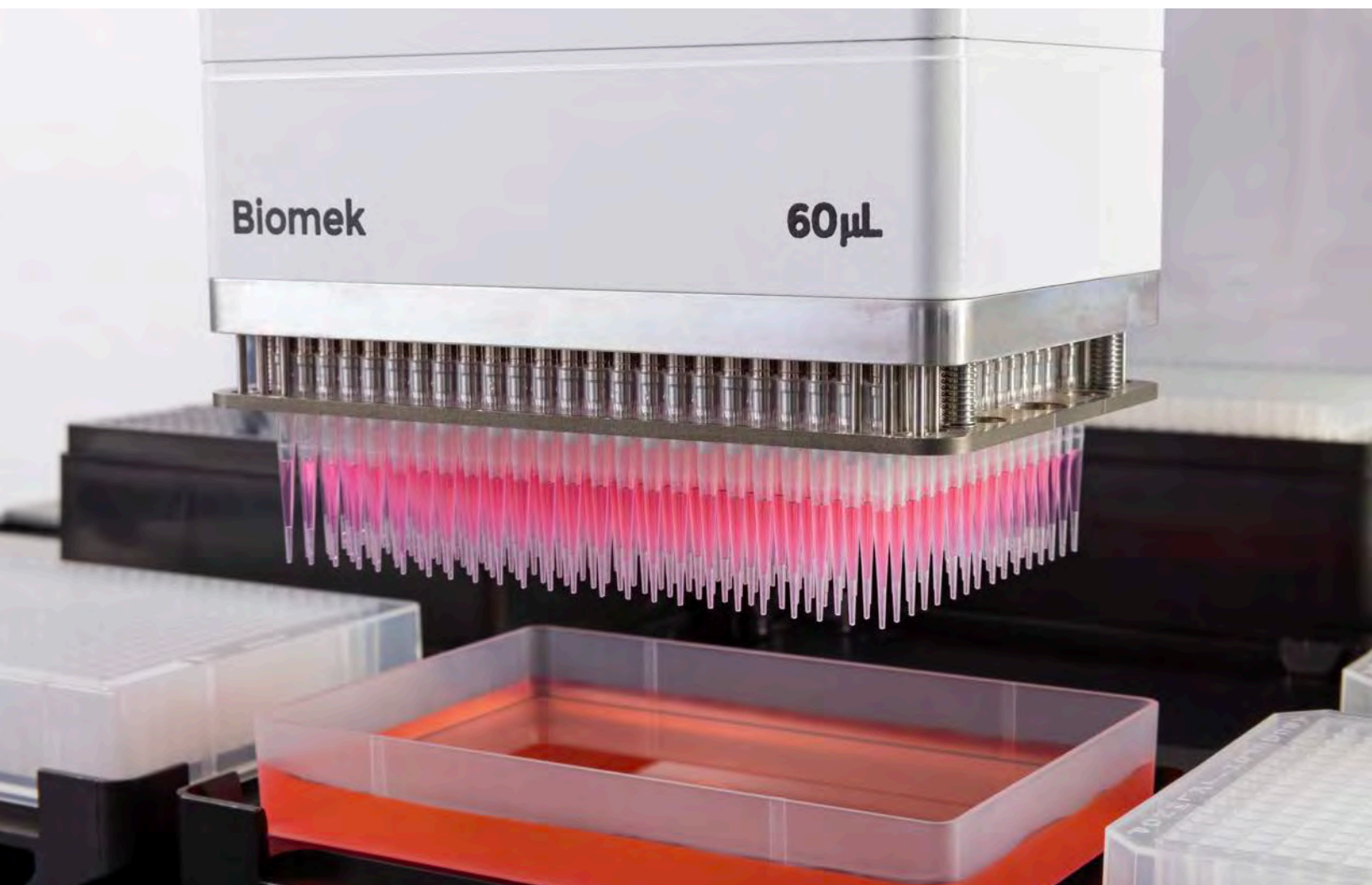
Black/conductive. Contain carbon that makes them electrically conductive. These tips can detect the liquid level when pipetting conducting liquids. They are available with filters and are used on the Biomek Span-8 Pipettor.

