

The MACSQuant® Analyzer 16

Colorful new possibilities in automated flow cytometry



Anous.



More colors, more applications

The MACSQuant[®] Analyzer 16 is engineered to expand the revolutionary automation of the MACSQuant Analyzer line of flow cytometers. It facilitates fully automated data acquisition in a 24-tube or 96-well plate format. The instrument's compact design is suited for basic research as well as advanced immune monitoring applications. The MACSQuant Analyzer 16 provides users with the flexibility and robustness that are required for the increasing demands of modern laboratories.

Automation

Load your samples, import your experiment settings, and leave the staining, data acquisition, and analysis to the MACSQuant Analyzer 16.

Expanded fluorescent channels

Amplify your data with a selection from 14 fluorescence channels and an average of 40% less sample volume required.

Simplicity

Minimize the learning curve with straightforward experiment setup and operation.

Multi-instrument alignment

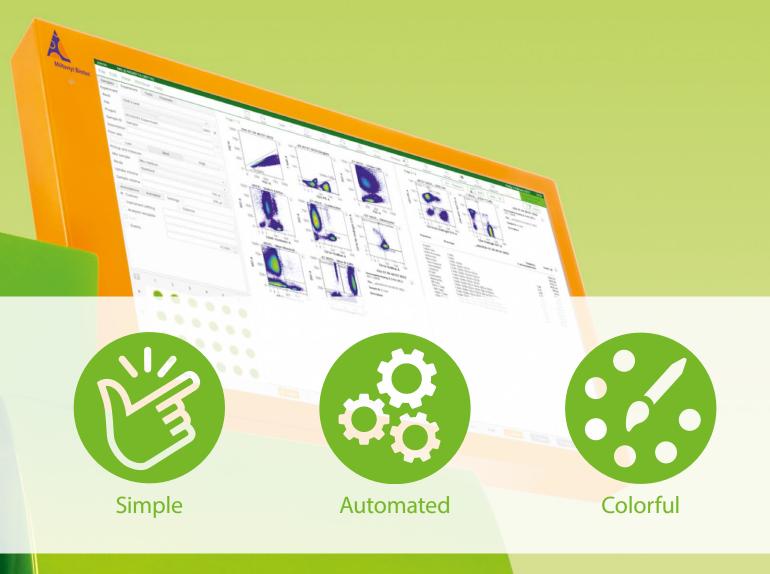
Using our Smart Gain software technology, users can harmonize data with collaborating labs to ensure reproducibility.

High quality

The MACSQuant Analyzer 16 was designed and developed within the established Quality Management System according to ISO 13485, providing you with an instrument of clinical quality.









Familiar small footprint with expanded fluorescence capabilities

As research questions become more complex, let the MACSQuant Analyzer 16 simplify your data acquisition. With expanded fluorescence capabilities, you can now maximize the number of parameters you can measure for each sample. Fluorescent channels are available with 405, 488, and 640 nm excitation lasers to cover a broad range of dyes and fluorescent proteins, enabling you to tell your story with more color. With the MACSQuant Analyzer 16, you can expand your immunophenotyping and functional applications to span a broad range of cell types, such as T cells, B cells, NK cells, dendritic cells, as well as tissue-specific, neural, and stem cell populations, and microorganisms. What would these new possibilities mean for your research?

Violat la car 105 vara	SSC	405/10 nm	
Violet laser 405 nm	V1	450/50 nm	VioBlue®/Vio® Bright V423
	V2	525/50 nm	VioGreen™
	V3	579/34 nm	BV 570
	V4	615/20 nm	Vio® Bright V600
	V5	667/30 nm	BV 650

	FSC	488/10 nm	
Blue laser 488 nm	B1	525/50 nm	FITC/Vio [®] Bright B515
	B2	579/34 nm	PE
	B3	615/20 nm	PE-Vio [®] 615
	B4	667/30 nm	PerCP
	B5	725/40 nm	PerCP-Vio [®] 700
	B 6	785/62 nm	PE-Vio [®] 770

Red laser 640 nm

R1	667/30 nm	APC/Vio [®] Bright R667
R2	725/40 nm	Vio® Bright R720
R3	785/62 nm	APC-Vio [®] 770

The features that make the difference

Large monitor

Integrated 15.6" monitor for simple touchscreen operation

Robotic needle arm

At the heart of automated sample mixing, processing, and autolabeling

Integrated MACS® Magnetic Cell Enrichment

Focus on your cells of interest and fortify your data

Precision volume uptake syringe

True volumetric cell counting without counting beads

Universal Reagent Rack

Flexible autolabeling options from glass or plastic vials

MACS MiniSampler Plus

Seamless sampling from single tubes, multiple tubes, or 96-well plates





Automated flow cytometry

The MACSQuant Analyzer 16 offers a plethora of built-in automation features to reduce hands-on time and operator variability.

