



The first built-for-purpose nanoscale flow cytometer capable of detecting nanoparticles at least as small as 40 nm* while simultaneously characterizing them in detail.



EXCITATION OPTICS

The CytoFLEX nano flow cytometer has the capacity for 12 simultaneous detection channels, broken out as 6 fluorescent channels, 5 side scatter channels and 1 forward scatter channel. Each instrument is available in one configuration with all lasers activated. The excitation lasers are listed in the table below.

LASER SPECIFICATIONS Colinear Laser Description

Laser	Wavelength	Power at Flow Cell
Violet	405 nm	120 mW
Blue	488 nm	50 mW
Yellow	561 nm	35 mW
Red	638 nm	100 mW

FLOW CELL

Fixed integrated optics and quartz flow cell design

FORWARD SCATTER DETECTION

Silicon photodiode with built-in 405/10 nm bandpass filter

BANDPASS FILTERS

Side scatter filters cannot be replaced, repositioned or removed. Fluorescence filters can be replaced, but can not be repositioned or removed. Please discuss custom filter placement with technical support.

Fluorescence	Scatter
447/60	405/10 (3) - VSSC1, VSSC2, VFSC
531/46	488/8 - BSSC
595/50	561/6 - YSSC
670/30	638/6 - RSSC
710/47	
792/64	

FLUORESCENCE AND SIDE SCATTER DETECTION

Fluorescent and side scatter light is delivered by fiber optics to Avalanche Photodiode detector arrays. Proprietary design ensures high performance, high efficiency, low-noise signal detection. Emission profiles are collected using reflective optics and single transmission bandpass filters.



QUALITY CONTROL

For detection channels off each of the lasers, CytExpert nano QC fluorescence pass/fail criteria is rCV <10% when using CytoFLEX nano QC Fluorospheres and QC Scatterspheres. QC process is enhanced with automated baseline monitor process.

FLUIDICS

Built-in diaphragm and maintenance-free piston pumps deliver both sheath and sample, respectively.

FLUIDIC CAPACITY

Fluidic cart with on-board fluid containers

10 L disposable sheath container

5 L disposable CytoFLEX Cleaner

SAMPLE FLOW RATES

Adjustable sample flow rates from approximately 1-6 μ L/min, in 1 μ L/min steps.

SAMPLE INPUT FORMATS

5 mL (12 x 75 mm) tubes, 1.5 mL micro-centrifuge sample tubes

VOLUMETRIC COUNTING ACCURACY

>90% - tested with 144nm QC Scatterspheres at all nominal sample flow rates.

MAINTENANCE

Automated sample and sheath baseline signal check and user defined cleaning and acceptance criteria.

Maintenance reminder can be set up in the software for some regular operations with custom cycle tim.e

User replaceable fluidic filters and sample line and probe

ELECTRONICS

NOMINAL ELECTRONIC ACQUISITION RATE

16,000 events per second with all configured parameters

RECOMMENDED ACQUISITION RATE

Up to 5,000 events per second with all configured parameters

DATA DISPLAY

Fully digital system with 7 decades data display

SIGNAL

Pulse area, height for every channel, width for one selectable channel

PERFORMANCE

ANALYSIS

SCATTER SENSITIVITY

Reflective optics with a single transmission bandpass filter in front of each detector

Violet (405 nm) Side Scatter Resolution (VSSC1, VSSC2): Able to detect scattered light from polystyrene particles at least as small as 40 nm* and as large as 1000 nm. Intended range of operation: VSSC1 - 40 nm to 150 nm, VSSC 2 - 80 nm to 1000 nm, when light scatter is equivalent to or greater than polystyrene.

CARRYOVER

<1.0% for both QC beads and biological material

FLUORESCENCE SENSITIVITY

Sensitivity is monitored by automated QC process using 500 nm CytoFLEX nano Multi-intensity Fluorospheres

DATA MANAGEMENT

SOFTWARE

The CytExpert nano software is a full-featured software package that controls the instrument's operation, collection of experiment data, and analysis of the results

If desired, data files can be exported as FCS files for analysis in Kaluza, Cytobank and other platforms

LANGUAGE

English and Chinese

OPERATING SYSTEM

Windows® 10 Enterprise LTSC 2019 x64 -bit

FCS FORMAT

FCS 3.0, FCS 3.1

MAXIMUM DATA FILE SIZE

25 million event per file with all parameters

RECOMMENDED MINIMUM COMPUTER SPECIFICATIONS

CPU: 10th Gen Intel $^{\circ}$ Core $^{\text{TM}}$ i7 (12MB Cache, 2.40 GHz)

Memory: 16 GB RAM or higher

Storage: 512 GB TB

Ethernet: Integrated 100M GB, Dual Ethernet ports

USB Port: >4 ports

Monitor: 32-inch monitor (2560x1440 resolution) or 24-inch

monitor (1920x1080 resolution)

COMPENSATION

Automatic full matrix compensation

Manual full matrix compensation

Novel Compensation Library: store fluorescent spillover values of dyes to easily determine the correct compensation matrix with new gain settings

Import/export compensation values or matrix between experiments

Absolute linear gain amplifications allows automatic compensation adjustment between experiments that use different gain settings

INSTALLATION REQUIREMENTS

DIMENSIONS (W X D X H)

FLOW CYTOMETER

59 cm x 50 cm x 44 cm

FLUIDICS CART

37 cm x 89 cm x 33 cm, cart can be placed on bench or floor

WEIGHT

CytoFLEX nano Flow Cytometer: 45 kg

Fluidics cart (with fluid): 22 kg

POWER SPECIFICATIONS

Voltage: AC 100-240V, 50/60Hz

Rated Power: 200VA

OPERATION ENVIRONMENT

Ambient Temperature: 15-27°C within ±2°C temperature variation during operation

Relative Humidity: 20-80% (noncondensing)

Altitude: 2,000 m (max)

COMPLIANCE AND SAFETY STANDARDS

EN 61010-1:2010+A1:2019

UL 61010-1:2012

CAN/CSA-C22.2 NO. 61010-1-12 + GI1 +GI2 (R2017) + A1

IEC 61010-2-081:2019

EN IEC 61010-2-081:2020

UL 61010-2-081:2019

CSA C22.2 NO. 61010-2-081:19 61010-1:2010, AMD1:2016

AS/NZS CISPR 11:2017+AMD 1:2020

ENVIRONMENTAL, HEALTH, AND SAFETY

EU RoHS 3(2015/863/EU)

EU waste electrical and electronic equipment 2012-19/EU

California Proposition 65

REACH 1907/2006

Packaging and Packaging Waste Directive 94/62/EC

California Seismic Regulation LAW

COMPLIANCE CERTIFICATION MARKS

cNTRLus Certification Mark

CB Certification Mark

CE Certification Mark

RCM Certification Mark

UKCA Certification Mark

*polystyrene when triggering on violet side scatter



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

© 2024 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.

