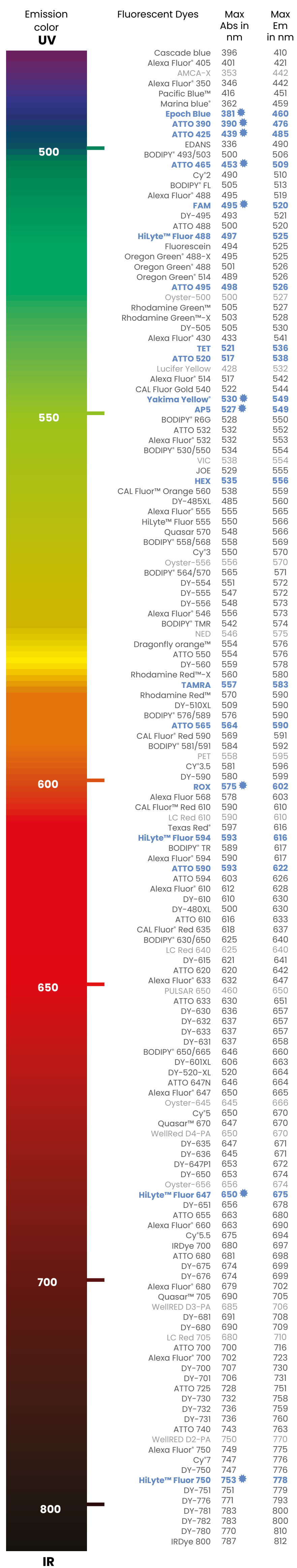
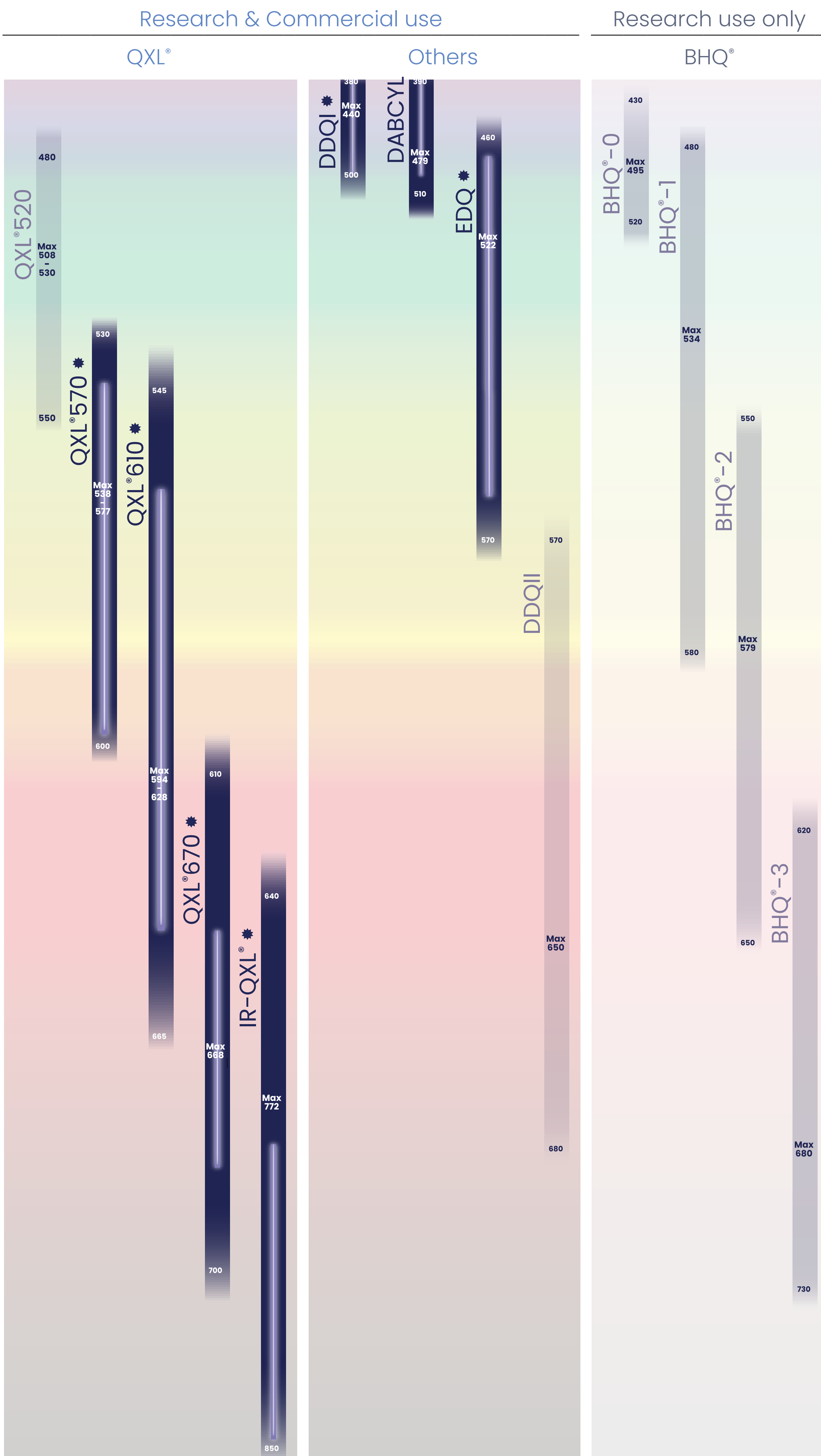