Hassle-free housekeeping

Automated start-up, calibration, cleaning, and shutdown ensure instrument robustness and high-quality data without cumbersome preparation.

Autolabeling

The missing step in a fully automated flow cytometry assay. Reduce the risk of pipetting error with the reliability of robotics using the automated pipetting arm, the MACS MiniSampler Plus, and our newly designed Universal Reagent Rack.

Integrated magnetic cell isolation

Using the integrated MACS Enrichment Column, you can magnetically enrich your target population to perform a deeper analysis of rare cell subsets. Removing the non-relevant events before your flow analysis makes your assay even more robust.

Automated compensation

Get rid of time-consuming manual compensation and let the MACSQuant Analyzer 16 do it for you.

Automated analysis

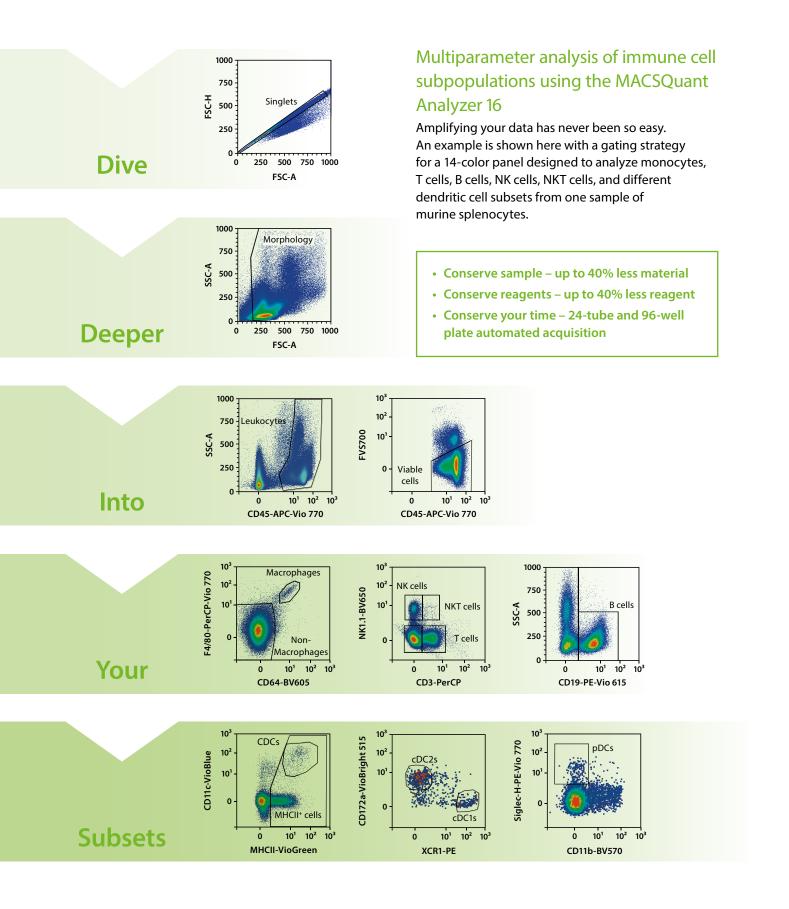
Use our Express Modes, unique add-ons for the MACSQuantify[™] Software, to automate measurement and analysis of your flow experiments ensuring simplicity and reproducible data analysis.





Get more with the MACSQuant Analyzer 16

More channels. More data. More subsets.





Harmonizing the world of flow cytometry is just an instrument setting away

Experience consistently reproducible results from day to day, instrument to instrument, and operator to operator. With the Smart Gain technology of the MACSQuantify Software, you can transfer your assay from one instrument to another, while passing on all necessary information to ensure an identical setup.