Wide Bore Tips



Wide Bore Tip

Biomek wide bore tips have a wider orifice that reduces resistance and shearing forces. These tips can be valuable when pipetting viscous liquids or delicate cell types or high molecular weight DNA when gentle pipetting is critical.



Size Information

The size of the tip determines the maximum volume of liquid that the tip can transfer. Here are a few examples:

Volume/Capacity Information

P30s		P30XL		P20		P50		P250		P1000	
Part No.	Capacity	Part No.	Capacity	Part No.	Capacity	Part No.	Capacity	Part No.	Capacity	Part No.	Capacity
B85739	30µL	B85753	50µL	B85764	80µL	B85881	90µL	B85903	230µL	B85940	1070µL
B85745	30µL	B85756	50µL	B85767	80µL	B85884	90µL	B85906	230µL	B85945	1070µL
B85749	25µL	B85760	40µL	B85771	40µL	B85888	50µL	B85911	190µL	B85955	1025µL

Placing Your Order

When you place a tip order, your order will be in case count.



P1000 Tips

1 tip rack (box) = 96 Tips
5 tip racks (box) = 1 Case

96 Tips x 5 Racks = 480 Tips
72 Cases = 1 Skid/Pallet



P20, P50, and P250 Tips

1 tip rack (box) = 96 Tips
10 tip racks (box) = 1 Case

96 Tips x 10 Racks = 960 Tips
90 Cases = 1 Skid/Pallet



P30 and P30XL Tips

1 tip rack (box) = 384 Tips
10 tip racks (box) = 1 Case

384 Tips x 10 Racks = 3840 Tips
90 Cases = 1 Skid/Pallet

Biomek i-Series Pipette Tips

Non-conductive tips for use with 60 µL Multichannel-384 Head.



30 µL Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85739	Biomek P30 Tips Non-Sterile			30 µL	384 tips/tip box, 10 tip boxes/case
B85745	Biomek P30 Tips Pre-Sterile	•		30 µL	384 tips/tip box, 10 tip boxes/case
B85749	Biomek P30 Tips Pre-Sterile with Barrier	•	•	25 µL	384 tips/tip box, 10 tip boxes/case



50 µL Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85753	Biomek P30XL Tips Non-Sterile			50 µL	384 tips/tip box, 10 tip boxes/case
B85756	Biomek P30XL Tips Pre-Sterile	•		50 µL	384 tips/tip box, 10 tip boxes/case
B85760	Biomek P30XL Tips Pre-Sterile with Barrier	•	•	40 µL	384 tips/tip box, 10 tip boxes/case



Biomek i-Series Pipette Tips

Non-conductive tips for use with 300 & 1200 μL Multichannel-96 Heads & Span-8 Pod.



80 μL Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85764	Biomek P20 Tips Non-Sterile			80 μL	96 tips/tip box, 10 tip boxes/case
B85767	Biomek P20 Tips Pre-Sterile	●		80 μL	96 tips/tip box, 10 tip boxes/case
B85771	Biomek P20 Tips Pre-Sterile with Barrier	●	●	40 μL	96 tips/tip box, 10 tip boxes/case



90 μL Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85881	Biomek P50 Tips Non-Sterile			90 μL	96 tips/tip box, 10 tip boxes/case
B85884	Biomek P50 Tips Pre-Sterile	●		90 μL	96 tips/tip box, 10 tip boxes/case
B85888	Biomek P50 Tips Pre-Sterile with Barrier	●	●	50 μL	96 tips/tip box, 10 tip boxes/case



230 μL Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85903	Biomek P250 Tips Non-Sterile			230 μL	96 tips/tip box, 10 tip boxes/case
B85906	Biomek P250 Tips Pre-Sterile	●		230 μL	96 tips/tip box, 10 tip boxes/case
B85911	Biomek P250 Tips Pre-Sterile with Barrier	●	●	190 μL	96 tips/tip box, 10 tip boxes/case



