

## 96-well plate compatibility with Real-Time thermocycler

		ABI GeneAmp® 5700	ABI Prism® SDS 7000	ABI Prism® SDS 7300	ABI Prism® SDS 7500	ABI Prism® SDS 7500 FAST	ABI Prism® SDS 7700	ABI Prism® SDS 7900HT	ABI Prism® SDS 7900 FAST	iCycler IQ®	My iQ	iQ5	Mx3000P®	Mx3005P®	MasterCycler® ep realplex	DNA Engine Opticon® 1	DNA Engine Opticon® 2	Chromo 4	MiniOpticon®	Quantica®	Reference
qPCR 96-well plate sub-skirted, high profile, natural (ABI systems)	25 (5 X 5)	✓ *	✓ *	✓ *	✓ *		✓ *	✓ *													RT-PL96-AB
qPCR 96-well plate sub-skirted, low profile, natural (ABI FAST systems)	25 (5 X 5)					✓			✓												RT-PL96-AF
qPCR 96-well plate non-skirted, high profile, natural	25 (5 X 5)	✓	✓	✓	✓		✓			✓	✓	✓	✓	✓	✓						RT-PL96-MQ
qPCR 96-well plate non-skirted, low profile, natural	25 (5 X 5)															✓	✓	✓	✓	✓	RT-PL96-OP
qPCR optical semi-dome 8-cap strip, natural	300	✓ *	✓ *	✓ *	✓ *		✓ *	✓ *		✓ **	✓ **	✓ **									RT-OCAP-300
qPCR optical flat 8-cap strip, natural	300	✓ *	✓ *	✓ *	✓ *		✓ *	✓ *		✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	RT-FLAT-300
qPCR opti-seals (optical, disposable, adhesive)	25	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	RT-OPSL-25
qPCR opti-seals (optical, disposable, adhesive)	100	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	RT-OPSL-100

(\*) When qPCR 96-well plates for ABI Systems, reference RT-PL96-AB are closed with semi-dome caps the height of caps and rim is the same, so the pressure of the lid is applied equally and a good closure is achieved, leading to minimum popping open of lids. When qPCR 96-well plates for ABI Systems, reference RT-PL96-AB are closed with flat caps a maximal signal is obtained as the optical area of a flat cap is bigger than of a domed cap. So depending on your need both cap types can be used.

(\*\*) Semi-dome caps should fit on iCycler IQ®, iQ5, My iQ® when the plate is properly closed. Our recommendation is to use flat caps.