

EpiQuik Red Taq 2x PCR Mix

(Catalog No. R12020)

Description

EpiQuik Red Taq 2x PCR Mix is a premixed solution containing everything needed for successful PCR reaction except specific primers and DNA template. The mix includes high-quality recombinant EpiQuik Red Taq Hot Thermostable DNA polymerase, nucleotides and magnesium in a PCR reaction buffer. The mix has a special additive - the inert red dye. This makes it very suitable for standard applications. Strong red color of the mix allows user to verify adequate mixing. Reaction products are ready for direct gel loading and the dye serves as a marker for electrophoresis progress monitoring. For the reaction set-up, add the PCR Mix (10 or 25 µl) to the primers, template and water for the total reaction volume of 20 or 50 µl. To activate the EpiQuik Red Taq Hot Thermostable DNA polymerase, it should be incubated at 95 – 97°C for 15 minutes as a first PCR step. This enzyme mix allows the PCR setup at ambient temperature without nonspecific annealing and extension.

Composition

- EpiQuik Red Taq DNA Polymerase (0.1 U/µl) in reaction buffer, 4 mM MgCl₂, 0.4 mM of each dNTPs (dATP, dCTP, dGTP, dTTP)
- Water (nuclease free) - 1 ml
- 1 ml of mix is sufficient for 100 reactions

Applications

- Ready-to-use mix reduces experimental variability
- Reduced pipetting steps, thereby less time-consuming set-up of PCR
- Set-up at room temperature
- Visualization of complete reaction mixing
- Direct loading of samples following amplification

Quality Data

Activity and stability tested at 20, 30 and 40 cycles of PCR reactions at 95°C. Free of detectable, nonspecific nucleases. Several cycles of freezing/thawing are allowed.

Safety Warnings and Precautions

This product is designed for research purposes and in vitro use only. According to common laboratory safety practice, it is recommended to wear protective clothing, gloves and safety glasses. Please refer to www.epigentek.com for Material Safety Data Sheet of the product. Some applications this product is used in may require a license which is not provided by the purchase of this product. Users should obtain the license if required.

Storage Conditions

Store at -20°C. Guaranteed stable for 1 year when properly stored.

Ordering Information

Products

EpiQuik Red Taq 2x PCR Mix

Size

1 ml

Cat. No.

R12020-1

Recommended PCR reaction mix:

Component	Quantity
EpiQuik Red Taq 2x PCR Mix	10 µl (1x)
Primer Forward	0.3-1 µM
Primer Reverse	0.3-1 µM
DNA Template	1-100 ng/µl
H ₂ O PCR Grade	Up to 20 µl
Total	20 µl

Recommended PCR cycles:

Cycle step	Temp.	Time	Cycles
Initial Denaturation	95°C	15 min	1
Denaturation	95°C	30-60 s	26-35
Annealing	50-68°C	30-60 s	26-35
Elongation	72°C	1-4 min	26-35
Final Elongation	72°C	5-10 min	1

IMPORTANT: Annealing temperature should be 2-6°C lower than the primer melting temperature.

This product is for research purposes only. Not intended for use in diagnostic procedures.