230 μL Wide Bore Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85926	Biomek P250 Wide Bore Tips Non-Sterile			230 μL	96 tips/tip box, 10 tip boxes/case
B85929	Biomek P250 Wide Bore Tips Pre-Sterile	●		230 μL	96 tips/tip box, 10 tip boxes/case
B85936	Biomek P250 Wide Bore Tips Pre-Sterile with Barrier	●	●	190 μL	96 tips/tip box, 10 tip boxes/case



Biomek i-Series Pipette Tips

Non-conductive tips for use with 300 & 1200 μL Multichannel-96 Heads & Span-8 Pod.



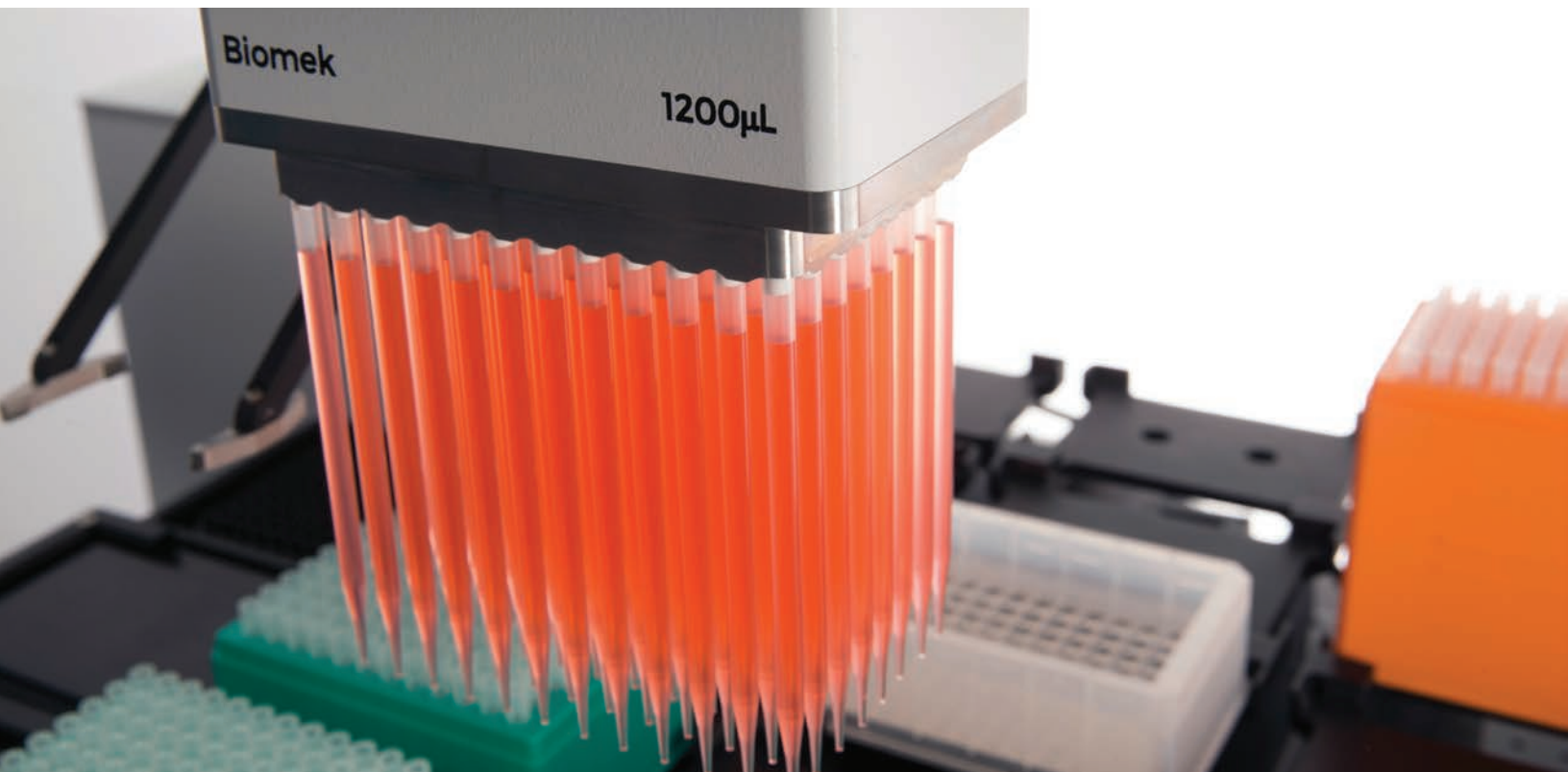
1070 μL Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85940	Biomek P1000 Tips Non-Sterile			1070 μL	96 tips/tip box, 5 tip boxes/case
B85945	Biomek P1000 Tips Pre-Sterile	•		1070 μL	96 tips/tip box, 5 tip boxes/case
B85955	Biomek P1000 Tips Pre-Sterile with Barrier	•	•	1025 μL	96 tips/tip box, 5 tip boxes/case



