

Standard PCR

Tailor-made solutions for a broad range of applications

Product	Cat.-Nr.	Convenience	Yield	Specificity	Fidelity	Application
Ready-to-Use Mixes / direct gel loading						
Red Load Taq Master	PCR-108	++++	++	++	+	<ul style="list-style-type: none"> • Routine PCR / optimized for minimal by-product formation • Plate based PCR and automated pipetting
Red Load Taq Master / high yield	PCR-106	++++	+++	+	+	<ul style="list-style-type: none"> • Routine PCR / optimized for high efficiency in a broad range of reaction conditions • Not recommended for plate based PCR and automated pipetting
Red Load Hot Start Master	PCR-109	++++	++	+++	+	<ul style="list-style-type: none"> • High specificity PCR / high sensitivity PCR • Plate based PCR and automated pipetting
Ready-to-Use Mixes						
Taq Master	PCR-102	+++	++	++	+	<ul style="list-style-type: none"> • Routine PCR / optimized for minimal by-product formation • Plate based PCR and automated pipetting
Taq Master / high yield	PCR-101	+++	+++	+	+	<ul style="list-style-type: none"> • Routine PCR / optimized for high efficiency in a broad range of reaction conditions • Not recommended for plate based PCR and automated pipetting
Hot Start Master	PCR-103	+++	++	+++	+	<ul style="list-style-type: none"> • High specificity PCR / high sensitivity PCR • Plate based PCR and automated pipetting
Core Kits						
Taq Core Kit	PCR-232	++	++	++	+	<ul style="list-style-type: none"> • Routine PCR / optimized for minimal by-product formation • Plate based PCR and automated pipetting
Taq Core Kit / high yield	PCR-231	++	+++	+	+	<ul style="list-style-type: none"> • Routine PCR / optimized for high efficiency in a broad range of reaction conditions • Not recommended for plate based PCR and automated pipetting
Hot Start Core Kit	PCR-233	++	++	+++	+	<ul style="list-style-type: none"> • High specificity PCR / high sensitivity PCR • Plate based PCR and automated pipetting
High Fidelity Core Kit	PCR-234	++	+++	++	+++	<ul style="list-style-type: none"> • High fidelity PCR / long range PCR up to 30 kb • Amplification of GC-rich, very long and other difficult templates
Pfu-X Core Kit	PCR-237	++	+++	++	++++	<ul style="list-style-type: none"> • Amplification with highest fidelity for critical applications • High speed amplification of difficult and long templates

Standard PCR

Tailor-made solutions for a broad range of applications

Thermophilic Polymerases						
Taq Pol	PCR-202	+	++	++	+	<ul style="list-style-type: none"> Routine PCR / optimized for minimal by-product formation Plate based PCR and automated pipetting
Taq Pol / high yield	PCR-201	+	+++	+	+	<ul style="list-style-type: none"> Routine PCR / optimized for high efficiency in a broad range of reaction conditions Incorporation of labeled or other modified nucleotides
Hot Start Pol	PCR-203	+	++	+++	+	<ul style="list-style-type: none"> High specificity PCR / high sensitivity PCR Plate based PCR and automated pipetting
High Fidelity Pol	PCR-204	+	+++	++	+++	<ul style="list-style-type: none"> High fidelity PCR / long range PCR up to 30 kb Amplification of GC-rich, very long and other difficult templates
Pfu-X Polymerase	PCR-205	+	+++	++	++	<ul style="list-style-type: none"> Amplification with highest fidelity for critical applications High speed amplification of difficult and long templates
Sequencing Pol	PCR-206	+	++	++	+	<ul style="list-style-type: none"> Incorporation of dNTPs and ddNTPs at equal rates DNA Cycle sequencing
Ready-to-Use Lyophilisates						
Red Load Taq Master Lyophilisate	PCR-151	+++++	++	++	+	<ul style="list-style-type: none"> Preloaded tubes and plates for routine PCR, room temperature stable Direct loading of the PCR reaction product onto the gel
Taq Master Lyophilisate	PCR-152	++++	++	++	+	<ul style="list-style-type: none"> Preloaded tubes and plates for routine PCR Room temperature stable
Hot Start Master Lyophilisate	PCR-153	++++	++	+++	+	<ul style="list-style-type: none"> Preloaded tubes and plates for high specificity / high sensitivity PCR Room temperature stable
Supplements						
Blue Loading Buffer	PCR-254-bl	-	-	-	-	<ul style="list-style-type: none"> Facilitates loading of DNA samples into wells of agarose and polyacrylamide gels For DNA fragments larger than 100 bp, contains bromophenol blue and xylene cyanol
Orange Loading Buffer	PCR-254-or	-	-	-	-	<ul style="list-style-type: none"> Facilitates loading of DNA samples into wells of agarose and polyacrylamide gels For small DNA fragments up to 200 bp, contains Orange G
PCR Optimization Kit	PCR-251	-	-	-	-	<ul style="list-style-type: none"> Easy optimization of the MgCl₂ concentration for difficult primer-template pairs Identification of optimal enzyme / buffer / enhancer combinations
PCR Additives Kit	PCR-252	-	-	-	-	<ul style="list-style-type: none"> Enhancer for GC-rich and other difficult templates Taq Stabilizer to increase amplification yield
PCR Control Kit	PCR-253	-	-	-	-	<ul style="list-style-type: none"> Quality and benchmark tests / positive control of PCR reactions / internal lab standard Includes primer and template for amplification of the -actin gene from human DNA