Your partner in a regulated environment

Further characterize and enumerate your cells in your cell manufacturing applications. The MACSQuant Analyzer 16 hardware and software are configured to support small- to large-scale manufacturing requirements and offer tools and support for GMP-compliant processes.

Compliance with 21 CFR Part 11

You can be sure that your data is suitable for submission to regulatory agencies with the optional 21 CFR Part 11 feature, which includes:

- Audit trail
- Analysis reports with e-signature
- User management system according to 21 CFR Part 11

Quality Management System according to ISO 13485

The MACSQuant Analyzer 16 was designed and developed within the established Quality Management System according to ISO 13485, providing you with an instrument of clinical quality.







Complete your automation loop with reliable reagents

In order to achieve reproducible and high-quality results, Miltenyi Biotec offers a complete flow cytometry solution including a dedicated range of reagents. We help you make sure that variations in your experiment are due to your sample and not due to unreliable antibodies or instruments.

REAfinity[™] Recombinant Antibodies – flow cytometry is in their genes

Miltenyi Biotec has introduced a portfolio of REAfinity Recombinant Antibodies that provide superior lot-to-lot consistency and purity compared to mouse or rat monoclonal and polyclonal antibodies. They diminish the need for FcR blocking and allow for analyses with only one single isotype control. They help you to generate high-quality data with no more background signal while saving efforts when setting up experiments.

Advantages of REAfinity Recombinant Antibodies:

- High lot-to-lot consistency
- One universal isotype
- No more background signal





Vio Dyes – brighter dyes for flow cytometry

When used in combination with our proprietary Vio and Vio Bright Dyes, REAfinity Antibodies provide superior mean fluorescence intensity and high stain indices. With the brightest dyes on the market, setting up complex multicolor experiments has never been so simple.

Ready-to-use kits

Use Miltenyi Biotec's range of ready-to-use, pretitrated kits and save valuable experiment set-up time and assay costs. Our kits have been validated to be used with the automatic labeling function of the MACSQuant Analyzer 16, which in combination with our Express Modes, give you true walk-away capacities. All you have to do is set up the experiment and come back to look at completely analyzed data.

Miltenyi Biotec Tested Panels (MBTPs)

Miltenyi Biotec also provides a list of pre-tested and validated antibody panels complete with gating strategy, staining protocols, and reagent list for easy and accurate flow cytometric analysis.

Customized solutions

Miltenyi Biotec's Custom Antibody Design Service enables researchers to benefit from personalized flow cytometry solutions. This service includes purified, functional-grade antibodies, single- and multicolor antibody conjugates, as well as multicolor antibody cocktails.



For more information about our products, visit:

- miltenyibiotec.com/reafinity
- miltenyibiotec.com/vio
- miltenyibiotec.com/MBTPs
- miltenyibiotec.com/customab



MACSQuant Analyzer 16 specifications

Optics					
Laser excitation	Spatially separated: 405 nm, 65 mW 488 nm, 50 mW 640 nm, 72 mW				
Emission detectors	FSC: 488/10 nm SSC: 405/10 nm	V1: 450/50 nm V2: 525/50 nm V3: 579/34 nm V4: 615/20 nm V5: 667/30 nm	B1: 525/50 nm B2: 579/34 nm B3: 615/20 nm B4: 667/30 nm B5: 725/40 nm B6: 785/62 nm	R1: 667/30 nm R2: 725/40 nm R3: 785/62 nm	
Fluorescence sensitivity and resolution	MESFs (CV <5%): FITC <110 PE <75 APC <100				
Scatter resolution	Scatter performance is optimized for resolving human peripheral blood lymphocytes, monocytes, and granulocytes				
Flow cell dimensions	$200 imes 250 \ \mu m$				
Minimal guaranteed particle size	500 nm				
Fluorescence detectors	Optimized with spectrally matched PMTs for all channels				
Laser spot size	15 × 45 μm				
Fluidics					
Minimal uptake volume ¹	1 μL (25 μl recommended for volumetric counting applications)				
Sample flow rate	25, 50, or 100 μ L/min or automate flow rate to maintain 500, 1,000, or 2,000 events/second				
Measurement speed ^{2,3}	<25 minutes per 96-well plate (5 µL measurement volume per well)				
Sample uptake	1–450 μL				
Maximal event rate	Up to 15,000 events/second				
System maintenance	Automated startup, PMT calibration, cleaning cycles, and shutdown				
Sample mixing	Aspiration				
Performance					
Absolute counts performance ^{2,4}	CV <7%				
Sample carryover ^{2,5}	0.01%				
Sample tube/plate	1.5 mL, 2 mL and 5 mL single tubes, up to 24 \times 5 mL single tubes, up to 96-well plate (U, V, flat bottom; deep-well)				
MACS Cell Enrichment Unit	For pre-analysis enrichment of rare cells				