1070 μL Wide Bore Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85971	Biomek P1000 Wide Bore Tips Non-Sterile			1070 μL	96 tips/tip box, 5 tip boxes/case
B85975	Biomek P1000 Wide Bore Tips Pre-Sterile	•		1070 μL	96 tips/tip box, 5 tip boxes/case
B85981	Biomek P1000 Wide Bore Tips Pre-Sterile with Barrier	•	•	1025 μL	96 tips/tip box, 5 tip boxes/case



Biomek i-Series Pipette Tips

Conductive tips for use with Span-8 Pod to enable Liquid Level Sensing.



80µL Conductive Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85775	Biomek P20 Conductive Tips Non-Sterile			80 µL	96 tips/tip box, 10 tip boxes/case
B85872	Biomek P20 Conductive Tips Pre-Sterile	●		80 µL	96 tips/tip box, 10 tip boxes/case
B85877	Biomek P20 Conductive Tips Pre-Sterile with Barrier	●	●	40 µL	96 tips/tip box, 10 tip boxes/case



90µL Conductive Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85892	Biomek P50 Conductive Tips Non-Sterile			90 µL	96 tips/tip box, 10 tip boxes/case
B85894	Biomek P50 Conductive Tips Pre-Sterile	●		90 µL	96 tips/tip box, 10 tip boxes/case
B85899	Biomek P50 Conductive Tips Pre-Sterile with Barrier	●	●	50 µL	96 tips/tip box, 10 tip boxes/case



230µL Conductive Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85915	Biomek P250 Conductive Tips Non-Sterile			230 µL	96 tips/tip box, 10 tip boxes/case
B85917	Biomek P250 Conductive Tips Pre-Sterile	●		230 µL	96 tips/tip box, 10 tip boxes/case
B85922	Biomek P250 Conductive Tips Pre-Sterile with Barrier	●	●	190 µL	96 tips/tip box, 10 tip boxes/case



1070µL Conductive Pipette Tips

Part No.	Description	Sterile	Filter	Tip Volume	Packaging
B85959	Biomek P1000 Conductive Tips Non-Sterile			1070 µL	96 tips/tip box, 5 tip boxes/case
B85961	Biomek P1000 Conductive Tips Pre-Sterile	●		1070 µL	96 tips/tip box, 5 tip boxes/case
B85965	Biomek P1000 Conductive Tips Pre-Sterile with Barrier	●	●	1025 µL	96 tips/tip box, 5 tip boxes/case



Biomek Span-8 Fixed Probes

For use with the Biomek Span-8 Pod.

