



Meet your lysing needs with unmatched reproducibility

The ImmunoPrep™ Reagent System.

Regardless of which workstation best fits your needs, Beckman Coulter's ImmunoPrep reagent system ensures that your lab can go from whole blood to lysis in just seconds. Designed specifically for use on the Q-Prep, TQ-Prep and FP 1000 workstations, this system provides the fastest no-wash, flow cytometry-ready preparation, while maintaining cellular integrity and unmatched reproducibility between samples. Benefits also include:

- Gentle lysing action to reduce cell stress and minimize cell shrinkage, while maintaining cellular morphology
- Minimizing debris or platelet adhesion to leukocytes
- Reduction of non-specific binding and excess manipulation of cells
- Prevention of selective cell loss due to centrifugation

Semi-automated performance at an incredible value

The Q-Prep™ Workstation.

When the Q-Prep Workstation was first introduced, it paved the way for improved sample preparation – virtually revolutionizing the traditional manual methods. Today, the Q-Prep remains a cost-effective choice for many lower-volume laboratories and offers a variety of features, including:

- Standardized cell preparation with excellent reproducibility
- The fastest patient results with a 35-second cell lyse process
- A reduction in labor cost, compared to manual methods

Service And Support You Can Count On

As a world leader in innovative service and support, Beckman Coulter ensures that your flow cytometry systems perform at peak efficiency throughout their lifetime.

Our global network of technical experts stands ready to assist with all system support needs. We provide customer support on-line, on the phone and on-site, so laboratories can maintain increased efficiency for years to come.

For more information on how a sample preparation system can improve the accuracy and efficiency of your lab's flow cytometry processes, contact your sales representative or visit www.beckmancoulter.com/cytomics today.



Simplify • Automate • Innovate

Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707. Australia, Gladstoneville (61) 2 9844 6000. Canada, Mississauga (1) 905 819 1234. China, Beijing (86) 10 6515 6028. Czech Republic, Prague (420) 1267 00 83 66. Hong Kong (852) 2814 7431, 2814 0481. France, Villepinte (33) 1 49 90 90 00. Germany, Krefeld (49) 2151 33 35. India, Mumbai (91) 22 3080 5101. Italy, Cassina de' Pecchi (Milano) (39) 02 953921. Japan, Tokyo (81) 3 6745 4704. Latin America (1) (305) 380 4709. Mexico, Mexico City (52) 55 560 57770. Netherlands, Mijdrecht (31) 297 230630. Puerto Rico (787) 747 3335. Singapore (65) 6339 3633. South Africa/Sub-Saharan Africa, Johannesburg (27) 11 805 2014. Sweden, Bromma (46) 8 564 85 900. Switzerland, Nyon 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) 714 953 5321. B2006-7430-LL-1.5K www.beckmancoulter.com © 2007 Beckman Coulter, Inc. PRINTED IN U.S.A.

BR-10281A

Enhanced Efficiency And Quality With An Increasing Workload

Standardized Sample Preparation Systems



Simplify • Automate • Innovate

Standardize And Improve Productivity From Start To Finish

When you need a more efficient and cost-effective way to manage your laboratory's flow cytometry sample preparation, turn to the world leader in progressive automation – Beckman Coulter. We offer not just one, but four state-of-the-art solutions that can meet your laboratory's needs and improve your entire flow cytometry process.

Whether your lab needs the cost-effective Q-Prep workstation, the automated benefits of the TQ-Prep and PrepPlus2 workstations or the integration of FP 1000 Cell Preparation System, Beckman Coulter has a solution that meets your sample preparation needs perfectly – and the expertise in automation that remains unmatched in the industry.

Achieve complete automation of pre-analytical processes

FP 1000 Cell Preparation System.

By combining the speed and efficiency of the PrepPlus 2 and the accuracy of the TQ-Prep into a single system, the FP 1000 delivers a whole new level of capability to the flow cytometry laboratory. Bar-code labeled specimens and daughter tubes are tracked and verified through the sample preparation process, reducing errors, re-runs and incorrect sample ID.

FP 1000 benefits:

- Full bi-directional LIS connectivity eliminates transcription errors and automates data transmission
- Streamlined workflow with electronic worklists for up to 4 FC 500 analyzers
- Improved sample processing efficiency - greater than 80 daughter tubes per hour (batch mode with tetraCHROME™ reagents)
- Cap piercing probe minimizes biohazard exposure and decreases occurrence of repetitive motion injuries
- Enhanced workflow with continuous processing - load and unload daughter carousels and samples without halting operation
- Decrease pre-analytical labor requirements - automate highly repetitive setup tasks including antibody/sample pipetting and lysis steps
- Full range of on-board reagents including, antibodies, lysing reagent, Flow-Count™, buffers and diluents

Standardize whole blood lysis, time after time

COULTER® TQ-Prep™ Workstation.

In response to the increasing demands on laboratories to perform immunophenotyping and CD4 measurements, the TQ-Prep workstation is the ideal way for your lab to meet its needs. Its direct compatibility with the Cytomics™ FC 500 Series and Epics™ XL-MCL™ flow cytometers ensures efficiency and cost savings, while delivering reliable operation for years to come. It also features:

- True walk-away operation from on-board incubation and lysis cycle processing
- Optimized cost-effectiveness and labor efficiency by reducing the number of repeat samples
- User-friendly interface with an icon-driven touch screen
- Maximum reliability with minimal maintenance
- Continuous real-time feedback allowing operators to follow the incubation and lysing process
- Precise and accurate reagent dispensing from an advanced syringe pump system to achieve standardized, reproducible sample preparation
- Sample throughput of up to 32 samples in only 24 minutes



Automate sample processing, while enhancing efficiency and precision

COULTER® PrepPlus 2 Workstation.

For laboratories that require the speed and efficiency of automated sample processing, Beckman Coulter's PrepPlus 2 is an ideal solution. By connecting the PrepPlus 2 workstation to a TQ-Prep system, your laboratory can gain the benefits of a single, comprehensive system that has the ability to automate your entire sample preparation process—resulting in a whole new level of lab productivity.

The PrepPlus 2—controlled by the TQ-Prep through its touch-screen—provides precision pipetting of reagents, patient samples, controls and calibrators into secondary tubes. Lysing is then automatically performed on the TQ-Prep, with minimal operator intervention.

The PrepPlus 2 offers a variety of advanced features that help keep your laboratory on a smooth course of efficiency, including:

- Enhanced precision through automated processing
- Increased sample volume with throughput of up to 90 daughter tubes per hour
- Flexible software programming of reagents, controls and pipetting parameters
- Sample flexibility that provides the ability to run whole blood and body fluids
- Improved safety with closed tube sampling, from a wide range of primary tubes
- Maintained whole blood cell suspensions with rotating specimen racks