FLUORESCENT DYES



QUENCHERS



Order

Oligonucleotides can be ordered via the web oligo configurators or on request. For more information, please contact us or visit our website. [www.eurogentec.com](http://www.eurogentec.com)

High performance

Access™ Dyes & Quenchers from Eurogentec are a class of high performance proprietary molecules with IP-friendly terms and conditions.

Usage

In bold blue: Dyes & Quenchers for commercial & research use  
In black: Dyes & Quenchers for research use only (RUO)  
In light grey: Not available at Eurogentec

Dark blue bands represent the indicative quenching range of recommended quenchers  
Gray bands represent the indicative quenching range of alternate quenchers  
Purple bars represent the optimal recommended quenching range  
Access™ Dyes & Quenchers

All wavelength values are in nm.  
Note: Spectral properties of a dye or a quencher may vary when coupled to another molecule

\* Perform a dye calibration for optimal results  
For complementary information, please refer to instrument manufacturer technical guide or contact us at [scientific.support@eurogentec.com](mailto:scientific.support@eurogentec.com)

YY = Yakima Yellow®  
DFO = Dragonfly Orange™  
TR = Texas Red™  
In blue = Recommended by Eurogentec  
In black = Available at Eurogentec  
In grey = Not available at Eurogentec

COMPATIBILITY CYCLER CHANNELS VS FLUORESCENT DYES

Thermocycler	Blue channel	Channel 1	Channel 2	Channel 3	Channel 4	Channel 5	Channel 6	Channel 7
GeneAmp® SDS 5700		FAM	AP5/YV/JOE/VIC/TET*	DFO*/TAMRA/NED				
ABI Prism® SDS 7000		FAM	AP5/YV/JOE/VIC/TET*	DFO*/TAMRA/NED				
ABI Prism® SDS 7700		FAM	AP5/YV/JOE/VIC/TET*	DFO*/TAMRA/NED	ROX			
ABI Prism® SDS 7900 HT		FAM	AP5/YV/JOE/VIC/TET*	DFO*/TAMRA/NED	ROX			
ABI Prism® SDS 7300		FAM	AP5/YV/JOE/VIC/TET*	DFO*/TAMRA/NED	ROX			
ABI Prism® SDS 7500		FAM	AP5/YV/JOE/VIC/TET*	DFO*/TAMRA/NED	ROX/TR			