Fixed Probes

Part No.	Description	Probe Length	Packaging
719543	Stainless Steel Fixed Probe*	60 mm	1/case
719808	Stainless Steel Fixed Probe*	60 mm	8/case
719544	Stainless Steel Fixed Probe*	100 mm	1/case
719810	Stainless Steel Fixed Probe*	100 mm	8/case
A34967	Teflon-Coated Fixed Probe	60 mm	1/case
A39379	Teflon-Coated Fixed Probe	60 mm	8/case
A34966	Teflon-Coated Fixed Probe	100 mm	1/case
A39377	Teflon-Coated Fixed Probe	100 mm	8/case
717931	Teflon-Coated Fixed Probe*	60 mm	1/case
719807	Teflon-Coated Fixed Probe*	60 mm	8/case
717933	Teflon-Coated Fixed Probe*	100 mm	1/case
719809	Teflon-Coated Fixed Probe*	100 mm	8/case
987873	Stainless Steel Piercing Probe*	120 mm	1/pack
987872	Stainless Steel Piercing Probe*	120 mm	4/pack
987870	Stainless Steel Piercing Probe*	120 mm	8/pack
A49675	Large Volume Tubing Kit†	—	—



* Requires Biomek Large Volume Tubing Kit (Part No. A49675).

† The Biomek Large Volume Tubing Kit includes tubing and other parts necessary to convert a system with low-volume tubing to large-volume tubing for use with large-volume syringes or stainless steel piercing probes.

100 mm and 120 mm probes are compatible with Biomek NX[®]/FX[®] and i-Series Span-8 Pods.

Teflon is a registered trademark of E.I. du Pont de Nemours and Company.



Biomek Micro- and Deep-Well Plates

For use with Biomek Automated Workstations.

Deep-well plates and microplates are used for sample preparation, sample storage, and fraction collection.

Biomek 96 Round Deep-Well Plates

Part No.	Description	Color	Well Bottom	Well Volume	Packaging
267001	Non-sterile, Polystyrene	Clear	Round	1 mL	24/case
267004	Sterile, Polystyrene	Clear	Round	1 mL	24/case
267006	Non-sterile, Polypropylene	Natural	Round	1 mL	24/case
267007	Sterile, Polypropylene	Natural	Round	1 mL	24/case



Biomek 96 Square Deep-Well Plates

Part No.	Description	Color	Well Bottom	Well Volume	Packaging
140504	Non-sterile, Polypropylene	Natural	Round	2 mL	24/case
609681	Sterile, Polypropylene	Natural	Round	2 mL	24/case



