



Crystal Buffer

Cat. No.	Amount
PCR-271S	5 x 1,2 ml
PCR-271	5 x 1,2 ml
PCR-271-10ML	10 ml
PCR-271-100ML	100 ml

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C or -20 °C

Shelf Life: 12 months

Form: Liquid

Concentration: 10x conc.

Description:

Crystal PCR Buffer guarantees robust and reliable amplification results in almost all PCR applications. The buffer contains a well-balanced ratio of potassium-, ammonium- and magnesium-ions to ensure high specificity and minimal by-product formation without the need of additional optimization steps.

Content:

Crystal PCR Buffer (green cap)

200 mM Tris-HCl, KCl, $(\text{NH}_4)_2\text{SO}_4$ and 20 mM MgCl_2 .

Optimization of MgCl_2 concentration:

A final Mg^{2+} concentration of 2.0 mM is recommended for most PCR applications. For an individual optimization of the reaction buffer add 25 mM MgCl_2 stock solution (PCR-266) as shown in the table below.

final MgCl_2 conc.	20 μl final assay volume	50 μl final assay volume
2 mM	-	-
3 mM	0.8 μl	2.0 μl
4 mM	1.6 μl	4.0 μl
5 mM	2.4 μl	6.0 μl