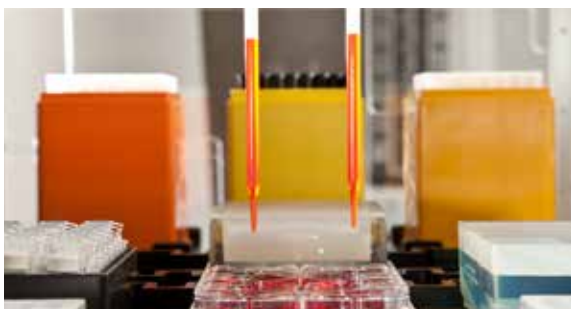




Ensure Biological Integrity of Samples

BIOMEK WIDE BORE PIPETTE TIPS

Biomek Wide Bore Pipette 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. Biomek Wide Bore Tips have been designed to address such sample preparation challenges with a pipetting orifice twice the size of standard pipette tips.¹



Comparison of Standard vs. Wide Bore Orifice Pipette Tip Diameters



* Enlarged for illustrative purposes only.

1 Ensure biological integrity of samples

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

2 Minimize DNA shearing

Avoid the risk of inadvertent DNA shearing, while preparing samples during nucleic acid extraction and purification protocols.

3 Improve sample throughput

Increase throughput when pipetting viscous reagent solutions, such as glycerol stocks, or bead mixtures.

4 Minimize cell lysis

Prevent inadvertent cell lysis when pipetting fragile cell suspensions, hybridomas, or embryoid bodies.

5 Reduce aggregate disruptions

Maintain the natural state of protein aggregates in therapeutic formulation studies and algal coagulated in biomass studies.

6 Prevent tip clogging

Avoid obstructions due to heavy particulate or fibrous sample matrices as encountered in biomass studies and agricultural research.

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

Improved Pipetting Efficiency

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 1 mL tips are commonly used for large volume transfers applied to nucleic acid extraction and purification protocols, “seed and feed” cellular applications, and general 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 added confidence in Biomek pipetting performance during method development, the non-conductive, Wide

Bore Tips offer the advantage of being clear (versus black, conductive) 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.

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 those found in biomass studies and agricultural research.

Ordering Information

Biomek Wide Bore Pipette Tips

	NON-STERILE	STERILE (PRE-STERILE)	STERILE WITH FILTER (BARRIER)	PACKAGING
For use with Biomek 4000, and Biomek NX [®] and FX [®] Span-8 Automated Workstations.*				
● P250	B01101	B01102	B01103	96 tips/tip box, 10 tip boxes/case
● P1000	B01112	B01113	B01114	96 tips/tip box, 5 tip boxes/case
For use with Biomek i-Series Automated Workstations.				
● 230/190 µL	B85926	B85929	B85936	96 tips/tip box, 10 tip boxes/case
● 1070/1025 µL	B85971	B85975	B85981	96 tips/tip box, 5 tip boxes/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.

Bio-certifications

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