

New products & catalog numbers from November 1st!



Endpoint PCR

Taq and Hot Start Polymerases come now with a re-optimized Crystal & Ruby Buffer system for reliable amplification results in PCR applications.



Thanks to a well-balanced ratio of potassium-, ammonium- and magnesium-ions **Crystal Buffer** ensures high specificity and minimal by-product formation. No more optimization steps are required.



Ruby Buffer additionally contains gel loading buffer and an inherent red dye. The red dye allows an easy visual control during PCR set-up and in combination with the density reagent the direct loading of the reaction product into the gel.

Product	Previous Cat.-No.	New Cat.-No.
Taq Polymerase	PCR-202	PCR-211
Hot Start Polymerase	PCR-203	PCR-212
Hot Start Polymerase Ab+	PCR-210	PCR-213
Taq Core Kit	PCR-232	PCR-214
Hot Start Core Kit	PCR-233	PCR-215
Hot Start Core Kit Ab+	PCR-239	PCR-216