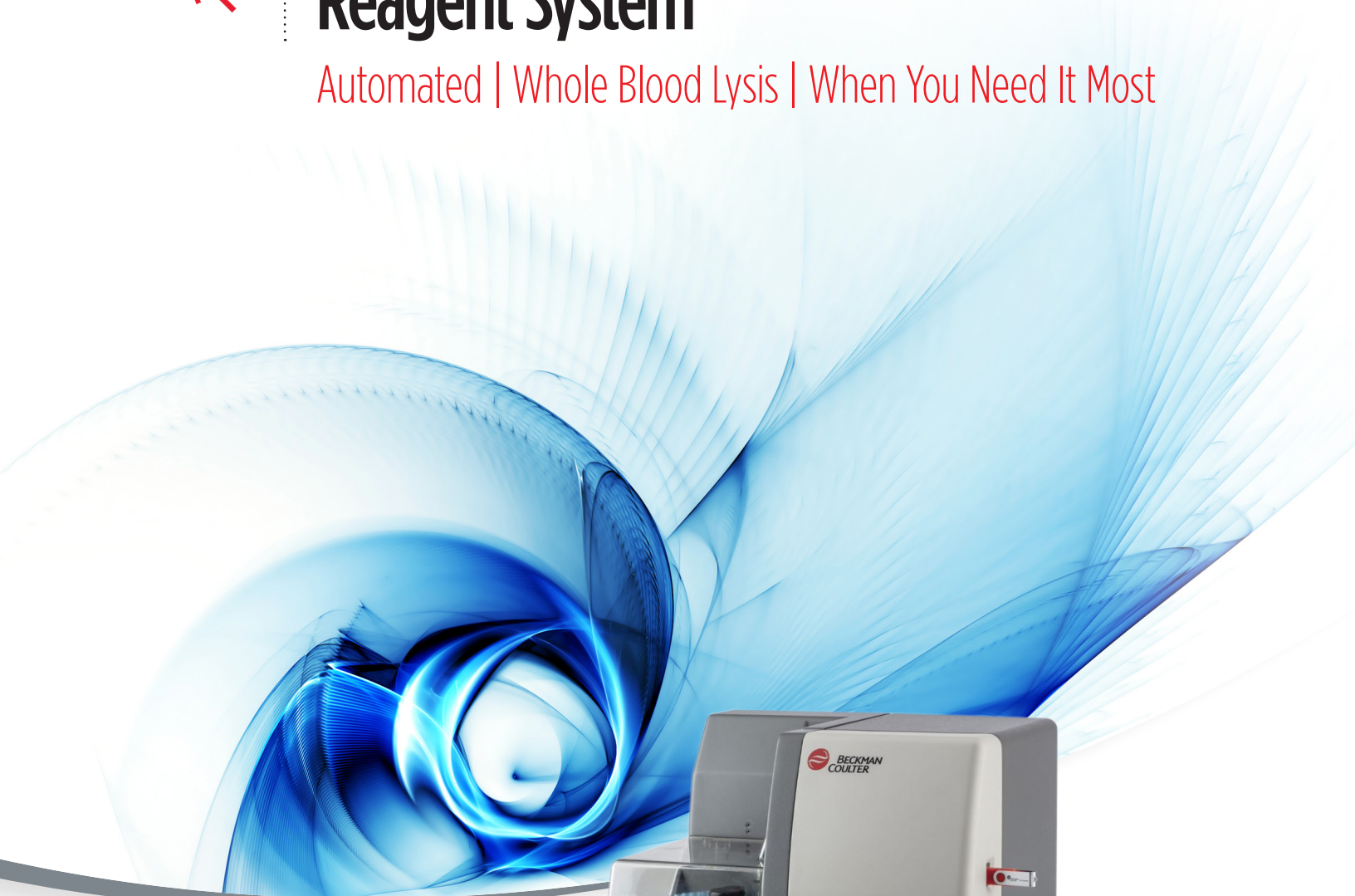




TQ-Prep Workstation & IMMUNOPREP Reagent System

Automated | Whole Blood Lysis | When You Need It Most



VISUALIZE
the possibilities.



The TQ-Prep Workstation and IMMUNOPREP Reagent System together provide rapid, no-wash whole-blood sample preparation. Automation provides inter-sample and inter-laboratory consistency and reproducibility. A standard 32-tube carousel allows walk-away processing and perfectly complements Beckman Coulter's Navios, Navios EX and DxFLEX flow cytometers from Beckman Coulter Life Sciences.

Proven Performance and Reliability

The IMMUNOPREP Reagent System with ready-to-use Cyto-Stat tetraCHROME Monoclonal Antibody Reagents provides fast no-wash sample preparations for flow cytometric analysis, while maintaining cellular integrity and morphology.

The TQ-Prep Workstation delivers whole blood lysis cell preparations for your flow cytometry needs time after time.

- True walk-away operation from on-board incubation and lysis cycle processing
- User-friendly interface with a larger icon-driven touchscreen
- Maximum reliability with minimal maintenance
- The vortex system mixing ensures reproducible homogenous cell lysis
- Reagent dispensing with syringe pump system consistently provides precise and accurate volumes
- Sample throughput of up to 32 samples in only 24 minutes
- Windows 10 IOT (Internet of things) Enterprise LTSC Operating System
- Data transfer via USB

TQ-Prep Upgraded Features

- Updated covers for front panel and reagent cover with improved fit
- Data privacy and security enhancements
- Updated safety certificates
- The touch screen interface has been replaced with a capacitive touch display interface
- Display icons have been resized for larger display



Beckman Coulter Life Sciences: 35 years of Innovations in SAMPLE PREPARATION

Sample Prep Automation has been provided by Beckman Coulter over the past 35 years



IMMUNOPREP Reagent System

The IMMUNOPREP Reagent System is a rapid, no-wash, standardized, whole blood preparation procedure maximizing sample-to-sample, and laboratory to laboratory reproducibility. The IMMUNOPREP reagent lysing action reduces cell stress and minimizes cell shrinkage while maintaining cellular integrity and morphology. Separation of leukocyte cell populations of interest from red blood cell ghosts and debris is maximized. Scatter characteristics are optimized for superior cell type resolution.



The IMMUNOPREP Reagent Kit is comprised of 3 ready-to-use reagents:

- Reagent A lyses the red blood cells
- Reagent B buffers the solution and stops the lysing process
- Reagent C fixes the cells

Performance Specifications

Accuracy of Delivery Volume of Distilled Water

- Volume delivery of syringe for Reagent A: 600 μ L \pm 5%
 - Lysing time: 8 seconds \pm 1 second
- Volume delivery of syringe for Reagent B: 265 μ L \pm 5%
 - Stabilizing time: 10 seconds \pm 1 second
- Volume delivery of syringe for Reagent C: 100 μ L \pm 5%
 - Fixative time: 10 seconds \pm 1 second

Performance Characteristics

- The precision (reproducibility) of the unit based on 32 replicate samples tested for CD4 have a coefficient of variation (CV) of 1.2%.

Order information

Product Name	Product Number
TQ-Prep Workstation	6605429
300 test IMMUNOPREP Reagent System	7546999



The device is not available for sale in all markets; please contact your Beckman Coulter representative for availability.

© 2023 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other product names and brands are properties of their respective owners.

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

2023-GBL-EN-101268-v2