



CFX Duet Real-Time PCR System

Powerful Two-Color Genomics Analysis, Accessible for Most Laboratories



The CFX Duet Real-Time PCR System makes uniform and precise quantitative PCR (qPCR) accessible for any lab. The CFX Duet provides the thermal performance and the proprietary, accurate optical shuttle system of our premium CFX Opus Real-Time PCR Systems for customers that run only singleplex or duplex assays.

The CFX Duet System utilizes CFX Maestro Software through an attached computer as its control and analysis system. The economical user can quickly get started with gene expression, genotyping, thermal gradient optimization, quantification, and even protein characterization using fluorescence resonance energy transfer (FRET). The two-color CFX Duet System is factory calibrated for common dyes such as SYBR® Green and EvaGreen®. This system is not compatible with the Br.io cloud platform.

Benefits

- Easy entry point into basic qPCR with 2-color system
- No calibration or passive reference dyes required
- Versatile capabilities for essential molecular biology and protein characterization
- Additional FRET LED channel, enabling protein melt analysis

- Time-saving experiment optimization with thermal gradient feature
- Optimal thermal performance equivalent to the premium CFX Opus Systems
- Continued support for and control of up to 4 CFX instruments with CFX Maestro Software, and Security Edition

Specifications

System

Sample block type	Fixed
Sample capacity	96 wells
Sample size	1–50 µl (10–50 µl recommended)
Dimensions (W x D x H)	33 x 56 x 36 cm (13 x 22 x 14 in.)
Weight	22 kg (48 lb)
Communications	USB 2.0 or above
Electrical approvals	IEC, CE

Optical Detection System

Excitation	3 filtered LEDs
Detection	3 filtered photodiodes
Range of excitation/emission wavelengths	450–580 nm
Scan time	
All channels	12 sec
Single channel	3 sec
Dynamic range	10 orders of magnitude
Sensitivity	Detects 1 copy of target sequence in human genomic DNA
Multiplex analysis	Up to 2 targets per well

continues

Specifications

96-Well Reaction Block

Heating and cooling method	Peltier
Lid heating	30–110°C
Temperature range	4–100°C
Maximum ramp rate	5°C/sec
Average ramp rate	3.3°C/sec
Thermal accuracy	±0.2°C
Thermal uniformity	±0.3°C (min – max ≤ 0.6°C), measured 10 sec after the block reaches the target temperature

Gradient	
Operational range	30–100°C
Programmable span	1–24°C

CFX Maestro Data Management and Analysis Software

Operating system	Windows 10 (64-bit)
Memory	Minimum of 1 GB
Data analysis modes	<p>PCR quantification with standard curve</p> <p>Melt curve analysis</p> <p>Gene expression analysis by relative quantity (ΔCq) or normalized expression ($\Delta\Delta Cq$)</p> <p>Multiple-file gene expression analysis for comparison of an unlimited number of quantification cycle (Cq) values</p> <p>Allelic discrimination</p> <p>Endpoint analysis</p> <p>Data analysis options include bar charts, box-and-whisker plots, dot plots, clustergrams, scatter plots, and volcano plots</p> <p>Statistics include <i>t</i>-test and analysis of variance (ANOVA)</p>
Data export	<p>Save, copy, and print all graphs and spreadsheets from right-click menu</p> <p>Export specified data in multiple formats</p> <p>Copy and paste into a Microsoft Excel, Word, or PowerPoint file</p> <p>Print directly or save as PDFs customizable reports containing run settings, data graphs, and spreadsheets</p> <p>Export to RDML</p>
Image export	<p>Export images at any pixel size and at a resolution up to 600 dpi</p> <p>Save images as .bmp, .jpg, or .png files</p>

Ordering Information

Catalog #	Description
Instrument	
12012625	CFX Duet Real-Time PCR System , includes power cord, USB cable
Software	
	All software for Windows unless otherwise noted.
12013758	CFX Maestro Software , includes USB installation drive, video quick guides, PDF instruction manual, license for qbase+ Software
12012832	CFX Maestro Software, Security Edition , 1 license
12013028	CFX Maestro Software, Security Edition , 5 licenses
12012833	CFX Maestro Software, Russian Edition
12012834	CFX Maestro Software, Chinese Edition
1845025	Precision Melt Analysis Software , includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit
Consumables	
HSP9655	Hard-Shell 96-Well PCR Plates , pkg of 50, low profile, thin wall, skirted, white shell/white wells
HSP9601	Hard-Shell 96-Well PCR Plates , pkg of 50, low profile, thin wall, skirted, white shell/clear wells
MSB1001	Microseal 'B' PCR Plate Sealing Film , pkg of 100, optically clear seals for PCR plates
TLS0851	0.2 ml 8-Tube PCR Strips without Caps , low profile, white
TLS0801	0.2 ml 8-Tube PCR Strips without Caps , low profile, clear
TCS0803	0.2 ml Flat PCR Tube 8-Cap Strips , optical, ultraclear
Reagents	
1845098	CFX Qualification Plate , 96-well
12010176	Reliance One-Step Multiplex RT-qPCR Supermix , 1 ml (1 x 1 ml), 200 x 20 μ l reactions
1725270	SsoAdvanced Universal SYBR® Green Supermix , 2 ml (2 x 1 ml), 200 x 20 μ l reactions
1725016	SsoAdvanced Universal Inhibitor-Tolerant SYBR® Green Supermix , 2 ml (2 x 1 ml), 200 x 20 μ l reactions
1725095	SingleShot SYBR® Green One-Step Kit for Cell Lysis and RT-qPCR , 100 x 50 μ l reactions
17005726	SEQuoia Complete Stranded RNA Library Prep Kit , 24 reactions
12011928	SEQuoia Dual Indexed Primers Set , 12 vials of unique dual indexed primers, 96 reactions

Search **CFX qPCR Systems** in the Apple App Store and Android Play Store for an augmented reality tour. Contact your local sales representative for an in-person demonstration.

Visit bio-rad.com/CFXDuet for more information.

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. EvaGreen is a trademark of Biotium, Inc.

SYBR is a trademark of Thermo Fisher Scientific Inc. Bio-Rad Laboratories, Inc. is licensed by Thermo Fisher Scientific Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

