

Compatibility of plastic plates and tubes with Real-Time PCR thermocyclers

		QuantStudio w/ standard (0.2 mL) 96-well blocks (1, 3, 5, 6&7 Flex/PRO, 12k Flex)	QuantStudio w/ FAST (0.1 mL) 96-well blocks (3, 5, 6&7 Flex/PRO, 12k Flex)	QuantStudio w/ 384-well blocks (5, 6&7 Flex/PRO, 12k Flex)	Agilent AriaMx & AriaDx	Analytik Jena qTOWER ³ & qTOWERiris	Bio-Rad iCycler iQ-iQ5 & MyiQ™	Bio-Rad CFX96 & CFX Opus/Connect 96 & CFX Duet	Bio-Rad CFX384 & CFX Opus/Connect 384	Qiagen QIAquant (2- & 5-plex)	Roche LightCycler® 480 (I & II)	Roche LightCycler® 96	ABI Prism® SDS 7500	ABI Prism® SDS 7900HT	ABI StepOne™ Plus	ABI Viia7	Regular block	FAST block	Regular block	FAST block	Regular block	FAST block	
qPCR 96-well plate type	Cover																						
Sub-skirted, high profile, frosted or white type (ABI systems) RT-PL96-AB or RT-PL96-ABW	Flat caps RT-FLAT-120	★	-	-	-	-	✓	-	-	-	-	-	-	-	★	-	-	★	-	-	★	-	
	Adhesive seals RT-OPSL-100	★	-	-	-	-	✓	-	-	-	-	-	★	-	★	-	-	★	-	-	★	-	
Sub-skirted, low profile, frosted or white type (ABI systems) RT-PL96-AF or RT-PL96-AFW	Adhesive seals RT-OPSL-100	-	★	-	★	-	-	-	-	-	-	-	-	★	-	★	★	-	-	-	-	★	
	Flat caps RT-FLAT-120	✓	-	-	-	✓	★	-	-	-	-	-	✓	-	-	-	-	✓	-	-	✓	-	
Non-skirted, high profile, natural or white type RT-PL96-MQ or RT-PL96-MQW	Adhesive seals RT-OPSL-100	✓	-	-	-	✓	★	-	-	-	-	-	✓	-	-	-	-	✓	-	-	✓	-	
	Flat caps RT-FLAT-120	-	✓	-	★	-	-	★	-	★	-	★	-	✓	-	✓	✓	-	-	-	-	-	
Non-skirted, low profile, frosted or white type RT-PL96-OP or RT-PL96-OPW	Adhesive seals RT-OPSL-100	-	✓	-	★	-	-	★	-	★	-	★	-	✓	-	✓	✓	-	-	-	-	-	
	Flat caps RT-FLAT-120	-	-	-	-	-	-	★	-	★	★	-	-	-	-	-	-	-	-	-	-	-	
Semi-skirted, low profile, white RT-PL96-LC480	Adhesive seals RT-OPSL-100	-	-	-	-	-	-	★	-	★	★	-	-	-	-	-	-	-	-	-	-	-	
	Flat caps RT-FLAT-120	-	-	-	-	-	-	★	-	★	★	-	-	-	-	-	-	-	-	-	-	-	
qPCR 384-well plate type	Cover																						
Skirted, low profile, white plate for ABI RT-PL384-ABW	Adhesive seals RT-OPSL-100	-	-	★	-	-	-	-	★	★	-	-	-	-	★	★	-	★	-	-	★	★	
	Adhesive seals RT-OPSL-100	-	-	-	-	-	-	-	★	-	★	-	-	-	-	-	-	-	-	-	-	-	
qPCR 8-tube strip type	Cover																						
Extra Robust, Regular Profile, Frosted, 0.2 ml RT-STR8-RF120	Flat caps RT-FLAT-120	✓	-	-	-	★	✓	-	-	-	-	✓	✓	-	✓	-	-	✓	-	-	✓	-	
	Flat caps RT-FLAT-120	✓	-	-	-	★	✓	-	-	-	-	✓	✓	-	✓	-	-	✓	-	-	✓	-	
Extra Robust, Low Profile, Frosted, 0.1 ml RT-STR8-LF120	Flat caps RT-FLAT-120	-	✓	-	✓	-	-	✓	-	★	▲	-	-	✓	-	✓	✓	-	-	-	✓	✓	
	Flat caps RT-FLAT-120	-	✓	-	✓	-	-	✓	-	★	▲	-	-	✓	-	✓	✓	-	-	-	✓	✓	

★ recommended for optimal compatibility | ▲ adaptor required (Ref: RT-PL96-LCAD) | ✓ compatible plastics | ✓ Refer to cyclers manufacturer instructions for required adaptor
 qPCR optical semi-domed 8-cap strips are also available (Ref: RT-OCAP-300: 300 cap strips) and can be used on several machines.