

**qPCR Buffer**

Optimized buffer system for real-time PCR

Cat. No.	Amount
PCR-279S	5 x 1,2 ml
PCR-279-10ML	10 ml
PCR-279-100ML	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:**

qPCR Buffer is optimized for use in real-time PCR applications. The buffer contains a well-balanced ratio of potassium-, ammonium- and magnesium-ions to ensure high specificity and minimal by-product formation in probe-based assays as well as in SybrGreen/EvaGreen-based assays. There is no additional optimization required for most assay set-ups.

Content:**qPCR Buffer (green cap)**200 mM Tris-HCl, KCl, $(\text{NH}_4)_2\text{SO}_4$ and 30 mM MgCl_2 .**Optimization of MgCl_2 concentration:**

MgCl_2 Solution - 25 mM (#PCR-266) is recommended for optimization of the final Mg^{2+} concentration. A concentration of 2 - 5 mM is required for optimal functionality. Lower concentrations give higher specificity, whereas higher concentrations generate a higher yield.