



Crystal Taq Master (2x)

Master mix for routine PCR
Ready-to-Use Mixes for PCR

Cat. No.	Amount
PCR-166S	4 x 1,25 ml (2x conc.)
PCR-166L	20 x 1.25 ml (2x conc.)
PCR-166XL	100 ml (2x conc.)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: Short term storage (up to 3 month) at 4 °C possible.

Shelf Life: 12 months

Form: liquid

Concentration: 2x conc.

Description:

Crystal PCR Master is a 2 x conc. ready-to-use master mix recommended for routine PCR applications (up to 4 kb fragment length), high throughput PCR or genotyping.

It contains all reagents required for PCR (except template and primer) in a well-balanced ratio to ensure high specificity and minimal by-product formation in almost all PCR applications without the need of additional optimization steps.

The mix guarantees robust and reliable amplification results with a minimum of pipetting steps, saves time and reduces the risk of contaminations.

The total PCR assay volume is freely adaptable to individual protocols or the requirements of automated pipetting systems.

Content:

Cat.No.	Master Mix	PCR-grade water	Assays x 50 µl
PCR-166S	4 x 1.25 ml	6 ml	200
PCR-166L	20 x 1.25 ml	2 x 12 ml	1000
PCR-166XL	100 ml	100 ml	4000

2 x concentrated PCR master mix containing Taq polymerase, nucleotides (dATP, dCTP, dGTP, dTTP), KCl, (NH₄)₂SO₄, MgCl₂, density reagent, enhancing and stabilizing additives.

Recommended PCR assay:

Before starting, vortex the master mix thoroughly to assure homogeneity.

component	stock conc.	20 µl assay	50 µl assay	final conc.
Crystal PCR Master	2x	10 µl	25µl	1x
Primer Mix or each primer	10 µM each primer	0.4-0.8 µl	1-2 µl	200-400 nM each primer
Template/sample DNA		< 10 ng	< 20 ng	
PCR-grade water		fill up to 20 µl	fill up to 50 µl	

Recommended cycling conditions:

Before cycling, vortex PCR tubes or plates to assure homogeneity and centrifuge briefly to remove bubbles.



Crystal Taq Master (2x)

Master mix for routine PCR
Ready-to-Use Mixes for PCR

Initial denaturation	95 °C	2 min	1x
Denaturation	95 °C	10 - 20 sec	25 - 30x
Annealing ¹⁾	50 - 68 °C	10 - 20 sec	
Elongation ²⁾	72 °C	20 sec - 4 min	

¹⁾The annealing temperature depends on the melting temperature of the primers used.

²⁾The elongation time depends on the length of the fragments to be amplified. A time of 1 min/kb is recommended.