

EpiQuik Taq 2x PCR Mix

(Catalog No. R12018)

Description

EpiQuik 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 Taq DNA polymerase, nucleotides and magnesium in a PCR reaction buffer. For the reaction set-up, add the PCR Mix (10 or 25 μ l) to the primers, template and water for the total reaction volume of 20 or 50 μ l.

Composition

- EpiQuik Taq DNA Polymerase(0.1 U/ μ l) 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

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.

Recommended PCR reaction mix:

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

Recommended PCR cycles:

Cycle step	Temp.	Time	Cycles
Initial Denaturation	95°C	3-5 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.

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 Taq 2x PCR Mix

Size

1 ml

Cat. No.

R12018-1

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