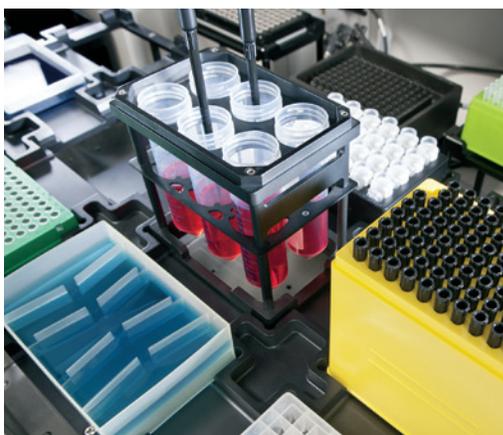


Optimize Large Volume Pipetting Transfers

Biomek Span-8 P1000 Pipette Tips

Biomek Span-8 P1000 Pipette Tips are now available in three designs: 1) non-conductive, 2) wide bore, and 3) conductive. The capability to transfer a full 1 mL of samples or reagents in a single pipetting step provides greater efficiency and timesavings during sample preparation. The P1000 tips are commonly used for large volume transfers applied to nucleic acid extraction and purification protocols, "seed and feed" cellular applications, and generic reagent pipetting. The overall length of the tips allows access to the full contents of long, narrow tubes, such as 12 x 75 mm and 16 x 100 mm test tubes, 15 and 50 mL centrifugation tubes, and 1 and 2 mL deep-well plates.



For the best in system performance, use only Beckman Coulter validated tips on Biomek workstations.

1 Faster sample preparation

Increase sample preparation times by transferring a full 1 mL of samples or reagents in a single pipetting action without having to split transfer volumes.

2 Observe pipetting transfers

Instill confidence during method development by visualizing pipetted volumes with non-conductive (clear) tips.

3 Ensure biological integrity of samples

Reduce flow-induced shear strain on native biological sample matrices with wide bore tips.

4 Process unknown sample volumes

Enable liquid level sensing with conductive tips for processing unknown sample volumes.

5 Optimize pipetting performance

Improve precision in critical dilution steps with single step transfers.

6 Minimize contamination risks

Eliminate the risk of contamination of samples, assay plates, and reagent reservoirs through the use of disposable pipette tips and by reducing the number of "fly-over" movements.

7 Free up deck space

Conserve valuable Biomek deck space due to less tip rack consumption.

Discovery
in motion.



Life Sciences

Non-conductive vs. Conductive Tips

For added confidence in Biomek pipetting performance during method development, non-conductive tips offer the advantage of being clear (versus black) to allow for visual inspection of aspirated volumes. With the clear tips, researchers are able to see how much volume they have aspirated into the tips, and the uniformity of aspirated volume across tips. Black, conductive P1000 Span-8 tips do not provide visual inspection, but do offer the advantage of enabling liquid level sensing of pipetted matrices on the Biomek NX^P and FX^P Span-8 instruments. The Biomek Span-8 pod incorporates this functionality to sense the level of liquid in a piece of labware and to track the volume of liquid as it is aspirated from, or dispensed to, the labware. This feature is invaluable when pipetting from labware with different liquid volumes and heights, and minimizes the amount of pipetted sample from wetting the outside of the tip and causing cross-contamination.

Ordering Information

P1000 Tips

Part No.	Description	Max Volume	Packaging
B01122	Biomek Span-8 P1000 Tips Non-sterile	1050 µL	96/rack, 5 racks/case
B01123	Biomek Span-8 P1000 Tips Pre-sterile	1050 µL	96/rack, 5 racks/case
B01124	Biomek Span-8 P1000 Tips Pre-sterile with Barrier	1025 µL	96/rack, 5 racks/case



P1000 Wide Bore Tips

Part No.	Description	Max Volume	Packaging
B01112	Biomek Span-8 P1000 Wide Bore Tips Non-sterile	1050 µL	96/rack, 5 racks/case
B01113	Biomek Span-8 P1000 Wide Bore Tips Pre-sterile	1050 µL	96/rack, 5 racks/case
B01114	Biomek Span-8 P1000 Wide Bore Tips Pre-sterile with Barrier	1025 µL	96/rack, 5 racks/case



P1000 Conductive Tips

Part No.	Description	Max Volume	Packaging
987935	Biomek Span-8 P1000 Conductive Tips Non-sterile	1050 µL	96/rack, 5 racks/case
987936	Biomek Span-8 P1000 Conductive Tips Pre-sterile	1050 µL	96/rack, 5 racks/case
987925	Biomek Span-8 P1000 Conductive Tips Pre-sterile with Barrier	1025 µL	96/rack, 5 racks/case



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.

Wide Bore (Clear) Tips

P1000 Span-8 Wide Bore tips are non-conductive (clear) tips, which are designed with a pipetting orifice twice the size, e.g. 1.4 mm, of standard P1000 tips. Biomek Wide Bore Tips minimize the inadvertent harm, which can be applied to native biological sample matrices during sample preparation. Loss of biological sample integrity can occur when pipetting fragile cell suspensions, hybridomas, embryoid bodies, algal coagulates, protein aggregates, and genomic materials due to the pipette tips' small internal orifice diameter and the speed of aspiration and dispense during pipetting.⁽¹⁾

In addition, Biomek Wide Bore Tips reduce sample preparation times by allowing faster pipetting speeds of viscous reagent solutions, such as glycerol stocks, or bead mixtures. The larger diameter bore also minimizes tip clogging when pipetting particulate-laden or fibrous sample matrices as found in biomass studies and agricultural research.

Bio-certifications

Biomek tips are certified to be free of RNase, DNase, DNA (human and mouse) / PCR inhibition, pyrogen / endotoxins, and trace metals.

⁽¹⁾ For additional information, request: Technical Information Bulletin, IB-17295A, *Analyzing the Effect of Pipette Tip Geometries on Fluid Velocity and Shear Strain Rate; Biomek Wide Bore vs. Standard Pipette Tips*, Robert D. Lund, Beckman Coulter Life Sciences, Indianapolis, IN.