¹ At every uptake, an additional excess volume is aspirated by the instrument. The excess volumes are calibration- and process-dependent and do not exceed ¹ 0 μL for fast, Standard, and Extended modes, and 20 μL for Screen mode.
² Referred value indicates the average of multiple experiments and can differ for individual sample materials.

 ² Referred value indicates the average of multiple experiments and can differ for individual sample materials.
³ The measurement speed is determined by measuring the time between the movement of the robotic needle arm into the first measured well, and its movement out of the last measured well. The measurements were carried out at High flow rate in Screen mode.
⁴ For counting performance, full 96-well plates were loaded with 200 µL/well of peripheral blood mononuclear cell (PBMC) suspension at a nominal concentration of 5,000 cells/µL. The uptake volume was set to 50 µL at Medium flow rate.
⁵ For carryover, full 96-well plates were loaded with 200 µL/well of PBMC suspension at a nominal concentration of 10,000 cells/µL in every other well ("SRC-wells"). Alternating wells were loaded with an equal volume of MACSQuant Running Buffer ("CO-wells"). The uptake volume was set to 100 µL and mean at Medium flow rate in Strand at Medium flow rate in Strand mode. measured at Medium flow rate in Standard mode. The carryover is defined by sum (CO-singlet count)/sum(SRC-singlet count) × 100%.

Support at your fingertips

MACSQuant Live Support

- Live support at your fingertips via MACSQuant Support portal
- Have your questions answered in real time by one of our experts

For more information, please visit:

- miltenyi.com/support
- miltenyi.com/customapplication-service
- miltenyi.com/training

Application and instrument support

- Technical and field application support for assay design and product advice
- Custom automation and express mode development

Instrument training

- Training at regional Miltenyi Innovation and Training Centers (MITC)
- Onsite training and assay development
- Online application resources

Service

- Comprehensive service options
- Globally distributed field service teams

Miltenyi Biotec – a company supporting your complete workflow

Miltenyi Biotec offers a complete portfolio of research tools that enables a stream-lined workflow, from reproducible sample preparation to a variety of downstream applications. MACS Solutions are the perfect companion for every step of your inspired research.



MACS Sample Preparation

The secret to the success of any experiment depends on the quality of the starting material. Our sample preparation portfolio has the tools you need to start smart. Our innovative instruments and reagents help you standardize your tissue preparations and get reproducible data.



MACS Cell Separation

Whether isolating cells in small-scale experiments or in high-throughput industrial settings – we offer manual, semi-automated, automated, and robotic integration solutions to meet your specific research demands.



MACS Antibodies

Miltenyi Biotec offers a huge range of fully validated antibodies suitable for multiple applications, such as flow cytometry, microscopy, and functional assays. The MACS Antibody Portfolio guarantees high performance consistency with innovative and proprietary recombinant antibody technology.



MACS Flow Cytometry

MACS Flow Cytometry provides best-in-class solutions for all your research needs. Instruments, reagents, kits, and software constitute a comprehensive portfolio to keep your finger on the pulse of advanced flow cytometry, cell sorting, and cell analysis.



MACS Imaging and Microscopy

Our growing MACS Imaging and Microscopy portfolio offers versatile, high-quality imaging solutions to study complex biological systems.



MACS Cell Culture and Stimulation

Our MACS Cell Culture and Stimulation portfolio comprises a specialized and versatile range of cell culture media and reagents for cell stimulation, expansion, and/or differentiation.



