

EONIS real-time PCR



Introduction

EONIS™ Q is a real-time PCR instrument with possibility to integrate into automation systems. The instrument is available in both 96- and 384-well formats. The EONIS™ Q has been designed to be robust, easy-to-use, affordable and meet the safety requirements of local regulations in target markets.

Product numbers	
EONIS™ Q96	2044-0020
EONIS™ Q384	2045-0020
PC	2044-8010



Hardware

Parameter	96 Well system	384 Well system
Sample size	5-100 µl (depending on consumables)	5-20 µl (depending on consumables)
Heating	Max. 5.5°C/s and ø 7°C/s	Max. 4°C/s and ø 3°C/s
Cooling	Max. 5.5°C/s and ø 4.5°C/s	Max. 2°C/s and ø 1.5°C/s
Adjustable temperature range	4°C to 99°C	4°C to 99°C
Max./Min. Gradient	40°C / 0.1 °C	24°C / 0.1 °C

Pc and software

Parameter	96 Well system	384 Well system
Control	PC with EONIS™ Q Software	
Operating system	Windows 10 or higher	

The software includes the following analysis methods:

- Absolute and relative quantification
- Delta-delta Ct-method
- Genotyping
- Allele discrimination
- PCR efficiency
- Melting curve (no HRM)
- Multi-gene and multi-plate analysis

Optics

Parameter	96 Well system	384 Well system
Measuring principal	Fibre optic shuttle system with 8fold scanner and color modules for excitation and emission filters	Fibre optic shuttle system with 16fold scanner and color modules for excitation and emission filters
Light source	4 combined LEDs RGBW	
Detector	PMT (Photo Multiplier)	

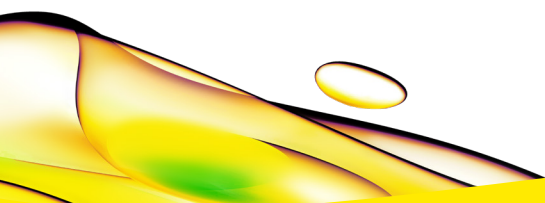
Parameters color modules (for 96 and 384 well)

Name	Excitation	Emission
Color module 1	465 nm \pm 15 nm	520 nm \pm 15 nm
Color module 2	510 nm \pm 15 nm	565 nm \pm 15 nm
Color module 3	530 nm \pm 15 nm	585 nm \pm 15 nm
Color module 4	560 nm \pm 15 nm	610 nm \pm 15 nm
Color module 5	625 nm \pm 10 nm	680 nm \pm 15 nm
Color module 6	625 nm \pm 10 nm	710 nm \pm 20 nm

Dimensions

Parameter	96 Well system	384 Well system
Weight	Approx. 38 Kg	
Dimensions (W x D x H)	310 mm x 345 mm x 613 mm	
Recommended footprint	310 mm x 477 mm (open drawer)	

For research use only. Not for use in diagnostic procedures.



revvity