

# Increase productivity and accuracy in your cell deposition.

MoFlo XDP CyClone

Blood Banking  
Capillary Electrophoresis  
Centrifugation  
**Flow Cytometry**  
Genomics  
Lab Automation  
Lab Tools  
Particle Characterization



Increase productivity with high-speed, high-precision cell deposition into any multi-well plate or slide format you choose.

The CyClone enables accurate cell deposition with the MoFlo XDP high-performance cell sorter.

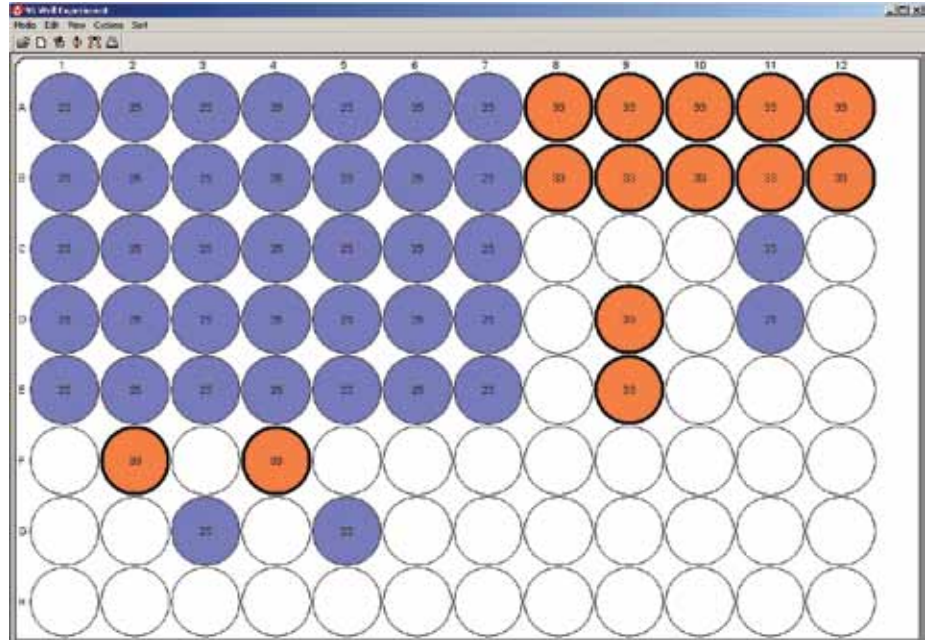
CyClone's precise positioning allows single cells to be sorted into standard or specialized plates and onto slides for your customized sort profiles.

Stable fluidics combined with the CyClone performance specifications; enable sorting into 1536-well plates on the MoFlo XDP.

# MoFlo XDP CyClone

CyClone sets the bar for performance, versatility and speed. From a one percent starting target population, CyClone can execute single cell deposition into a 96-well plate in less than 50 seconds with 99.9 percent purity. The CyClone arm enables single cell sorting into 1536-well plates and onto the smaller surface area of a slide. This ensures the more common 96-well and 384-well plate sorting to be more precise.

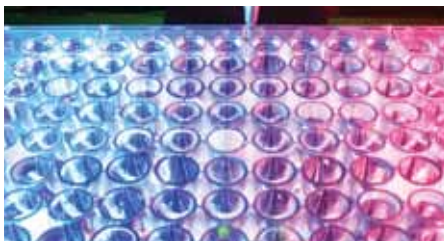
- Purity greater than 99.9 percent.
- Single cell accuracy is assured with a proprietary algorithm for “one-and-one only” cell deposition.
- Deposition positioning accuracy of  $\pm 43.541 \mu\text{m}$  on the y-axis and  $\pm 126.27 \mu\text{m}$  the x-axis.
- Well-to-well position speed of 10 cm per second.
- The retractable arm makes the switch between cloning and traditional tube sorting fast and easy.



Summit Software user interface for the CyClone has the capability to customize sort setups for multi-well plate or slide deposition.

## CyClone Specifications

<b>Supported Receptacle</b>	Standard multi-well plate formats including 6, 12, 24, 48, 96, 384, 1536; formats on standard custom microscope slides
<b>Maximum Density of Discrete Droplet Deposition</b>	1818 rows by 952 columns for a total of 1.73 million objects on an Omni tray (80 mm x 120 mm)
<b>Cell Deposition Accuracy</b>	Standard deviation of $\pm 126.27 \mu\text{m}$ on the x-axis; $\pm 43.541 \mu\text{m}$ on the y-axis
<b>Maximum Well-to-Well Speed</b>	10 cm per second



Standard 96-well plate



96-well deep plate

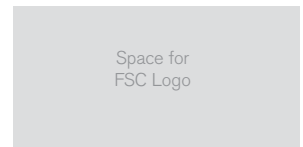


Standard 1536-well plate

[www.coulterflow.com](http://www.coulterflow.com)



For Research Use Only. Not for use in diagnostic procedures. Optional parameters available outside U.S.  
 MoFlo XDP is a class 1 laser.  
 Beckman Coulter, the stylized logo and MoFlo are registered trademarks of Beckman Coulter, Inc.  
 CyClone is a trademark of Beckman Coulter, Inc.



Printed with soy ink

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at [www.beckmancoulter.com](http://www.beckmancoulter.com)

B2010-11847-DG-2.5

© 2010 Beckman Coulter, Inc.

PRINTED IN U.S.A.