

Two-way communication,
in one **powerful solution.**

Instrument Manager* Version 8.05 with Bidirectional LIS Connectivity for FC 500 Systems

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



Powerful features deliver the information you need:

- Test requests and patient demographics are automatically sent to the FC 500
- Worklists can be automatically created
- The workstation delivers robust graphics review capabilities with histograms that can be transmitted to the LIS
- Track samples throughout the testing process, including all bar-coded daughter tubes
- Use Beckman Coulter's FP 1000 cell preparation system or the TQ-Prep™ and PrepPlus™ 2 workstations for added simplicity and efficiency

Improved safety measures give you peace of mind:

- Manual data entry is eliminated, reducing the risk of error and improving patient safety
- The FC 500 CXP software automatically verifies sample position and test
- Results are password protected
- The software helps users meet patient safety goals from accrediting agencies such as CAP and JCAHO

Enhanced data management saves time and gives you more control:

- Automatically release test results that meet specific user-defined criteria
- Create conditional statements in order to flag patient results that are outside of user-defined ranges or to flag Quality Control violations
- Create specific decision rules for particular patients or doctors

Innovative design means software is easy to use:

- Installation is simple
- Report templates are customizable
- Minimal operator intervention is required
- Current Beckman Coulter customers with Instrument Manager can download the software for free

Imagine your flow cytometer automatically downloading test requests and patient demographics from the LIS. No need to enter the data manually. No need for any other instrument.

Now imagine worklists that are automatically created for you. Powerful data management capabilities that are just a click away. Sample identification that improves patient safety.

Beckman Coulter delivers this and more through the Instrument Manager (IM) v.8.05 Data Management Package.

When you add this capability to your FC 500, you establish a path for true bidirectional information flow between your instrument and the LIS.

* *Instrument Manager* is a registered trademark of Data Innovations, Inc.

Bidirectional LIS Connectivity for FC 500 Systems

Simplicity. Efficiency. Features that make your day easier.

Instrument Manager Version 8.05 Bidirectional Data Management for the FC 500 System. Welcome to a new level of flow cytometry testing.



Please call for ordering information.



Simplify • Automate • Innovate

Australia, Gladesville (61) 2 9844 6000 Canada, Mississauga (1) 905 819 1234 China, Beijing (86) 10 6515 6028
Czech Republic, Prague (420) 267 00 83 66 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 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 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, Brea, CA (1) 800 352 3433, (1) 714 993 5321

B2007-7778-PO-2K

www.beckmancoulter.com

© 2007 Beckman Coulter, Inc.

DIAG-PRINTED IN U.S.A.