MACS Molecular Analysis

MACS Technology has been perfectly adapted for molecular applications for fast and sensitive analysis of cells at any subcellular level. Explore our products for high-yield mRNA, protein, and organelle isolation, cDNA synthesis, and stable cellular transfection.

miltenyibiotec.com/mq16



Germany/Austria

Miltenyi Biotec B.V. & Co. KG Friedrich-Ebert-Straße 68 51429 Bergisch Gladbach Germany Phone +49 2204 8306-0 Fax +49 2204 85197 macsde@miltenyi.com

USA/Canada

Miltenyi Biotec, Inc. 2303 Lindbergh Street Auburn, CA 95602, USA Phone 800 FOR MACS Phone +1 866 811 4466 Fax +1 877 591 1060 macsus@miltenyi.com

Australia

Miltenyi Biotec Australia Pty. Ltd. Unit 11, 2 Eden Park Drive Macquarie Park, NSW 2113 Australia Phone +61 2 8877 7400 Fax +61 2 9889 5044 macsau@miltenvi.com

Benelux

Miltenyi Biotec B.V. Dellaertweg 9C 2316 WZ Leiden The Netherlands macsnl@miltenyi.com Customer service for: The Netherlands Phone 0800 4020120 Fax 0800 4020100

Belgium Phone 0800 94016 Fax 0800 99626 Luxembourg Phone 800 24971 Fax 800 24984

China

Miltenyi Biotec Technology & Trading (Shanghai) Co., Ltd. Room A401, 4/F No. 1077, Zhangheng Road Pudong New Area 201203 Shanghai, P.R. China Phone +86 21 6235 1005-0 Fax +86 21 6235 0953 macscn@miltenyi.com.cn

France

Miltenyi Biotec SAS 10 rue Mercoeur 75011 Paris, France Phone +33 1 56 98 16 16 macsfr@miltenyi.com

Hong Kong Miltenyi Biotec Hong Kong Ltd. Unit 301, Lakeside 1 No. 8 Science Park West Avenue Hong Kong Science Park Pak Shek Kok, New Territories Hong Kong Phone +852 3751 6698 Fax +852 3619 5772 macshk@miltenyi.com.hk

Italy

Miltenyi Biotec S.r.l. Via Paolo Nanni Costa, 30 40133 Bologna Italv Phone +39 051 6 460 411

Fax +39 051 6 460 499 macsit@miltenyi.com

Japan

Miltenyi Biotec K.K. NEX-Eitai Building 5F 16-10 Fuyuki, Koto-ku Tokyo 135-0041, Japan Phone +81 3 5646 8910 Fax +81 3 5646 8911 macsjp@miltenyi.com

Nordics and Baltics

Miltenyi Biotec Norden AB Medicon Village Scheeletorget 1 223 81 Lund Sweden macsse@miltenyi.com **Customer service for:** Sweden Phone 0200 111 800 Fax +46 280 72 99 Denmark

Phone 80 20 30 10 Fax +46 46 280 72 99 Norway, Finland, Iceland, and Baltic countries Phone +46 46 280 72 80 Fax +46 46 280 72 99

Singapore

Miltenyi Biotec Asia Pacific Pte Ltd. 438B Alexandra Road, Block B Alexandra Technopark #06-01 Singapore 119968 Phone +65 6238 8183 Fax +65 6238 0302 macssg@miltenyi.com

South Korea

Miltenyi Biotec Korea Co., Ltd. Donggeuk 7F, 562 Nonhyeon-ro Gangnam-gu Seoul 06136, South Korea Phone +82 2 555 1988 Fax +82 2 555 8890 macskr@miltenyi.com

Spain

Miltenyi Biotec S.L. C/Luis Buñuel 2 Ciudad de la Imagen 28223 Pozuelo de Alarcón (Madrid) Spain Phone +34 91 512 12 90 Fax +34 91 512 12 91 macses@miltenyi.com

Switzerland

Miltenyi Biotec Swiss AG **Gibelinstrasse 27** 4500 Solothurn Switzerland Phone +41 32 623 08 47 Fax +49 2204 85197 macsch@miltenyi.com

United Kingdom

Miltenyi Biotec Ltd. Almac House, Church Lane Bisley, Surrey GU24 9DR, UK Phone +44 1483 799 800 Fax +44 1483 799 811 macsuk@miltenyi.com

www.miltenyibiotec.com

30-118-919.03

Miltenyi Biotec provides products and services worldwide. Visit www.miltenyibiotec.com/local to find your nearest Miltenyi Biotec contact.

Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for therapeutic or diagnostic use. MACS, MACSQuant, MACSQuantify, REAfinity, Vio, VioBlue, VioGreen, and the Miltenyi Biotec logo are registered trademarks or trademarks of Miltenyi Biotec and/or its affiliates in various countries worldwide.