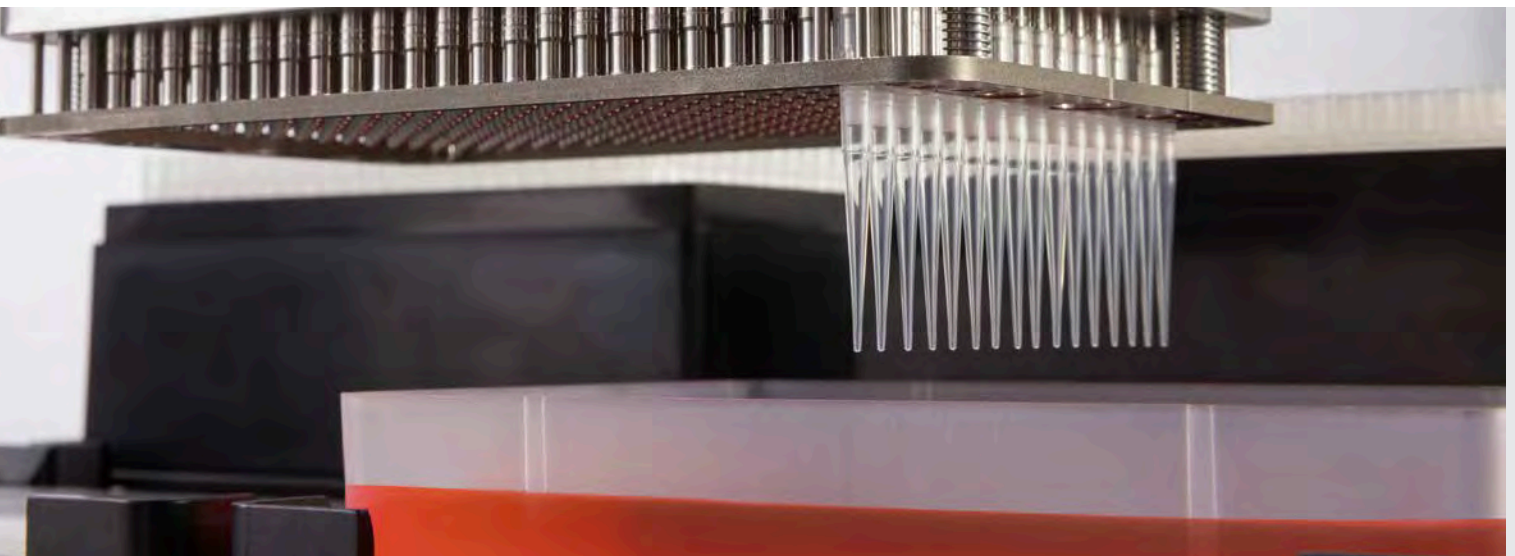
Biomek 96-Well Microplate

Part No.	Description	Color	Well Bottom	Well Volume	Packaging
609844	Non-sterile, Polystyrene	Clear	Flat	300 μ L	100/case
609801	Non-sterile, Polypropylene	Natural	Conical	200 μ L	25/case



For optimum sealing, use Biomek Aluminum Foil Seals (Part No. 538619).

For Biomek 96 Round Deep-Well Plates, you can also use Biomek Cap Mats (Part No. 267002 or 267005).



Biomek Modular Reservoir System

For use with Biomek Automated Workstations.

Reservoirs are essential accessories used in laboratories for temporarily storing reagents or samples during pipetting applications.

Full Reservoirs

(Span-8 Pod Use Only)

Part No.	Description	Capacity	Packaging
372784	Sterile, Polypropylene	150 mL	12/case
534678	Sterile, Polypropylene, Nonpyrogenic	150 mL	12/case



Half Reservoirs

Part No.	Description	Capacity	Packaging
372786	Sterile, Polypropylene	75 mL	24/case
534681	Sterile, Polypropylene, Nonpyrogenic	75 mL	24/case
372794	Sterile, Polypropylene, Divided by Width	37 mL/section	24/case



Quarter Reservoirs

Part No.	Description	Capacity	Packaging
372790	Sterile, Polypropylene	40 mL	48/case
534680	Sterile, Polypropylene, Nonpyrogenic	40 mL	48/case
372792	Sterile, Polypropylene, Divided by Width	38 mL (19 mL/section)	48/case
372788	Sterile, Polypropylene, Divided by Length	38 mL (19 mL/section)	48/case