Quantstudio® 3.5, 6, 7		FAM	AP5/YV/HEX*/JOE/VIC/TET*	DFO*/TAMRA/cy3/NED	ROX/TR	cy5/HL647		
Via7		FAM	AP5/YV/HEX*/JOE/VIC/TET*	DFO*/TAMRA/cy3/NED	ROX/TR	cy5/HL647		
StepOne®		FAM	AP5/YV/HEX*/JOE/VIC/TET*	ROX				
StepOnePlus®		FAM	AP5/YV/HEX*/JOE/VIC/TET*	DFO*/TAMRA/cy3/NED	ROX			
Cycler IQ™		FAM	AP5/YV/HEX/cy3/TET	DFO*/TAMRA/cy3/NED	ROX/TR	cy5		
My IQ™		FAM						
iQ5		FAM	AP5/YV/HEX/JOE/TET	DFO*/TAMRA/cy3/NED	ROX/TR	cy5		
CFX 96™		FAM	AP5/YV/HEX/JOE/TET	ROX/TR	cy5			
MiniOpticon™		FAM	AP5/YV/HEX/TET*					
DNA Engine Opticon™1		FAM						
DNA Engine Opticon™2		FAM						
Chromo 4™		FAM	AP5/YV/HEX/TAMRA/VIC/TET					
Mx3000P (choice of 4 filters)		FAM	TET	YY*/HEX/JOE/VIC	cy3	DFO/TAMRA/NED		TR/ROX
Mx3005P (choice of 5 filters)	Epoch Blue / Alexa Fluor® 350	FAM	TET	YY*/HEX/JOE/VIC	cy3	DFO/TAMRA/NED		cy5
Mx4000™ (choice of 4 filters)		FAM	TET	YY*/HEX/JOE/VIC	cy3	DFO/TAMRA/NED		TR/ROX
Aria™ (choice of up to 6 filters)	ATTO 425	FAM	HEX	cy3	ROX	cy5		cy5
Mastercycler™ep realplex2		FAM	AP5/YV/HEX/JOE/VIC/TET					
Mastercycler™ep realplex4		FAM	AP5/YV/HEX/TET/JOE/VIC	DFO*/TAMRA	ROX			
LightCycler™ 1.5		FAM	AP5/YV/HEX/JOE/VIC/TET	cy5				
LightCycler™ 2		FAM	AP5/HEX/YV/JOE*/VIC	ROX/TR/IC Red 610	ATTO 620/IC Red 640	cy5/IC Red 670		ATTO 680/IC Red 705
LightCycler 480™ I	ATTO 425	FAM	AP5/YV/HEX/JOE*/VIC	ROX/TR/IC Red 610	ATTO 620/IC Red 640	cy5		
LightCycler 480™ II	ATTO 425	FAM	AP5/YV/HEX/JOE/VIC	ROX/TR/IC Red 610	ATTO 620/IC Red 640	cy5 / cy5.5 / IC Red 705		
LightCycler 96		FAM	AP5/YV/HEX/VIC	TR/IC Red 610	cy5			
Smartcyler™1		FAM	AP5/YV/cy3/JOE/VIC/TET	TR	cy5			
Smartcyler™2		FAM	AP5/cy3/YY*/JOE*/TET	TR/ROX	cy5			
Rotor-Gene™ 2000 / 3000		FAM	AP5/YV/JOE/VIC/TET	TAMRA/ROX/cy3.5/TR				
Rotor-Gene™ 6000	Epoch Blue / Alexa Fluor® 350	FAM	AP5/YV/HEX/JOE/VIC/TET	TAMRA/ROX/cy3.5/TR			ATTO 680	
PikoReal		FAM	AP5/HEX/YY	ROX/TR	cy5			
Qtower		FAM	AP5/YV/HEX/JOE/VIC	DFO/TAMRA/NED	ROX/TR/cy3.5*	cy5		cy5.5