

Consistent and Accessible Quantitative PCR Analysis for Any Laboratory

Powerful Genomic Analysis Platform for Singleplex and Duplex Assays

The Bio-Rad™ 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 who run only singleplex and duplex assays.

The CFX Duet System utilizes CFX Maestro Software through an attached computer as its control and analysis system. CFX Maestro Software makes experiment setup and analysis easy and provides customized graphing and reports without exporting to other programs. 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

- No calibration or passive reference dyes required
- Intuitive and robust software
- Versatile capabilities for essential molecular biology and protein characterization
- Time-saving experiment optimization with thermal gradient feature
- Optimal thermal performance equivalent to the premium CFX Opus Systems

Ready Out of the Box

Our consistent distance scanning shuttle technology comes factory calibrated for common fluorophores and does not require passive reference dyes, so even labs new to qPCR will be ready to go upon receiving the system.



Ease of Use

The CFX Duet System uses the same software as our premium CFX Opus Systems. CFX Maestro Software makes setup and analysis easy. Postanalysis data management can be done within the software, which makes customized graphing and basic statistical analysis available without having to export files.

Essential Molecular Analysis Is Accessible

Since this two-color system does not require passive reference dyes, it can perform genotyping and many gene expression analyses. Quantification of up to two targets is possible and an additional FRET mode enables protein melt analysis for basic protein characterization experiments.

Efficient Optimization and Thermal Management

The CFX Duet System has thermal gradient functions across the eight rows of the thermal block. Even with this two-channel system, you can optimize thermal parameters in fewer runs. The CFX Duet System has the same thermal uniformity and precision as our premium CFX Opus Systems, so every lab can have access to its robust thermal capabilities.

Ordering Information

Catalog #	Description
Instrument	
12016265	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
12012832	CFX Maestro Software, Security Edition , 1 license
12013028	CFX Maestro Software, Security Edition , 5 licenses
12012833	CFX Maestro Software, Russian Edition , 1 license
12012834	CFX Maestro Software, Chinese Edition , 1 license
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 µl reactions
1725270	SsoAdvanced Universal SYBR® Green Supermix , 2 ml (2 x 1 ml), 200 x 20 µl reactions
1725016	SsoAdvanced Universal Inhibitor-Tolerant SYBR® Green Supermix , 2 ml (2 x 1 ml), 200 x 20 µl reactions
1725095	SingleShot SYBR® Green One-Step Kit for Cell Lysis and RT-qPCR , 100 x 50 µl 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](https://www.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. © 2023 Bio-Rad Laboratories, Inc.

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](https://www.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** 82 080 007 7373 **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