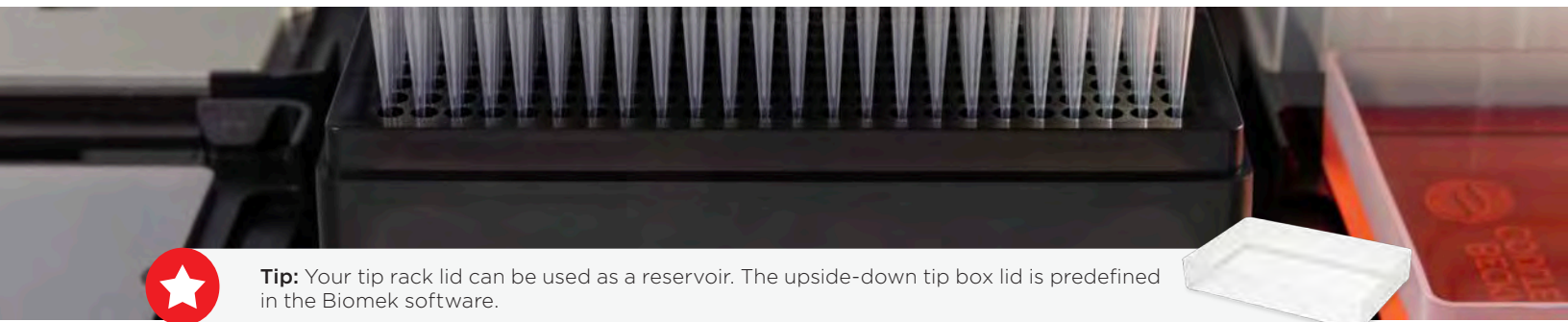


Frame For Reservoirs

Part No.	Description	Capacity	Packaging
372795	—	—	1 each



Each frame can accommodate up to two (2) half reservoirs or four (4) quarter reservoirs. Full reservoirs do not require a reservoir frame.



Tip: Your tip rack lid can be used as a reservoir. The upside-down tip box lid is predefined in the Biomek software.

Biomek Labware and Accessories

For use with Biomek Automated Workstations.

Tube racks provide organized storage to various sizes of glass, polystyrene and polypropylene tubes. Tube rack inserts ensure a proper fit of the tubes.

Biomek Tube Rack

Part No.	Description	Color	Packaging
373661	24-Position	Black	1 each



This 24-Position Tube Rack holds various sizes of glass, polystyrene, and polypropylene tubes, Microfuge tubes, and cryovials. Use inserts to ensure proper tube fit.

Biomek Tube Rack Inserts

Part No.	Description	Color	Packaging
373699	10 mm Diameter	Black	25/case
373696	11 mm Diameter (Securely holds 1.5 mL Microfuge tube flip cap in place)	White	25/case
373698	12 mm Diameter	Red	25/case
373697	13 mm Diameter	Blue	25/case



1.5 mL Microcentrifuge Tubes

Part No.	Description	Color	Packaging
356090	Polypropylene	Clear	500/case
356091	Polypropylene	Blue	500/case
356092	Polypropylene	Green	500/case
356093	Polypropylene	Yellow	500/case
356094	Polypropylene	Orange	500/case
357448	Polypropylene	Natural	500/case



Microcentrifuge Tubes, with 11 mm inserts (Part No. 373696), fit into the Biomek 24-Position Tube Rack (Part No. 373661).

Biomek Labware and Accessories Continued

For use with Biomek Automated Workstations.

Biomek Cap Mats

Cap mats are made from research-grade silicone to tightly seal 96-round deep-well plates.

Part No.	Description	Color	Packaging
267002	Non-sterile	Natural	12/case
267005	Sterile	Natural	12/case



