

Flexible, automated
sample handling.

MoFlo™ XDP SmartSampler

Chemistry
Lab Automation
Information Systems
Molecular Diagnostics
Immunodiagnosics
Centrifugation
Disease Management
Hematology
Hemostasis
Flow Cytometry
Primary Care



Designed around the way you work, with an open platform and automated features to increase productivity.

SmartSampler is the latest product innovation for sample introduction on the MoFlo XDP high-performance cell sorter, streamlining sample handling operations normally performed manually by the instrument operator.

SmartSampler easily accommodates tube sizes from 0.5 to 50 mL for your custom formats. Other features include agitation and temperature control, as well as single push button operation for starting or stopping a run.



MoFlo XDP SmartSampler

From small volumes to large, your sample requirements change frequently. Now, use a sample input system that saves both time and money.

Flexible Sample Input

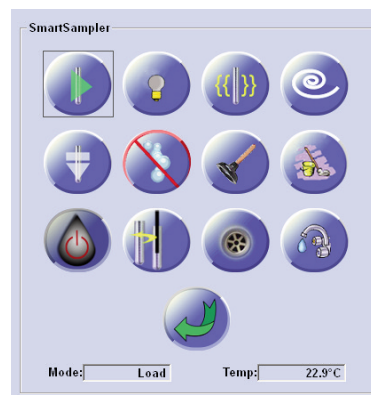
- Unique small-tube format reduces sample dilution requirements.
- Large-tube format for longer continuous sorts.
- Agitation keeps cells in homogeneous solution, critical for even distribution of rare events.
- Temperature control maintains viability for cell re-culture.
- End-of-sample (air) detection enables an entire sample run with minimal waste of source material and reagents.
- Automated backflush and cleaning for dense or sticky samples reduces clogs and facilitates high-speed sorting.
- Front panel button and LED displays promote one touch operation for high throughput and automated operation.
- The Lexan chamber seals the sample environment for superior control and operator safety.
- Filtration removes potential aerosols as small as 0.3 μm in exhausted air at 99.995 percent efficiency, protecting the operator.

aXcess Control Panel for SmartSampler Screen



SmartSampler menu button

SmartSampler abbreviated menu



SmartSampler full menu

SmartSampler Specifications

| | |
|---------------------|-----------------------------|
| Tube Sizes | 0.5 – 50 mL |
| Nozzle Sizes | 70 – 200 μm |
| Temperature Control | -4 – 40 °C |
| Sample Pressure | 5 – 85 psi |
| Dimensions | 18 cm W x 10 cm D x 44 cm H |

For Research Use Only. Not for use in diagnostic procedures.

Australia, Gladesville (61) 2 9844-6000 Canada, Mississauga (1) 905 819 1234 China, Beijing (86) 10 6515 6028
 Czech Republic, Prague (420) 272 01 73 32 Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707
 France, Villepinte (33) 1 49 90 90 00 Germany, Krefeld (49) 2151 33 35 Hong Kong (852) 2814 7431
 India, Mumbai (91) 22 3080 5101 Italy, Cassina de' Pecchi, Milan (39) 02 953921 Japan, Tokyo (81) 3 5530 8500 Korea, Seoul (82) 2 404 2146
 Latin America (1) (305) 380 4709 Mexico, Mexico City (001) 52 55 9183 2800 Netherlands, Mijdrecht (31) 297 230630 Puerto Rico (787) 747 3335
 Singapore (65) 6339 3633 South Africa/Sub-Saharan Africa, Johannesburg (27) 11 805 2014/5 Spain, Madrid (34) 91 3836080
 Sweden, Bromma (46) 8 564 85 900 Switzerland, Nyon (41) 0800 850 810 Taiwan, Taipei (886) 2 2378 3456 Turkey, Istanbul (90) 216 309 1900
 UK, High Wycombe (44) 01494 441181 USA, Fullerton, CA (1) 800 742 2345