

Compatibility cycler channels vs fluorescent dyes

Thermocycler	UV-A/Blue Channel	Channel 1	Channel 2	Channel 3	Channel 4	Channel 5	Channel 6	Channel 7
Real-Time PCR Thermocyclers								
AriaMx & AriaDx (up to 6 filters)	ATTO 425	FAM	YY*/HEX/JOE/ATTO 532*	ATTO 550*/TAMRA/Cy ³	ATTO 590*/ROX/TR	ATTO 647N/Cy ⁵		
qTOWER ³		FAM	YY/HEX/JOE/VIC	DFO/TAMRA/Alexa546/NED	ROX/TR/Cy ^{3.5}	ATTO 647N*/Alexa633/Cy ⁵	Cy ^{5.5}	
qTOWERiris	ATTO 425/ATTO 390*	FAM	YY/HEX/JOE/TET/VIC	ATTO 550*/TAMRA/DFO*	ATTO 590/ROX/TR/Cy ^{3.5}	ATTO 633/Cy ⁵	Cy ^{5.5} /ATTO 665	
ABI Prism [®] SDS 7500		FAM	YY*/HEX*/JOE/TET*/VIC	DFO*/TAMRA/Cy ³ /NED	ROX/TR	ATTO 647N*/HL647/Cy ⁵		
ABI Prism [®] SDS 7900HT		FAM	YY/JOE/TET/VIC	DFO*/TAMRA/NED	ROX			
QuantStudio™ 1		FAM	YY*/HEX/JOE/TET/VIC	ROX/TR/JUN				
QuantStudio™ 3		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy ³ *	ROX/ATTO 590*/TR*			
QuantStudio™ 5 (384-W.) & 6 Flex/Pro		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy ³ *	ROX/ATTO 590*/TR*	ATTO 647N*/HL647*/Cy ⁵		
QuantStudio™ 5 (96-W.) & 7 Flex/Pro & 12k Flex		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy ³ *	ROX/ATTO 590*/TR*	ATTO 647N*/HL647*/Cy ⁵	ATTO 680*/Cy ^{5.5} *	
StepOnePlus		FAM	YY*/HEX*/JOE/TET*/VIC	DFO*/TAMRA/Cy ³ */NED	ROX			
ViiA7		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy ³ *	ROX/ATTO 590*/TR*	ATTO 647N*/HL647*/Cy ⁵	ATTO 680*/Cy ^{5.5} *	
CFX Connect & CFX Duet		FAM	YY*/HEX/TET*/JOE*					
CFX Opus & CFX96 Touch		FAM	YY*/HEX/TET*/JOE*	TR/ROX*/ATTO 590*	ATTO 647N*/Cy ⁵	ATTO 700*/Cy ^{5.5} *		
CFX384 Touch		FAM	YY*/HEX/TET*/JOE*	TR/ROX*/ATTO 590*	ATTO 647N*/Cy ⁵			
iCycler iQ™		FAM	YY*/HEX/Cy ³ /JOE/TET	ATTO 550*/DFO*/TAMRA/Cy ³	ROX/TR/ATTO 590*	ATTO 647N*/Cy ⁵		
iQ™5		FAM	YY*/HEX/Cy ³ */JOE/TET	ATTO 550*/DFO*/TAMRA/Cy ³	ROX/TR/ATTO 590*	ATTO 647N*/Cy ⁵		
QIAquant 96 & 384 5plex		FAM	YY*/HEX/JOE/VIC	ROX/TR/ATTO 590*	ATTO 647N*/Cy ⁵	Cy ^{5.5} /ATTO 680*		
QIAquant 96 2plex		FAM	YY*/HEX/JOE/VIC					
Rotor-Gene [®] 6000 & Q	ATTO 390*	FAM	YY/HEX/JOE/TET/VIC	ROX/TR/Cy ^{3.5} /ATTO 590*	ATTO 647N*/Cy ⁵ /Alexa633	ATTO 680*/Alexa680		
LightCycler [®] 480 I	ATTO 425*/Cyan 500	FAM	YY*/HEX/JOE/VIC	ROX*/TR*/LC610	ATTO 620/LC640	ATTO 633*/Cy ⁵		
LightCycler [®] 480 II	ATTO 425*/Cyan 500	FAM	YY*/HEX/JOE/VIC	ROX*/TR*/LC610	ATTO 620/LC640	ATTO 633*/Cy ⁵ /Cy ^{5.5}		
LightCycler [®] 96		FAM	YY*/HEX/VIC	TR/ROX*/ATTO 590*/LC610	ATTO 647N*/Cy ⁵			
LightCycler [®] PRO	ATTO 425*/Cyan 500	FAM	YY*/HEX/JOE/R6G/VIC	ROX*/TR*/ATTO 590*/LC610	ATTO 633*/LC640/JA 270	ATTO 647N*/Cy ⁵	Cy ^{5.5} /ATTO 700*	
Mic		FAM	YY/HEX/JOE/VIC	ATTO 590/ROX/TR	ATTO 647N/HL647/Cy ⁵			
Digital PCR Platforms								
Naica [®] Prism3		FAM	HEX/Cy ³ /VIC	ATTO 647N*/Cy ⁵				
Naica [®] Prism6		FAM	YY	ATTO 550/TAMRA	ROX	ATTO 647N*/Cy ⁵	ATTO 700/Cy ^{5.5}	
Nio & Nio+		FAM	YY/HEX*/ATTO 532*/JOE*	ATTO 550/TAMRA/Cy ³	ROX/TR*/ATTO 590*	ATTO 647N*/Cy ⁵	Cy ^{5.5} /ATTO 700*	DY-521-XL
QX One ddPCR		FAM	YY*/HEX/VIC	ATTO 647N*/Cy ⁵	Cy ^{5.5}			
QX100 & QX200 ddPCR		FAM	YY*/HEX/VIC					
QX600 ddPCR		FAM	YY*/HEX/VIC	ATTO647N*/Cy ⁵	Cy ^{5.5}	ROX	ATTO 590	
QuantStudio™ Absolute Q		FAM	HEX/YY*/VIC	ATTO 550*/TAMRA*/ABY	ROX/ATTO 590*/TR*	ATTO 633*/Cy ⁵ /JUN		
QIAcuity Four/Eight (5plex)		FAM	YY*/HEX/JOE/VIC	TAMRA/ATTO 550	ROX/TR/ATTO 590*	ATTO 647N*/Cy ⁵		
QIAcuity One (2plex)		FAM	YY*/HEX/JOE/VIC					
Digital LightCycler [®]	ATTO 425/Cyan 500	FAM	HEX/VIC	TR/CFR610/LC610	ATTO 647N/Cy ⁵ /JA270/LC640	Cy ^{5.5}		
DQ24 Plus	ATTO 425	FAM	YY*/HEX/JOE/VIC/SUN	ROX/TR	ATTO 647N*/Cy ⁵ /Q670	Cy ^{5.5} /Quasar 705		



For complementary information, please refer to instrument manufacturer technical guide or contact us at scientific.support@eurogentec.com

* For optimal results in multiplex assays, a dye calibration may be required to avoid fluorescence cross-talk

YY = Yakima Yellow[®]
DFO = Dragonfly Orange™

TR = Texas Red[®]
R6G = Rhodamine-6G

In bold = Recommended by Eurogentec
In regular black = Available at Eurogentec
In grey = Not available at Eurogentec