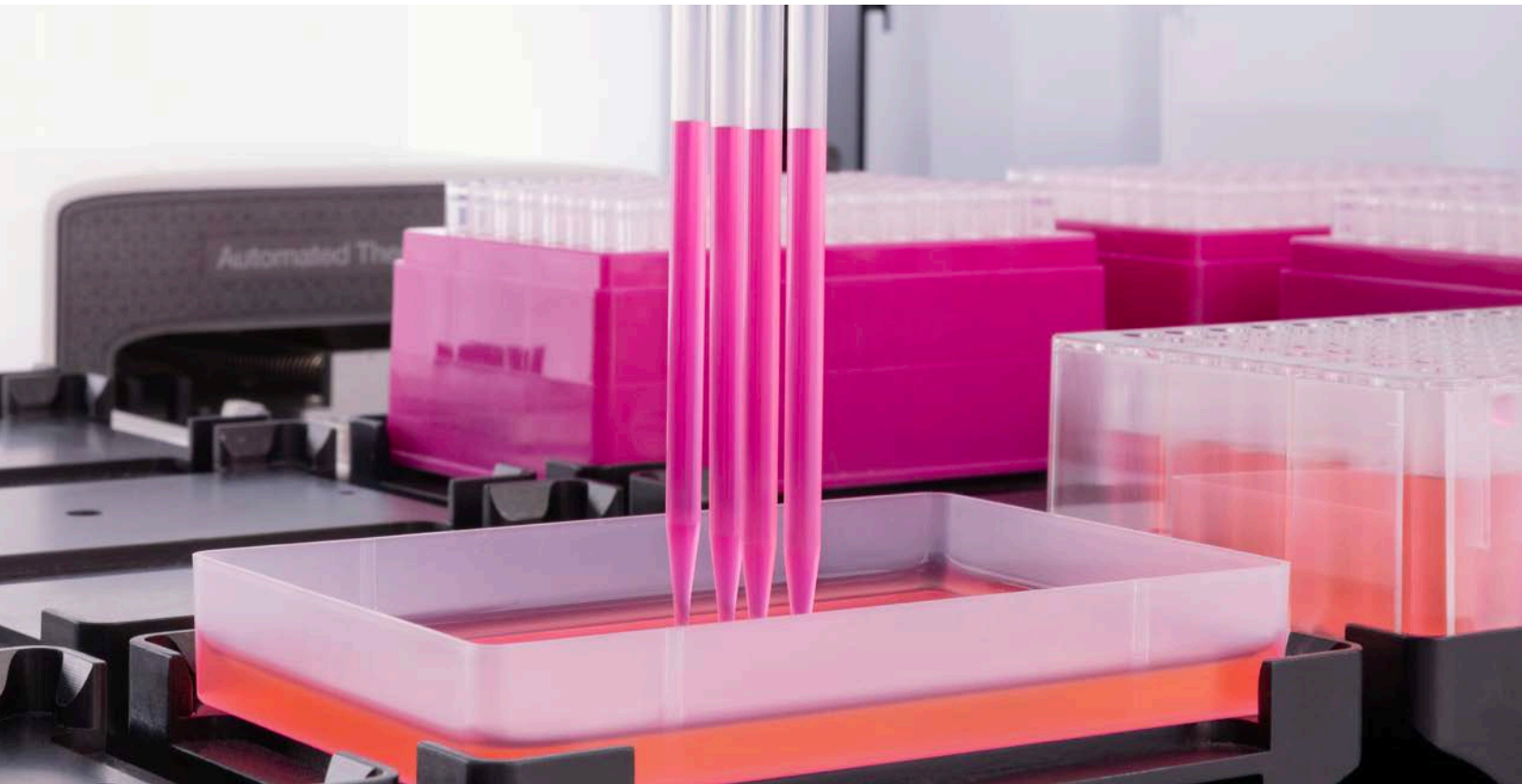
Biomek Aluminum Foil Seals

Aluminum foil seals are used to tightly seal 96- and 384-format plates.

Part No.	Description	Packaging
538619	Aluminum Foil Seals	100/case
538618	Roller for Aluminum Foil Seals	1 each



Use roller (Part No. 538618) to ensure complete plate sealing.



Beckman Coulter Life Sciences & Sustainable Innovation

Plastic circularity.

Beckman Coulter Life Sciences in partnership with Polycarbin, are working to divert lab plastics away from landfills and incinerators, keeping materials within the supply chain for recycling and reuse. Please contact your Beckman Coulter Life Sciences sales representative for more details about our recycling solution.

Environmental Awareness & Sustainability



Biomek tips, racks, and rack covers are molded from polypropylene resin to provide chemical resistance during sample preparation. The Society of the Plastics Industry (SPI) identifies polypropylene with an internationally recognized resin identification code number of 5 for recycling purposes.



Biomek tip packaging is comprised of 100% recyclable corrugated fiberboard.

POLYCARBIN



Accelerating Answers to Your Important Questions

With a relentless mission to empower those seeking answers to life's important scientific questions, Beckman Coulter Life Sciences is a preferred laboratory partner providing automation and innovation solutions for centrifugation, flow cytometry, genomics, particle analysis and liquid handling workflows.

With a legacy dating back to 1935, our technologies reduce manual laboratory processes and can provide greater speed, accuracy, and advanced analytics to help you accelerate answers in your lab.

To order, visit beckman.com or contact your local representative.



Biomek Automated Workstations are not intended or validated for use in the diagnosis of disease or other conditions.

© 2024 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
2024-GBL-EN-106312-v1

